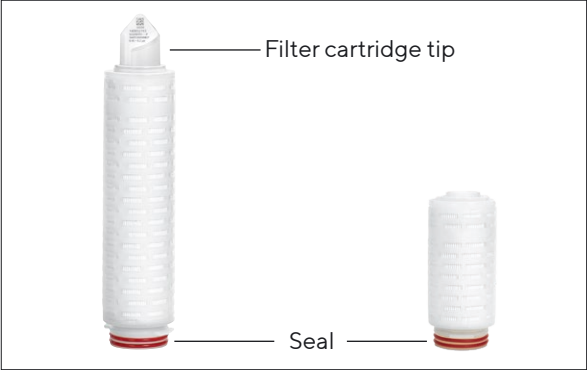




Instructions for Use

Standard and Mini Filter Cartridges



1. Validity

These instructions apply for the following product lines:

- Aerosart®
- Aquasart® Plus | Plus XL | PS
- Sartobev® PS
- Sartobran® P
- Sartoclean® CA | GF
- Sartocool® PS
- Sartofluor® GA | LG | HR
- Sartoguard GF | PES
- Sartolon®
- Sartopore® 2 | 2 XLG | 2 XLI | 2 XLM | Air | Platinum | Platinum HB
- Sartopore® Evo
- Sartopure® GA | GF Plus | IND | PP3
- Sartovent®
- Vinosart® PS

2. Safety Information

There are no inherent dangers associated with Standard and Mini filter cartridges. However, specific application scenarios may pose risks in connection with the technical systems in which the filter cartridges are operated. The warning and safety instructions in the instructions for use for the system in which the standard and mini filter cartridges are used must therefore be observed. This document only applies in combination with the system’s instructions for use.

3. Labeling

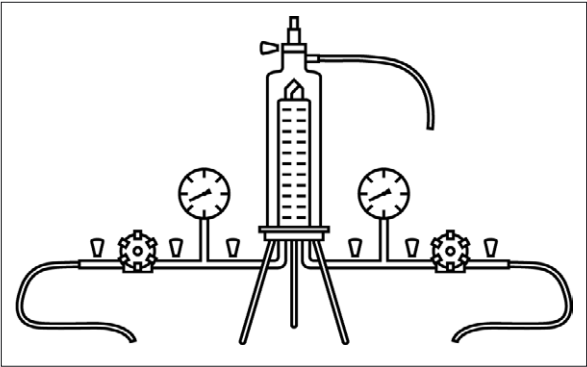
Information on filter cartridge tip on the label of the plastic bag and the outer packaging:

- Product designation
- Pore size or rejection rate
- Batch number
- Individual item number
- Order no.

Additional information on the label of the plastic bag and the outer packaging:

- Expiry date
- Batch number

4. Typical Filtration Setup with Stainless Steel Housing



5. Inserting Filter Cartridges into Housing

- Cut open the plastic bag at the bottom end on the seal side.
- Push the filter cartridge far enough out of the plastic bag so that the seals can be well moistened.
- Moisten the seals with water or the liquid to be filtered.
- Hold the filter cartridge firmly at the bottom end and insert into the underside of the housing by turning gently.
- Remove the plastic bag.
- Mount the housing dome on the base.

6. Moistening the Filter Unit

Filter units must be moistened before the integrity test.

Filter	Liquid for moistening
Hydrophilic filters	Aqueous solution
PTFE membrane filter	Liquid with surface tension of < 28 dyn/cm

- Close all valves.
- Open the vent valve until liquid comes out of it with no air bubbles.
- When the liquid emerges from the vent valve, stop filling.

- Close the vent valve.
- Set a differential pressure of 0.3 bar.
- Rinse for 3 – 5 minutes.
- To empty, open the vent valve and the discharge valve until no more liquid comes out of the housing.
- Close the valves.

7. Autoclaving the Filter Unit

- Sterilization conditions: Up to 134 °C, 2 bar, 30 min.
- Place the filter unit in an autoclavable sterilization bag.
 - Once the sterilization temperature has been reached, continue autoclaving the filter element for a further 30 minutes.
 - After autoclaving, remove the filter unit from the sterilization bag.

8. Performing In-Line Steam Sterilization

- Sterilization conditions: Up to 134 °C, 0.3 bar, 20 min. Perform in-line steam sterilization of the filter cartridge in the filtration direction.
- Throughout the entire procedure, apply differential pressure of max 0.5 bar, 121 °C or max. 0.3 bar, 134 °C.
 - Open the vent valve and the drain valve a little or use a steam separator.
 - Let the steam flow in slowly, by opening the inlet valve slowly.
 - If steam is escaping from all valves, leave at least one valve slightly open on the sterile side so that the steam can flow or use a steam separator. Close the other valves.

Error Prevention	Procedure
Preventing vacuum condensation	Blanket the filter system with compressed air on the inlet and outlet side following steam sterilization. The compressed air should be approx. 200 mbar higher than the previously applied sterilization pressure.
Cooling off filter cartridge	Flush system with hydrophilic filters with water or buffer until the system reaches ambient temperature. Hydrophobic filters can be dried with flowing air.

9. Performing Sanitization with Hot Water

- Open the vent valve and leave open during filling.
- Heat water to 85 °C.
- Completely fill the housing with water and allow to circulate for 30 minutes.
- Turn off the hot water supply.

10. Performing Integrity Tests

- Requirement: Filter unit is moistened, sterilized and cooled to room temperature.
- Moisten the filter unit again.
 - Perform the integrity test with an automatic test device, e.g., with Sartocheck®. Observe the Validation Guide and QAC for the IT values.

11. Performing Liquid Filtration

- Fully vent the housing. To do so, open the vent valve until some liquid escapes.
- Close the vent valve again.
- Set the filtration target pressure.
- Perform the filtration.

12. Returning Used Filter Elements

- In the event of a complaint, the product can be returned to Sartorius. Returned products must be clean, decontaminated by sterilization, and packaged. Products contaminated with hazardous materials, e.g., biological or chemical substances that pose a risk to human health, will not be accepted for disposal.
- Contact your local Sartorius contact person for instructions on how to return parts (see www.sartorius.com).
 - Pack the product for return.

13. Disposal

- The product must be disposed of properly by disposal facilities. The product must be decontaminated by sterilization for this purpose.
- Dispose of the product in accordance with local government regulations.
 - Dispose of the packaging in accordance with local government regulations.

14. Application Support

If needed, you can request the corresponding documents with further technical data, information on how to perform product integrity tests, or specific information about additional applications from your local Sartorius contact person.

Instructions for Use

Maxicaps®, Midicaps®, and Capsules



1. Validity

These instructions apply for the following product lines:

- Aerosart®
- Aquasart® Plus | Plus XL | PS
- Sartobev® PS
- Sartobran® P
- Sartoclean® CA | GF
- Sartocool® PS
- Sartofluor® GA | LG | HR
- Sartoguard GF | PES
- Sartolon®
- Sartopore® 2 | 2 XLG | 2 XLI | 2 XLM | Air | Platinum | Platinum HB
- Sartopore® Evo
- Sartopure® GA | GF Plus | IND | PP3
- Sartovent®
- Vinosart® PS

2. Safety Information

There are no inherent risks associated with Maxicaps®, Midicaps®, and capsules. However, specific application scenarios may pose risks in connection with the technical systems in which Maxicaps®, Midicaps® or capsules are operated. The warning and safety instructions in the instructions for use for the system in which the Maxicaps®, Midicaps®, or capsules are used must therefore be observed. This document only applies in combination with the system’s instructions for use.

3. Labeling

Information on the housing of the Maxicaps®, Midicaps®, and the capsules:

- Product designation
- Pore size or rejection rate
- Batch number
- Individual item number
- Order no.
- Operating parameters
- Direction of filtration
- Expiry date

Information on the label of the plastic bag | the outer packaging:

- Product designation
- Pore size or rejection rate
- Expiry date
- Batch number
- Order no.

4. Moistening the Filter Unit

Filter units must be moistened before the integrity test.

Filter	Liquid for moistening
Hydrophilic filters	Aqueous solution

PTFE membrane filter	Liquid with surface tension of < 28 dyn/cm
----------------------	--

- Close all valves.
- Open the vent valve until liquid comes out of it with no air bubbles.
- When the liquid emerges from the vent valve, stop filling.
- Close the vent valve.
- Set a differential pressure of 0.3 bar.
- Rinse for 3–5 minutes.
- To empty, open the vent valve and the discharge valve until no more liquid comes out of the housing.
- Close the valves. Ensure that the valves of Maxicaps® and Midicaps® are not closed against pressure exceeding 2 bar.

5. Autoclaving | Gamma Irradiating the Filter Unit

Depending on the filter unit used, only one sterilization method can be used: gamma irradiation or autoclaving. Products cannot be autoclaved after gamma irradiation.

Autoclaving

- Requirement: hydrophilic membrane filters must be moistened (exception is Sartopore® Platinum)
- Sterilization conditions: Up to 134 °C, 2 bar, 30 min.
- Loosely connect the stainless steel adapter with the filter unit connection.
 - Place the filter unit with the stainless steel adapter in an autoclavable sterilization bag.
 - Once the sterilization temperature has been reached, continue autoclaving the filter element for a further 20 minutes.
 - After autoclaving, remove the filter unit from the sterilization bag and tighten the stainless steel adapter by hand.

Gamma Irradiation

- Use a gamma-compatible capsule.
- Perform gamma irradiation of the capsule at max. 50 kGy.

6. Performing Integrity Tests

- The filter unit has been moistened, sterilized, and cooled to room temperature.
- Perform the integrity test with an automatic test device, e.g., with Sartocheck®. Observe the Validation Guide and QAC for the IT values.

7. Performing Liquid Filtration

- Fully vent the housing.
- Open the vent valve until the liquid to be filtered escapes.
- Close the vent valve again.
- Set the filtration target pressure.
- Perform the filtration.

8. Returning Used Filter Elements

- In the event of a complaint, the product can be returned to Sartorius. Returned products must be clean, decontaminated by sterilization, and packaged. Products contaminated with hazardous materials, e.g., biological or chemical substances that pose a risk to human health, will not be accepted for disposal.
- Contact your local Sartorius contact person for instructions on how to return parts (see www.sartorius.com).
 - Pack the product for return.

9. Disposal

- The product must be disposed of properly by disposal facilities. The product must be decontaminated by sterilization for this purpose
- Dispose of the product in accordance with local government regulations.
 - Dispose of the packaging in accordance with local government regulations.

10. Application Support

If needed, you can request the corresponding documents with further technical data, information on how to perform product integrity tests, or specific information about additional applications from your local Sartorius contact person.

Sartorius Stedim Biotech GmbH
August-Spindler-Str. 11
37079 Goettingen
Phone +49 551 308 0
www.sartorius.com

© 2025
Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
ML | Publication No.: SPK9077-e250302
DIR: 3030658-000-02
Status: 03 | 2025

List of Sartorius material numbers applying to EPA-FIFRA

16702–27	5102507H4	5107307H1-SS
16702–28	5102507H4--EP	5107307H1-SSIT
16703–79	5102607H2–E	5107307H2-FF
16TPB002-05	5102607H2–EIVG	5107307H2-OO
4-FLEX	5102707H1	5107307H2-OOIT
5101507H9	5102707H2	5107307H2-SO
5101507H9--B	5102707H3	5107307H2-SOIT
5102107H1	5102807H1	5107307H2-SS
5102107H2	5102807H3	5107307H2-SSIT
5102107H2–E	5105307H9-FF-A	5107307H3-FF
5102507H1	5105307H9-OO-1	5107307H3-OO
5102507H1--IVG	5105307H9-OO-A	5107307H3-OOIT
5102507H1–E	5105307H9-SO-1	5107307H3-SO
5102507H1–ELE	5105307H9-SO-A	5107307H3-SOIT
5102507H1–X	5105307H9-SS-A	5107307H3-SS
5102507H2	5105307H9-SSVPA	5107307H3-SSIT
5102507H2--IVG	5105307H9-EOO-A	5152507T0--EA
5102507H2–E	5105307H9-ESO-A	5152507T1--C
5102507H2–EIVG	5105307H9-ESS-A	5152507T1--EA
5102507H2–ELE	5107307H1-FF	5152507T1–E
5102507H3	5107307H1-OO	5152507T1–E-C
5102507H3--IVG	5107307H1-OOIT	5152507T2--C
5102507H3–E	5107307H1-SO	5152507T2--EA
5102507H3–X	5107307H1-SOIT	5152507T2–E

5152507T2-E-C	5181507T7--B	5182107T1--A-K
5152507T3--EA	5181507T7--GPN	5182107T1--GA
5152507T3-E-C	5181507T7-E-B	5182107T1--GF
5152707T1-E	5181507T7-EGPN	5182107T1--LG
5152707T1-E-C	5181507T8--B	5182107T1-FULG
5152707T2-E	5181507T8--MDB	5182158T1--A-K
5152707T2-E-C	5181507T8-E-B	5182158T1-FULG
5152807T1-E	5181507T9--B	5182506T0--GA
5152807T1-E-C	5181507T9--EAB	5182506T1--GA
5152807T2-E	5181507T9-E-B	5182506T1-EGA
5152807T2-E-C	5181507T9-EK-B	5182506T2--GA
5172507T1--HR	5181507T9-V-B	5182506T3--GA
5172507T2--HR	5181507T9-25-B	51825070-TCA
5172507T3--HR	5181507T9-25BXB	5182507T0--GA
5181207T7--B	5181807T6	5182507T0--GAG
5181207T8--B	5181807T6--D	5182507T0-EGA
5181207T8--MTB	5181807T6--K	5182507T0-TCA
5181207T8--STB	5181807T6--EAA	5182507T0-TGA
5181207T9--B	5181807T7--B	5182507T0-TLE
5181207T9--MTB	5181807T7--KSH	5182507T0-VGA
5181307T4-OO-B	5181807T8--B	5182507T1--F2
5181307T4-SO-B	5181807T8-E-B	5182507T1--F3
5181307T4-SS-1	5181807T9--B	5182507T1--GA
5181307T4-SS-B	5182106T1-FULG	5182507T1--GAG
5181307T5-OO-D	5182107T0--A-K	5182507T1--LG
5181407T7--B	5182107T0--GF	5182507T1--LGG

5182507T1--SC	5182507T3--EGA	5182707T3--LG
5182507T1--EGA	5182507T3--ELG	5182707T3-FSLG
5182507T1--ELG	5182507T3--VGA	5182807T1--LG
5182507T1--ESG	5182558T1--GA	5182807T1--EKO
5182507T1--VGA	5182558T1--LG	5182807T1--ELG
5182507T1--XGA	5182558T3--LG	5182807T2--LG
5182507T1--XLG	5182603T2--LGE	5182807T2--EKO
5182507T1-FSLGG	5182606T3--GA	5182807T2--ELG
5182507T1A--LG	5182607T1--GA	5182807T3--LG
5182507T2--GA	5182607T2--LG	5185307T0--OO-1
5182507T2--GAG	5182607T2--LGE	5185307T0--SO-1
5182507T2--HR	5182607T2--EGA	5185307T0--SO-V
5182507T2--LG	5182607T3--GA	5185307T0--SS-V
5182507T2--SC	5182607T3--VGA	5185307T7--FF
5182507T2--EGA	5182658T1--LGE	5185307T7--FF-A
5182507T2--EGAG	5182658T2--LGE	5185307T7--FH
5182507T2--ELG	5182706T1--FULG	5185307T7--FH-A
5182507T2--VGA	5182707T0--GA	5185307T7--FO
5182507T2--XGA	5182707T1--GA	5185307T7--FO-A
5182507T2--FSLG	5182707T1--LG	5185307T7--HH-A
5182507T2B--GA	5182707T1--KFSG	5185307T7--NN-A
5182507T3--F2	5182707T1--FSLG	5185307T7--OO-1
5182507T3--GA	5182707T1--FULG	5185307T7--OO-A
5182507T3--GAG	5182707T2--GA	5185307T7--RR-A
5182507T3--LG	5182707T2--LG	5185307T7--SO-1
5182507T3--SC	5182707T2--FSLG	5185307T7--SO-A

5185307T7-SS-A	5185307T9-OOOAA	5195307A0G-SS-V
5185307T7-SSP1A	5185307T9-SO-1	5195307A7G-FF-1
5185307T7-WW-A	5185307T9-SO-A	5195307A7G-FF-A
5185307T8-FF-A	5185307T9-SS	5195307A7G-FO-1
5185307T8-NN-A	5185307T9-SS-A	5195307A7G-FO-A
5185307T8-OO	5185507T1--GA	5195307A7G-OO-1
5185307T8-OO-1	5187307T1-BB	5195307A7G-OO-A
5185307T8-OO-A	5187307T1-FF	5195307A7G-SO-1
5185307T8-OO-K	5187307T1-SB	5195307A7G-SO-A
5185307T8-OOOAA	5187307T1-SS	5195307A7G-SS-1
5185307T8-RR-A	5187307T2-SS	5195307A7G-SS-A
5185307T8-SO-1	5191307A4G-OO-1	5195307A8G-FF-1
5185307T8-SO-A	5191307A4G-OO-B	5195307A8G-FF-A
5185307T8-SS	5191307A4G-SO-1	5195307A8G-FO-1
5185307T8-SS-1	5191307A4G-SO-B	5195307A8G-FO-A
5185307T8-SS-A	5191307A4G-SS-1	5195307A8G-OO-1
5185307T8-SSCCA	5191307A4G-SS-B	5195307A8G-OO-A
5185307T8-WW-A	5195307A0G-FF-1	5195307A8G-SO-1
5185307T8-ZZ-A	5195307A0G-FF-V	5195307A8G-SO-A
5185307T9-FF-1	5195307A0G-FO-1	5195307A8G-SS-1
5185307T9-FF-A	5195307A0G-FO-V	5195307A8G-SS-A
5185307T9-FH-A	5195307A0G-OO-1	5195307A9G-FF-1
5185307T9-HF-A	5195307A0G-OO-V	5195307A9G-FF-A
5185307T9-NN-A	5195307A0G-SO-1	5195307A9G-FO-1
5185307T9-OO-1	5195307A0G-SO-V	5195307A9G-FO-A
5185307T9-OO-A	5195307A0G-SS-1	5195307A9G-OO-1

5195307A9G-OO-A	5202506A3	5202805A3
5195307A9G-SO-1	5202506A4	5202806A2
5195307A9G-SO-A	5202506F1--BE	5202806A3
5195307A9G-SS-1	5202506F2--BE	5202807D3
5195307A9G-SS-A	5202506F3--BE	5202807D3---G
5197307A1G-OO	5202506F3---BEC	5202B06F3--BE
5197307A1G-OOIT	5202506F4--BE	5202B06F3--BEC
5197307A1G-SO	5202507D3	5212107A3
5197307A1G-SOIT	5202507D4	5212507A1
5197307A1G-SS	5202566D3---BEC	5212507A2
5197307A1G-SSIT	5202605F2--BE	5212507A3
5197307A2G-OO	5202605F3--BE	5212507A3---G
5197307A2G-OOIT	5202606F1--BE	5212707A3
5197307A2G-SO	5202606F2--BE	5212807A3
5197307A2G-SOIT	5202606F3--BE	5212807A3---G
5197307A2G-SS	5202606F3---BEC	5212807A3-E
5197307A2G-SSIT	5202705A2	5231306D4-SS-B
5202106F1--BE	5202705A3	5231306D5-OO-B
5202505A1	5202705F1--BE	5231307H4-OO
5202505A2	5202705F2--BE	5231307H4-OO-B
5202505A3	5202705F3--BE	5231307H4-OOAEB
5202505A4	5202705F3---BEC	5231307H4-ODDPB
5202505F3--BE	5202706F1--BE	5231307H4-OOPTB
5202505F3--BEC	5202706F2--BE	5231307H4-SO-B
5202506A1	5202706F3--BE	5231307H4-SOAEB
5202506A2	5202706F3---BEC	5231307H4-SS-B

5231307H4-SSVPB	5231807H7--B	5232507H1--PH
5231307H5-NO-B	5231807H8--B	5232507H1-EIVG
5231307H5-NO-V	5231807H9--B	5232507H1-EP
5231307H5-OO	5232106D1--P	5232507H2--BH
5231307H5-OO-B	5232107H1--IVG	5232507H2--HB
5231307H5-OO-D	5232107H1--P	5232507H2--IVG
5231307H5-OO-V	5232107H2--P	5232507H2--P
5231307H5-ODDPB	5232158H1--P	5232507H2-EIVG
5231307H5-OOPTB	5232506D0--P	5232507H2-EP
5231358G0-SS-V	5232506D1--IVG	5232507H3--IVG
5231358H0-SO-1	5232506D1--P	5232507H3--P
5231358H5-OO-D	5232506D1--P-H	5232507H3--PH
5231506D7--B	5232506D2--IVG	5232507H3-EIVG
5231506D8--B	5232506D2--P	5232507H3-EP
5231506D9	5232506D3--IVG	5232558H0--P
5231506D9--B	5232506D3--P	5232558H1--P
5231507H7	5232506D4--IVG	5232558H2--P
5231507H7--B	5232506D4--P	5232558H3--P
5231507H7-MRUM	5232507G1-PH-2O	5232606D2--P
5231507H8	5232507G2-PH-2O	5232607G1-PH-2O
5231507H8--B	5232507G3-PH-2O	5232607H1--P
5231507H8-V-B	5232507H0--P	5232607H2--P
5231507H9	5232507H0-VP	5232607H3--P
5231507H9--B	5232507H0-V-P	5232706D1--P
5231558H8--B	5232507H1--IVG	5232706D2--P
5231558H9--B	5232507H1--P	5232706D3--P

5232707G1--PH	5235306D0-SS-V	5235307H0-SO-1
5232707G2-PH-2O	5235306D7-FF-A	5235307H0-SO-V
5232707H1--P	5235306D7-OO-1	5235307H0-SS-V
5232707H2--P	5235306D7-OO-A	5235307H4-ODDPB
5232707H3--P	5235306D7-ODPA	5235307H5-ODDPB
5232758H1--P	5235306D7-ODDPB	5235307H7-FF-A
5232758H2--P	5235306D7-SO-1	5235307H7-FH-A
5232758H3-EP	5235306D7-SO-A	5235307H7-FO-A
5232806D1--IVG	5235306D7-SS-A	5235307H7-HH-A
5232806D1--P	5235306D7-WW-A	5235307H7-NN-A
5232806D2--P	5235306D8-FF-A	5235307H7-OO
5232806D3--P	5235306D8-OO-1	5235307H7-OO-1
5232807H1--IVG	5235306D8-OO-A	5235307H7-OO-A
5232807H1--P	5235306D8-SO-1	5235307H7-ODPA
5232807H2--P	5235306D8-SO-A	5235307H7-SO
5232807H3--IVG	5235306D8-SS-A	5235307H7-SO-1
5232807H3--P	5235306D9-NN-A	5235307H7-SO-A
5232858H1--P	5235306D9-OO-1	5235307H7-SO-K
5232858H2--P	5235306D9-OO-A	5235307H7-SODPA
5232858H3--P	5235306D9-SO-1	5235307H7-SS-A
5232858H4--P	5235306D9-SO-A	5235307H7-SSDPA
5235306D0-FF-V	5235306D9-SS-A	5235307H7-WW-A
5235306D0-OO-1	5235307H0-FF-V	5235307H8-FF-A
5235306D0-OO-V	5235307H0-HH-V	5235307H8-NO-A
5235306D0-SO-1	5235307H0-OO-1	5235307H8-OO
5235306D0-SO-V	5235307H0-OO-V	5235307H8-OO-1

5235307H8-OO-A	5235358H0-OO-1	5237306D2-ODX
5235307H8-RO-A	5235358H0-SO-1	5237306D2-OOIT
5235307H8-RR-A	5235358H0-SO-V	5237306D2-SB
5235307H8-SO	5235358H0-SS-V	5237306D2-SO
5235307H8-SO-1	5235358H7-OO-1	5237306D2-SS
5235307H8-SO-A	5235358H7-OO-A	5237306D3-BB
5235307H8-SO-K	5235358H7-SO-1	5237306D3-FF
5235307H8-SODPA	5235358H7-SO-A	5237306D3-OO
5235307H8-SS	5235358H7-SS-A	5237306D3-SB
5235307H8-SS-A	5235358H8-OO-1	5237306D3-SO
5235307H8-SSDPA	5235358H8-OO-A	5237306D3-SS
5235307H8-WW-A	5235358H8-SO-1	5237307G3H-OO2O
5235307H8-ZO-A	5235358H8-SS-A	5237307G3H-SS2O
5235307H8-ZZ-A	5235358H9-OO-1	5237307H1-BB
5235307H9-FF	5235358H9-OO-A	5237307H1-FF
5235307H9-FF-A	5235358H9-SO-1	5237307H1-OO
5235307H9-NN-A	5235358H9-SS-A	5237307H1-OOIT
5235307H9-OO-1	5237306D1-BB	5237307H1-SB
5235307H9-OO-A	5237306D1-FF	5237307H1-SO
5235307H9-OOCEA	5237306D1-OO	5237307H1-SOIT
5235307H9-RR-A	5237306D1-SB	5237307H1-SS
5235307H9-SO-1	5237306D1-SO	5237307H1-SSDP
5235307H9-SO-A	5237306D1-SS	5237307H1-SSIT
5235307H9-SODPA	5237306D2-BB	5237307H2-BB
5235307H9-SS-A	5237306D2-FF	5237307H2-FF
5235307H9-SSDPA	5237306D2-OO	5237307H2-OO

5237307H2-OOIT	5237358H3-SO	5302505A3--SC
5237307H2-SB	5237358H3-SS	5302505A4
5237307H2-SO	5238306D1-OO	5302505A4--EA
5237307H2-SOIT	5238306D2-OO	5302505A4--KI
5237307H2-SS	5238306D3-OO	5302506A1
5237307H2-SSDP	5238307H1-OO	5302506A2
5237307H2-SSIT	5238307H1-SO	5302506A2--G
5237307H3-BB	5238307H1-SS	5302506A3
5237307H3-FF	5238307H1-SY	5302506A3--G
5237307H3-OO	5238307H1-YY	5302506A3--EA
5237307H3-ODDX	5238307H2-OO	5302506A3--KI
5237307H3-OOIT	5238307H2-SO	5302506A3--SC
5237307H3-SB	5238307H2-SS	5302506A4
5237307H3-SO	5238307H2-SY	5302506A4--G
5237307H3-SOIT	5238307H2-YY	5302506A4--EA
5237307H3-SS	5238307H3-OO	5302506A4--KI
5237307H3-SSIT	5238307H3-SO	5302706A3
5237358H1-FF	5238307H3-SS	5302805A3
5237358H1-OO	5238307H3-SY	5302805A3--G
5237358H1-SO	5238307H3-YY	5302805A3--EA
5237358H1-SS	5302105A1--A-K	5302806A1
5237358H2-FF	5302505A1	5302806A2
5237358H2-OO	5302505A2	5302806A3
5237358H2-SO	5302505A3	5302806A3--G
5237358H2-SS	5302505A3--EA	5302806A3--EA
5237358H3-OO	5302505A3--KI	5302806A4

5302806A4--EA	5401307A5-OO-B	5412507A4--EB
5312505A3	5402807A1	5412507A4--SC
5312505A3-E	5412107A1	5412707A1--EB
5312505A4	5412107A1--A-K	5412707A2
5312524A1--EB	5412107A1--FW	5412707A3
5312524A2	5412107A1--GF	5412807A1
5312524A2--EB	5412107A2	5412807A2
5312524A3	5412507A1	5412807A2--EA
5312524A3--G	5412507A1--EB	5412807A2--EB
5312524A3--EB	5412507A1--XEB	5412807A3
5312524A4	5412507A1-E	5412807A3--G
5312524A4--EB	5412507A1-XEB	5412807A3--EA
5312724A3	5412507A2	5412807A3--EB
5312805A3	5412507A2--EA	5412807A3-E
5312824A2	5412507A2--EB	5412807A3-EK
5312824A3	5412507A2-E	5412807A4
5312824A4	5412507A3	5415307A7--BRK
5382507H2	5412507A3--G	5422105A1--A-K
5382507H2--G	5412507A3--EA	5422105A1--GF
5382507H3	5412507A3--EB	5422106A1--A-K
5382507H3--G	5412507A3--SC	5422106A1--GF
5382507H4	5412507A3-E	5422505A1
5382507H4--G	5412507A3-V	5422505A1--EB
5382558K3	5412507A4	5422505A2
5382807H3	5412507A4--G	5422505A2--EA
5382858K3	5412507A4--EA	5422505A3

5422505A3--G	5422705A3--EB	5441307G4-SS-B
5422505A3--EA	5422706A3	5441307G4-SSVPB
5422505A3--EB	5422706A3--EB	5441307G4G-OO-1
5422505A3--FW	5422805A1	5441307G4G-OO-B
5422505A3--JK	5422805A2	5441307G4G-SO-1
5422505A3--KI	5422805A3	5441307G4G-SO-B
5422505A3--SC	5422805A3--G	5441307G4G-SS-1
5422505A3-E	5422805A3--EA	5441307G4G-SS-B
5422505A4	5422806A1	5441307G4G-SSVPB
5422505A4--EB	5422806A2	5441307H4-NO-B
5422505A4--KI	5422806A2--EA	5441307H4-NOIN
5422506A1	5422806A3	5441307H4-OO
5422506A1--EB	5422806A3--G	5441307H4-OO-B
5422506A2	5422806A3--EA	5441307H4-SO-B
5422506A2--EA	5422806A4	5441307H4-SS-B
5422506A2--EB	5441306A5-OOEIK	5441307H4-SSVP
5422506A2-E	5441306G4-OO-B	5441307H4-SSVPB
5422506A3	5441306G5-OO	5441307H4G-OO-1
5422506A3--G	5441306G5-OO-1	5441307H4G-OO-B
5422506A3--EA	5441306G5-OO-B	5441307H4G-OOOXQ
5422506A3--EB	5441307A4G-OO-B	5441307H4G-OOSD
5422506A3--SC	5441307A4G-SO-B	5441307H4G-OOTCB
5422506A4	5441307A4G-SS-B	5441307H4G-SO-1
5422506A4--EB	5441307G4-NO-B	5441307H4G-SO-B
5422506A4--SC	5441307G4-OO-B	5441307H4G-SOTC1
5422705A3	5441307G4-SO-B	5441307H4G-SOTCB

5441307H4G-SS-1	5441307I4G-SSVPB	5441507H7--B
5441307H4G-SS-B	5441307IA-OO-B	5441507H7--IOB
5441307H4G-SSVPB	5441358K4-OO-B	5441507H7-E-B
5441307H5-OO	5441358K4-SO-B	5441507H8
5441307H5-OO-1	5441358K4-SS-B	5441507H8--B
5441307H5-OO-B	5441358K4G-OO-1	5441507H9
5441307H5-OOPIB	5441358K4G-OO-B	5441507H9--B
5441307H5G-OO-1	5441358K4G-SO-1	5441507H9-E-B
5441307H5G-OO-B	5441358K4G-SO-B	5441507H9-25-B
5441307H5G-OOSD	5441358K5-OO	5441558K7--B
5441307H5G-OOSD1	5441358K5-OO-B	5441558K8--B
5441307H5G-OOSDK	5441358K8G-SS-B	5441558K9
5441307H5G-OOTCB	5441358K9G-SS-A	5441558K9--B
5441307H9G-OO-A	5441358M4-OO-B	5441558M8-25-B
5441307I4-NO-B	5441358M4-SO-B	5441558M9-25-B
5441307I4-OO-B	5441358M4-SS-B	5441807H7-E-B
5441307I4-SO-B	5441358M4G-OO-1	5441807H8--B
5441307I4-SS-B	5441358M4G-OO-B	5441807H9--B
5441307I4-SSVP	5441358M4G-SO-1	5442106G1
5441307I4-SSVPB	5441358M4G-SO-B	5442107H1
5441307I4G-OO-1	5441358M4G-SS-1	5442107H2
5441307I4G-OO-B	5441358M4G-SS-B	5442107H3
5441307I4G-SO-1	5441506G7--B	5442158K1--GF
5441307I4G-SO-B	5441506G8--B	5442506G0
5441307I4G-SS-1	5441506G9--B	5442506G1
5441307I4G-SS-B	5441507H7	5442506G1--IVG

5442506G2	5442507H1--HB	5442558K2
5442506G2--IVG	5442507H1--IVG	5442558K2-E
5442506G2-E	5442507H1--SI	5442558K3
5442506G3	5442507H1-E	5442558K3--IVG
5442506G3--IVG	5442507H2	5442558K3-E
5442506G4	5442507H2--G	5442558M0
5442507A1--HF	5442507H2--HB	5442558M1
5442507A1--HFG	5442507H2--IVG	5442558M1--EB
5442507A2--HF	5442507H2-E	5442558M2
5442507A2--HFG	5442507H3	5442558M2--EB
5442507A3--HF	5442507H3--G	5442558M3
5442507A3--HFG	5442507H3--AO	5442558M3--EB
5442507G0	5442507H3--IVG	5442558M4
5442507G1	5442507H3-E	5442606G2-EAB
5442507G1--G	5442507H3-EIVG	5442607G2--ABG
5442507G2	5442507H4	5442607H1
5442507G2--G	5442507I1	5442706G1
5442507G2-E	5442507I1-E	5442706G2
5442507G3	5442507I2	5442706G3
5442507G3--G	5442507I2-E	5442706G4
5442507H0	5442507I3	5442707H1
5442507H0--G	5442507I3-E	5442707H1-VSI
5442507H0--HB	5442558K0	5442707H2
5442507H0-E	5442558K1	5442707H2--IVG
5442507H1	5442558K1--IVG	5442707H3
5442507H1--G	5442558K1-E	5442758K1

5442758K2	5445306G7-SS-A	5445306G9G-SO-1
5442806G3--IVG	5445306G7G-OO-1	5445306G9G-SO-A
5442807H1	5445306G7G-OO-A	5445306G9G-SS-1
5442807H1--IVG	5445306G7G-SO-1	5445306G9G-SS-A
5442807H2	5445306G7G-SO-A	5445307A0G-OO-V
5442807H3	5445306G7G-SS-1	5445307A0G-SO-V
5442858K1	5445306G7G-SS-A	5445307A0G-SS-V
5442858K2	5445306G8-FF-A	5445307A7G-OO-A
5442858K3	5445306G8-HH-A	5445307A7G-SO-A
5445306A7-HHEIK	5445306G8-OO-A	5445307A7G-SS-A
5445306G0-FF-V	5445306G8-SO-A	5445307A8G-OO-A
5445306G0-OO-V	5445306G8-SS-1	5445307A8G-SO-A
5445306G0-SO-V	5445306G8-SS-A	5445307A8G-SS-A
5445306G0-SS-V	5445306G8G-OO-1	5445307A9G-OO-A
5445306G0G-OO-1	5445306G8G-OO-A	5445307A9G-SO-A
5445306G0G-OO-V	5445306G8G-SO-1	5445307A9G-SS-A
5445306G0G-SO-1	5445306G8G-SO-A	5445307G0-FF-V
5445306G0G-SO-V	5445306G8G-SS-1	5445307G0-FO-V
5445306G0G-SS-1	5445306G8G-SS-A	5445307G0-OO-V
5445306G0G-SS-V	5445306G9-FF-A	5445307G0-SO-V
5445306G7-FF-A	5445306G9-OO-A	5445307G0-SS-V
5445306G7-FH-A	5445306G9-SO	5445307G0G-OO-1
5445306G7-FO-A	5445306G9-SO-A	5445307G0G-OO-V
5445306G7-HH-A	5445306G9-SS-A	5445307G0G-SO-1
5445306G7-OO-A	5445306G9G-OO-1	5445307G0G-SO-V
5445306G7-SO-A	5445306G9G-OO-A	5445307G0G-SS-1

5445307G0G-SS-V	5445307G9-SO-A	5445307H7-SO-A
5445307G7-FF-A	5445307G9-SS-A	5445307H7-SS-A
5445307G7-FO-A	5445307G9G-OO-1	5445307H7-EOO-A
5445307G7-OO-A	5445307G9G-OO-A	5445307H7-ESS-A
5445307G7-SO-A	5445307G9G-SO-1	5445307H7G-FF-1
5445307G7-SS-A	5445307G9G-SO-A	5445307H7G-FF-A
5445307G7G-OO-1	5445307G9G-SS-1	5445307H7G-FH-A
5445307G7G-OO-A	5445307G9G-SS-A	5445307H7G-FO-A
5445307G7G-SO-1	5445307H0-FF-V	5445307H7G-HH-1
5445307G7G-SO-A	5445307H0-OO-V	5445307H7G-HH-A
5445307G7G-SS-1	5445307H0-SF-V	5445307H7G-OO-1
5445307G7G-SS-A	5445307H0-SO-V	5445307H7G-OO-A
5445307G8-FF-A	5445307H0-SS-V	5445307H7G-SO-1
5445307G8-FO-A	5445307H0G-OO	5445307H7G-SO-A
5445307G8-OO-A	5445307H0G-OO-1	5445307H7G-SS-1
5445307G8-SO-A	5445307H0G-OO-V	5445307H7G-SS-A
5445307G8-SS-A	5445307H0G-SO-V	5445307H8-FF-A
5445307G8G-OO-1	5445307H0G-SO-1	5445307H8-OO-A
5445307G8G-OO-A	5445307H0G-SO-V	5445307H8-SF-A
5445307G8G-SO-1	5445307H0G-SS-1	5445307H8-SO-A
5445307G8G-SO-A	5445307H0G-SS-V	5445307H8-SS-A
5445307G8G-SS-1	5445307H7-FF-A	5445307H8-WW-A
5445307G8G-SS-A	5445307H7-FH-A	5445307H8-EOO-A
5445307G9-FF-A	5445307H7-FO-A	5445307H8-ESO-A
5445307G9-FO-A	5445307H7-HH-A	5445307H8-ESS-A
5445307G9-OO-A	5445307H7-OO-A	5445307H8G-OO-1

5445307H8G-OO-A	5445307I0G-OO-1	5445307I8G-FF-1
5445307H8G-SO-1	5445307I0G-OO-V	5445307I8G-FF-A
5445307H8G-SO-A	5445307I0G-SO-1	5445307I8G-FO-1
5445307H8G-SS-1	5445307I0G-SO-V	5445307I8G-FO-A
5445307H8G-SS-A	5445307I0G-SS-1	5445307I8G-OO-1
5445307H9-FF-A	5445307I0G-SS-V	5445307I8G-OO-A
5445307H9-OO-A	5445307I7-FF-A	5445307I8G-SO-1
5445307H9-SO-A	5445307I7-FO-A	5445307I8G-SO-A
5445307H9-SS-A	5445307I7-OO-A	5445307I8G-SS-1
5445307H9G-OO-1	5445307I7-SO-A	5445307I8G-SS-A
5445307H9G-OO-A	5445307I7-SS-A	5445307I9-FF-A
5445307H9G-SO-1	5445307I7G-FF-1	5445307I9-FO-A
5445307H9G-SO-A	5445307I7G-FF-A	5445307I9-OO-A
5445307H9G-SS-1	5445307I7G-FO-1	5445307I9-SO-A
5445307H9G-SS-A	5445307I7G-FO-A	5445307I9-SS-A
5445307H0G-SO	5445307I7G-OO-1	5445307I9G-FF-1
5445307H0G-SS	5445307I7G-OO-A	5445307I9G-FF-A
5445307I0-FF-V	5445307I7G-SO-1	5445307I9G-FO-1
5445307I0-FO-V	5445307I7G-SO-A	5445307I9G-FO-A
5445307I0-OO-V	5445307I7G-SS-1	5445307I9G-OO-1
5445307I0-SO-V	5445307I7G-SS-A	5445307I9G-OO-A
5445307I0-SS-V	5445307I8-FF-A	5445307I9G-SO-1
5445307I0G-FF-1	5445307I8-FO-A	5445307I9G-SO-A
5445307I0G-FF-V	5445307I8-OO-A	5445307I9G-SS-1
5445307I0G-FO-1	5445307I8-SO-A	5445307I9G-SS-A
5445307I0G-FO-V	5445307I8-SS-A	5445358K0-FF-V

5445358K0-OO-V	5445358K7G-SO-1	5445358KOG-OO
5445358K0-SO-V	5445358K7G-SO-A	5445358M0-FF-V
5445358K0-SS-V	5445358K7G-SS-1	5445358M0-FO-V
5445358KOG-OO-1	5445358K7G-SS-A	5445358M0-OO-V
5445358KOG-OO-V	5445358K8-FF-A	5445358M0-SO-V
5445358KOG-SO	5445358K8-OO-A	5445358M0-SS-V
5445358KOG-SO-1	5445358K8-SO-A	5445358M0-OO-V
5445358KOG-SO-V	5445358K8-SS-A	5445358M0-SO-V
5445358KOG-SS	5445358K8G-OO-1	5445358M0-SS-V
5445358KOG-SS-1	5445358K8G-OO-A	5445358M0G-FF-1
5445358KOG-SS-V	5445358K8G-SO-1	5445358M0G-FF-V
5445358K7-FF-A	5445358K8G-SO-A	5445358M0G-FO-1
5445358K7-FH-A	5445358K8G-SS-1	5445358M0G-FO-V
5445358K7-FO-A	5445358K8G-SS-A	5445358M0G-OO-1
5445358K7-HH-1	5445358K9-FF	5445358M0G-OO-V
5445358K7-HH-A	5445358K9-FF-A	5445358M0G-SO-1
5445358K7-OO-A	5445358K9-OO-1	5445358M0G-SO-V
5445358K7-SO-A	5445358K9-OO-A	5445358M0G-SS-1
5445358K7-SS-A	5445358K9-SO-A	5445358M0G-SS-V
5445358K7G-FF-1	5445358K9-SS-A	5445358M7-FF-A
5445358K7G-FF-A	5445358K9G-OO-1	5445358M7-FO-A
5445358K7G-FH-A	5445358K9G-OO-A	5445358M7-OO-A
5445358K7G-FO-A	5445358K9G-SO-1	5445358M7-SO-A
5445358K7G-HH-A	5445358K9G-SO-A	5445358M7-SS-A
5445358K7G-OO-1	5445358K9G-SS-1	5445358M7-FF-A
5445358K7G-OO-A	5445358K9G-SS-A	5445358M7-OO-A

5445358M7-SO-A	5445358M8G-SO-A	5447306G1G-OO
5445358M7-SS-A	5445358M8G-SS-1	5447306G1G-SO
5445358M7G-FF-1	5445358M8G-SS-A	5447306G1G-SS
5445358M7G-FF-A	5445358M9-FF-A	5447306G2-FF
5445358M7G-FO-1	5445358M9-FO-A	5447306G2-OO
5445358M7G-OO-1	5445358M9-OO-A	5447306G2-SO
5445358M7G-OO-A	5445358M9-SO-A	5447306G2-SS
5445358M7G-SO-1	5445358M9-SS-A	5447306G2G-OO
5445358M7G-SO-A	5445358M9-OO-A	5447306G2G-SO
5445358M7G-SS-1	5445358M9-SO-A	5447306G2G-SS
5445358M7G-SS-A	5445358M9-SS-A	5447306G3-FF
5445358M8-FF-A	5445358M9G-FF-1	5447306G3-OO
5445358M8-FO-A	5445358M9G-FF-A	5447306G3-SO
5445358M8-OO-A	5445358M9G-FO-1	5447306G3-SS
5445358M8-SO-A	5445358M9G-FO-A	5447306G3-SSAC
5445358M8-SS-A	5445358M9G-OO-1	5447306G3G-OO
5445358M8-OO-A	5445358M9G-OO-A	5447306G3G-SO
5445358M8-SO-A	5445358M9G-SO-1	5447306G3G-SS
5445358M8-SS-A	5445358M9G-SO-A	5447307G1-FF
5445358M8G-FF-1	5445358M9G-SS-1	5447307G1-OO
5445358M8G-FF-A	5445358M9G-SS-A	5447307G1-OOIT
5445358M8G-FO-1	5447306G1-FF	5447307G1-SO
5445358M8G-FO-A	5447306G1-OO	5447307G1-SOIT
5445358M8G-OO-1	5447306G1-SO	5447307G1-SS
5445358M8G-OO-A	5447306G1-SS	5447307G1-SSAC
5445358M8G-SO-1	5447306G1G-SSAC	5447307G1-SSIT

5447307G1G-OO	5447307G3-SSIT	5447307H2-OOIT
5447307G1G-OOIT	5447307G3G-OO	5447307H2-SO
5447307G1G-SO	5447307G3G-OOIT	5447307H2-SOHB
5447307G1G-SOIT	5447307G3G-SO	5447307H2-SOIT
5447307G1G-SS	5447307G3G-SOIT	5447307H2-SS
5447307G1G-SSIT	5447307G3G-SS	5447307H2-SSAC
5447307G2-FF	5447307G3G-SSIT	5447307H2-SSHB
5447307G2-OO	5447307H1-FF	5447307H2-SSIT
5447307G2-OOIT	5447307H1-OO	5447307H2G-OO
5447307G2-SO	5447307H1-OOAC	5447307H2G-SO
5447307G2-SOIT	5447307H1-OOIT	5447307H2G-SOIT
5447307G2-SS	5447307H1-SF	5447307H2G-SS
5447307G2-SSAC	5447307H1-SO	5447307H2G-SSIT
5447307G2-SSIT	5447307H1-SOIT	5447307H3-FF
5447307G2G-OO	5447307H1-SS	5447307H3-OO
5447307G2G-OOIT	5447307H1-SSAC	5447307H3-OOAC
5447307G2G-SO	5447307H1-SSIT	5447307H3-OOIT
5447307G2G-SOIT	5447307H1G-FO	5447307H3-SO
5447307G2G-SS	5447307H1G-OO	5447307H3-SOIT
5447307G2G-SSIT	5447307H1G-SO	5447307H3-SS
5447307G3-OO	5447307H1G-SOIT	5447307H3-SSAC
5447307G3-OOIT	5447307H1G-SS	5447307H3-SSIT
5447307G3-SO	5447307H1G-SSIT	5447307H3G-OO
5447307G3-SOIT	5447307H2-FF	5447307H3G-SO
5447307G3-SS	5447307H2-OO	5447307H3G-SOIT
5447307G3-SSAC	5447307H2-OOAC	5447307H3G-SS

5447307H3G-SSIT	5447307I2G-SO	5447358K1G-FO
5447307H3GMSS	5447307I2G-SOIT	5447358K1G-OO
5447307I1-OO	5447307I2G-SS	5447358K1G-SO
5447307I1-OOAC	5447307I2G-SSIT	5447358K1G-SOIT
5447307I1-OOIT	5447307I3-OO	5447358K1G-SS
5447307I1-SO	5447307I3-OOAC	5447358K1G-SSIT
5447307I1-SOIT	5447307I3-OOIT	5447358K2-FF
5447307I1-SS	5447307I3-SO	5447358K2-OO
5447307I1-SSAC	5447307I3-SOIT	5447358K2-OOIT
5447307I1-SSIT	5447307I3-SS	5447358K2-SO
5447307I1G-OO	5447307I3-SSAC	5447358K2-SOIT
5447307I1G-OOIT	5447307I3-SSIT	5447358K2-SS
5447307I1G-SO	5447307I3G-OO	5447358K2-SSAC
5447307I1G-SOIT	5447307I3G-OOIT	5447358K2-SSIT
5447307I1G-SS	5447307I3G-SO	5447358K2G-FO
5447307I1G-SSIT	5447307I3G-SOIT	5447358K2G-SO
5447307I2-OO	5447307I3G-SS	5447358K2G-SOIT
5447307I2-OOAC	5447307I3G-SSIT	5447358K2G-SS
5447307I2-OOIT	5447358K1-FF	5447358K2G-SSIT
5447307I2-SO	5447358K1-OO	5447358K3-FF
5447307I2-SOIT	5447358K1-OOIT	5447358K3-OO
5447307I2-SS	5447358K1-SO	5447358K3-OOIT
5447307I2-SSAC	5447358K1-SOIT	5447358K3-SO
5447307I2-SSIT	5447358K1-SS	5447358K3-SOIT
5447307I2G-OO	5447358K1-SSAC	5447358K3-SS
5447307I2G-OOIT	5447358K1-SSIT	5447358K3-SSIT

5447358K3G-SO	5447358M3G-SO	5448307G1G-SY
5447358K3G-SS	5447358M3G-SOIT	5448307G1G-YY
5447358K3G-SSIT	5447358M3G-SS	5448307G2G-OO
5447358M1-OO	5447358M3G-SSIT	5448307G2G-SO
5447358M1-SO	5448306G1G-OO	5448307G2G-SS
5447358M1-SS	5448306G1G-SO	5448307G2G-SY
5447358M1G-OO	5448306G1G-SS	5448307G2G-YY
5447358M1G-OOIT	5448306G2G-OO	5448307G3G-OO
5447358M1G-SO	5448306G2G-SO	5448307G3G-SO
5447358M1G-SOIT	5448306G2G-SS	5448307G3G-SS
5447358M1G-SS	5448306G3G-OO	5448307G3G-SY
5447358M1G-SSIT	5448306G3G-SO	5448307G3G-YY
5447358M2-OO	5448306G3G-SS	5448307H1G-OO
5447358M2-SO	5448307A1G-OO	5448307H1G-SO
5447358M2-SS	5448307A1G-SO	5448307H1G-SS
5447358M2G-OO	5448307A1G-SS	5448307H1G-SSAC
5447358M2G-OOIT	5448307A2G-OO	5448307H1G-SSPZ
5447358M2G-SO	5448307A2G-SO	5448307H1G-SY
5447358M2G-SOIT	5448307A2G-SS	5448307H1G-YY
5447358M2G-SS	5448307A3G-OO	5448307H2G-OO
5447358M2G-SSIT	5448307A3G-SO	5448307H2G-SO
5447358M3-OO	5448307A3G-SS	5448307H2G-SS
5447358M3-SO	5448307G1G-OO	5448307H2G-SSPZ
5447358M3-SS	5448307G1G-OOAC	5448307H2G-SY
5447358M3G-OO	5448307G1G-SO	5448307H2G-YY
5447358M3G-OOIT	5448307G1G-SS	5448307H3G-OO

5448307H3G-SO	5448358K3G-SS	5462507G3
5448307H3G-SS	5448358K3G-SY	5462558M1
5448307H3G-SSPZ	5448358K3G-YY	5462558M2
5448307H3G-SY	5448358M1G-OO	5462558M3
5448307H3G-YY	5448358M1G-SO	5465307G0G-FF-1
5448307I1G-OO	5448358M1G-SS	5465307G0G-FF-B
5448307I1G-SO	5448358M1G-SY	5465307G0G-FF-V
5448307I1G-SS	5448358M1G-YY	5465307G0G-FO-1
5448307I2G-OO	5448358M2G-OO	5465307G0G-FO-B
5448307I2G-SO	5448358M2G-SO	5465307G0G-FO-V
5448307I2G-SS	5448358M2G-SS	5465307G0G-OO-1
5448307I3G-OO	5448358M2G-SY	5465307G0G-OO-B
5448307I3G-SO	5448358M2G-YY	5465307G0G-OO-V
5448307I3G-SS	5448358M3G-OO	5465307G0G-SO-1
5448358K1G-OO	5448358M3G-SO	5465307G0G-SO-B
5448358K1G-SO	5448358M3G-SS	5465307G0G-SO-V
5448358K1G-SS	5448358M3G-SY	5465307G0G-SS-1
5448358K1G-SY	5448358M3G-YY	5465307G0G-SS-B
5448358K1G-YY	5461307G4-OO-B	5465307G0G-SS-V
5448358K2G-OO	5461307G4-SO-B	5465307G7G-OO-A
5448358K2G-SO	5461307G4-SS-B	5465307G8G-FF-1
5448358K2G-SS	5461358M4-OO-B	5465307G8G-FF-A
5448358K2G-SY	5461358M4-SO-B	5465307G8G-FO-1
5448358K2G-YY	5461358M4-SS-B	5465307G8G-FO-A
5448358K3G-OO	5462507G1	5465307G8G-OO-1
5448358K3G-SO	5462507G2	5465307G8G-OO-A

5465307G8G-SO-1	5465358M0G-SS-1	5467307G1G-SO
5465307G8G-SO-A	5465358M0G-SS-B	5467307G1G-SOIT
5465307G8G-SS-1	5465358M0G-SS-V	5467307G1G-SS
5465307G8G-SS-A	5465358M7G-SO-A	5467307G1G-SSIT
5465307G9G-FF-1	5465358M8G-FF-1	5467307G2G-OO
5465307G9G-FF-A	5465358M8G-FF-A	5467307G2G-OOIT
5465307G9G-FO-1	5465358M8G-FO-1	5467307G2G-SO
5465307G9G-FO-A	5465358M8G-FO-A	5467307G2G-SOIT
5465307G9G-OO-1	5465358M8G-OO-1	5467307G2G-SS
5465307G9G-OO-A	5465358M8G-OO-A	5467307G2G-SSIT
5465307G9G-SO-1	5465358M8G-SO-1	5467307G3G-OO
5465307G9G-SO-A	5465358M8G-SO-A	5467307G3G-OOIT
5465307G9G-SS-1	5465358M8G-SS-1	5467307G3G-SO
5465307G9G-SS-A	5465358M8G-SS-A	5467307G3G-SOIT
5465358M0G-FF-1	5465358M9G-FF-1	5467307G3G-SS
5465358M0G-FF-B	5465358M9G-FF-A	5467307G3G-SSIT
5465358M0G-FF-V	5465358M9G-FO-1	5467358M1G-OO
5465358M0G-FO-1	5465358M9G-FO-A	5467358M1G-SO
5465358M0G-FO-B	5465358M9G-OO-1	5467358M1G-SS
5465358M0G-FO-V	5465358M9G-OO-A	5467358M2G-OO
5465358M0G-OO-1	5465358M9G-SO-1	5467358M2G-SO
5465358M0G-OO-B	5465358M9G-SO-A	5467358M2G-SS
5465358M0G-OO-V	5465358M9G-SS-1	5467358M3G-OO
5465358M0G-SO-1	5465358M9G-SS-A	5467358M3G-SO
5465358M0G-SO-B	5467307G1G-OO	5467358M3G-SS
5465358M0G-SO-V	5467307G1G-OOIT	5468307G1

5468307G1G-OO	5468358M3G-SO	5472507F4-E
5468307G1G-SO	5468358M3G-SS	5472558G1
5468307G1G-SS	5468358M3G-SY	5472558G2
5468307G1G-SY	5468358M3G-YY	5472558G3
5468307G1G-YY	5471307F4-OO-B	5475307F0-OO-V
5468307G2G-OO	5471307F4-SO-B	5475307F0-SO-V
5468307G2G-SO	5471307F4-SS-B	5475307F0-SS-V
5468307G2G-SS	5471307F4G-OO-B	5475307F0G-OO-1
5468307G2G-SY	5471307F4G-SS-1	5475307F0G-OO-V
5468307G2G-YY	5471307F4G-SS-B	5475307F0G-SO-1
5468307G3G-OO	5471358G4-OO-B	5475307F0G-SO-V
5468307G3G-SO	5471358G4-SO-B	5475307F0G-SS-1
5468307G3G-SS	5471358G4-SS	5475307F0G-SS-V
5468307G3G-SY	5471358G4-SS-B	5475307F7-OO-A
5468307G3G-YY	5471358G4G-OO-B	5475307F7-SO-A
5468358M1G-OO	5471358G4G-SS-1	5475307F7-SS-A
5468358M1G-SO	5471358G4G-SS-B	5475307F7G-OO-1
5468358M1G-SS	5471558G0-25-V	5475307F7G-OO-A
5468358M1G-SY	5471558G9-25-B	5475307F7G-SO-1
5468358M1G-YY	5472507F1	5475307F7G-SO-A
5468358M2G-OO	5472507F1--G	5475307F7G-SS-1
5468358M2G-SO	5472507F2	5475307F7G-SS-A
5468358M2G-SS	5472507F2--G	5475307F8-OO-A
5468358M2G-SY	5472507F2-E	5475307F8-SO-A
5468358M2G-YY	5472507F3	5475307F8-SS-A
5468358M3G-OO	5472507F3-E	5475307F8G-OO-1

5475307F8G-OO-A	5475358G7G-OO-1	5475358G9G-SS-A
5475307F8G-SO-1	5475358G7G-OO-A	5477307F1-OO
5475307F8G-SO-A	5475358G7G-SO-1	5477307F1-SO
5475307F8G-SS-1	5475358G7G-SO-A	5477307F1-SS
5475307F8G-SS-A	5475358G7G-SS-1	5477307F1-SSAC
5475307F9-OO-A	5475358G7G-SS-A	5477307F1G-OO
5475307F9-SO-A	5475358G8-FF-A	5477307F1G-OOIT
5475307F9-SS-A	5475358G8-OO-A	5477307F1G-SO
5475307F9G-OO-1	5475358G8-SO-A	5477307F1G-SOIT
5475307F9G-OO-A	5475358G8-SS-A	5477307F1G-SS
5475307F9G-SO-1	5475358G8G-OO-1	5477307F1G-SSIT
5475307F9G-SO-A	5475358G8G-OO-A	5477307F2-OO
5475307F9G-SS-1	5475358G8G-SO-1	5477307F2-SO
5475307F9G-SS-A	5475358G8G-SO-A	5477307F2-SS
5475358G0-OO-V	5475358G8G-SS-1	5477307F2-SSAC
5475358G0-SO-V	5475358G8G-SS-A	5477307F2G-OO
5475358G0-SS-V	5475358G9-OO-1	5477307F2G-OOIT
5475358G0G-OO-1	5475358G9-OO-A	5477307F2G-SO
5475358G0G-OO-V	5475358G9-SO-A	5477307F2G-SOIT
5475358G0G-SO-1	5475358G9-SS-1	5477307F2G-SS
5475358G0G-SO-V	5475358G9-SS-A	5477307F2G-SSIT
5475358G0G-SS-1	5475358G9G-OO-1	5477307F3-OO
5475358G0G-SS-V	5475358G9G-OO-A	5477307F3-SO
5475358G7-OO-A	5475358G9G-SO-1	5477307F3-SS
5475358G7-SO-A	5475358G9G-SO-A	5477307F3-SSAC
5475358G7-SS-A	5475358G9G-SS-1	5477307F3G-OO

5477307F3G-OOIT	5477358G3G-OO	5478358G2G-OO
5477307F3G-SO	5477358G3G-OOIT	5478358G2G-SO
5477307F3G-SOIT	5477358G3G-SO	5478358G2G-SS
5477307F3G-SS	5477358G3G-SOIT	5478358G2G-SY
5477307F3G-SSIT	5477358G3G-SS	5478358G2G-YY
5477358G1-OO	5477358G3G-SSIT	5478358G3G-OO
5477358G1-SO	5478307F1G-OO	5478358G3G-SO
5477358G1-SS	5478307F1G-SO	5478358G3G-SS
5477358G1G-OO	5478307F1G-SS	5478358G3G-SY
5477358G1G-OOIT	5478307F1G-SY	5478358G3G-YY
5477358G1G-SO	5478307F1G-YY	5481307G4-SS-B
5477358G1G-SOIT	5478307F2G-OO	5481358M4-SS-B
5477358G1G-SS	5478307F2G-SO	5482507G1
5477358G1G-SSIT	5478307F2G-SS	5482507G2
5477358G2-OO	5478307F2G-SY	5482507G3
5477358G2-SO	5478307F2G-YY	5482507G3--G
5477358G2-SS	5478307F3G-OO	5482507G4
5477358G2G-OO	5478307F3G-SO	5482558M1
5477358G2G-OOIT	5478307F3G-SS	5482558M2
5477358G2G-SO	5478307F3G-SY	5482558M3
5477358G2G-SOIT	5478307F3G-YY	5485307G0-FF-V
5477358G2G-SS	5478358G1G-OO	5485307G0-FO-V
5477358G2G-SSIT	5478358G1G-SO	5485307G0-OO-1
5477358G3-OO	5478358G1G-SS	5485307G0-OO-V
5477358G3-SO	5478358G1G-SY	5485307G0-SO-1
5477358G3-SS	5478358G1G-YY	5485307G0-SO-V

5485307G0-SS-V	5485358M8-OO-1	5487307G3-SS
5485307G7-OO-A	5485358M8-OO-A	5487358M1-BB
5485307G7-SO-A	5485358M8-SO-1	5487358M1-OO
5485307G8-FF-A	5485358M8-SO-A	5487358M1-SB
5485307G8-FO-A	5485358M8-SS-A	5487358M1-SO
5485307G8-OO-1	5485358M9-FF-A	5487358M1-SS
5485307G8-OO-A	5485358M9-FO-A	5487358M2-BB
5485307G8-SO-1	5485358M9-OO-1	5487358M2-OO
5485307G8-SO-A	5485358M9-OO-A	5487358M2-SB
5485307G8-SS-A	5485358M9-SO-1	5487358M2-SO
5485307G9-FF-A	5485358M9-SO-A	5487358M2-SS
5485307G9-FO-A	5485358M9-SS-A	5487358M3-BB
5485307G9-OO-1	5487307G1-BB	5487358M3-OO
5485307G9-OO-A	5487307G1-OO	5487358M3-SB
5485307G9-SO-1	5487307G1-SB	5487358M3-SO
5485307G9-SO-A	5487307G1-SO	5487358M3-SS
5485307G9-SS-A	5487307G1-SS	5488307G1-OO
5485358M0-FF-V	5487307G2-BB	5488307G1-SO
5485358M0-FO-V	5487307G2-OO	5488307G1-SS
5485358M0-OO-1	5487307G2-SB	5488307G1-SY
5485358M0-OO-V	5487307G2-SO	5488307G1-YY
5485358M0-SO-1	5487307G2-SS	5488307G2-OO
5485358M0-SO-V	5487307G3-BB	5488307G2-SO
5485358M0-SS-V	5487307G3-OO	5488307G2-SS
5485358M8-FF-A	5487307G3-SB	5488307G2-SY
5485358M8-FO-A	5487307G3-SO	5488307G2-YY

5488307G3-OO	5491307H4-SOHBB	5492507H1-S
5488307G3-SO	5491307H4-SS-B	5492507H2
5488307G3-SS	5491307H4-SSHB1	5492507H2--HB
5488307G3-SY	5491307H4-SSHBB	5492507H2-S
5488307G3-YY	5491307H4-SSVPB	5492507H3
5488358M1-OO	5491307H4G-NO-1	5492507H3--HB
5488358M1-SO	5491307H4G-NO-B	5492507H3-S
5488358M1-SS	5491307H4G-OO-1	549307H8-OO-A
5488358M1-SY	5491307H4G-OO-B	549307H9-OO-A
5488358M1-YY	5491307H4G-OOHB1	5495307H0-FF-V
5488358M2-OO	5491307H4G-OOHBB	5495307H0-OO-V
5488358M2-SO	5491307H4G-SO-1	5495307H0-OOHBV
5488358M2-SS	5491307H4G-SO-B	5495307H0-SO-V
5488358M2-SY	5491307H4G-SOHB1	5495307H0-SOHBV
5488358M2-YY	5491307H4G-SOHBB	5495307H0-SS-V
5488358M3-OO	5491307H4G-SS-1	5495307H0-SSHBV
5488358M3-SO	5491307H4G-SS-B	5495307H0G-OO-1
5488358M3-SS	5491307H4G-SSHB1	5495307H0G-OO-V
5488358M3-SY	5491307H4G-SSHBB	5495307H0G-OOHB1
5488358M3-YY	5491307H4G-SSVPB	5495307H0G-OOHBV
5491307H4-NO-B	5491307HC-FF-V	5495307H0G-SO-1
5491307H4-OO-B	5491307HC-OO-V	5495307H0G-SO-V
5491307H4-OOHB1	5491307HC-OO-V	5495307H0G-SOHB1
5491307H4-OOHBB	5491507H9-25-B	5495307H0G-SOHBV
5491307H4-SO-B	5492507H1	5495307H0G-SS-1
5491307H4-SOHB1	5492507H1--HB	5495307H0G-SS-V

5495307H0G-SSHB1	5495307H7G-SS-1	5495307H9-SO-A
5495307H0G-SSHBV	5495307H7G-SS-A	5495307H9-SOHBA
5495307H7-FF-A	5495307H7G-SSHB1	5495307H9-SS-A
5495307H7-HH-A	5495307H7G-SSHBA	5495307H9-SSHBA
5495307H7-OO-A	5495307H8-FF-A	5495307H9G-OO-1
5495307H7-OOHBA	5495307H8-OO-A	5495307H9G-OO-A
5495307H7-SO-A	5495307H8-OOHBA	5495307H9G-OOHB1
5495307H7-SOHBA	5495307H8-SO-A	5495307H9G-OOHBA
5495307H7-SS-A	5495307H8-SOHBA	5495307H9G-SO-1
5495307H7-SSHBA	5495307H8-SS-A	5495307H9G-SO-A
5495307H7G-FF-1	5495307H8-SSHBA	5495307H9G-SOHB1
5495307H7G-FF-A	5495307H8G-OO-1	5495307H9G-SOHBA
5495307H7G-FH-1	5495307H8G-OO-A	5495307H9G-SS-1
5495307H7G-FH-A	5495307H8G-OOHB1	5495307H9G-SS-A
5495307H7G-FO-1	5495307H8G-OOHBA	5495307H9G-SSHB1
5495307H7G-FO-A	5495307H8G-SO-1	5495307H9G-SSHBA
5495307H7G-HH-1	5495307H8G-SO-A	5495307H0G-OOHB1
5495307H7G-HH-A	5495307H8G-SOHB1	5495307H0G-SOHB1
5495307H7G-OO-1	5495307H8G-SOHBA	5497307H1-OO
5495307H7G-OO-A	5495307H8G-SS-1	5497307H1-OOIT
5495307H7G-OOHB1	5495307H8G-SS-A	5497307H1-SO
5495307H7G-OOHBA	5495307H8G-SSHB1	5497307H1-SOIT
5495307H7G-SO-1	5495307H8G-SSHBA	5497307H1-SS
5495307H7G-SO-A	5495307H9-FF-A	5497307H1-SSIT
5495307H7G-SOHB1	5495307H9-OO-A	5497307H1G-OO
5495307H7G-SOHBA	5495307H9-OOHBA	5497307H1G-OOHB

5497307H1G-OOIT	5497307H2G-SS	5498307H1G-SO-1
5497307H1G-SO	5497307H2G-SSHB	5498307H1G-SOHB
5497307H1G-SO-1	5497307H2G-SSHB1	5498307H1G-SOHB1
5497307H1G-SOHB	5497307H2G-SSIT	5498307H1G-SS
5497307H1G-SOHB1	5497307H3-OO	5498307H1G-SSHB
5497307H1G-SOIT	5497307H3-OOIT	5498307H1G-SSHB1
5497307H1G-SOIT1	5497307H3-SO	5498307H2G-OO
5497307H1G-SS	5497307H3-SOIT	5498307H2G-OOHB
5497307H1G-SSHB	5497307H3-SS	5498307H2G-SO
5497307H1G-SSHB1	5497307H3-SSIT	5498307H2G-SO-1
5497307H1G-SSIT	5497307H3G-OO	5498307H2G-SOHB
5497307H2-OO	5497307H3G-OOHB	5498307H2G-SOHB1
5497307H2-OOIT	5497307H3G-OOIT	5498307H2G-SS
5497307H2-SO	5497307H3G-SO	5498307H2G-SS-1
5497307H2-SOIT	5497307H3G-SO-1	5498307H2G-SSHB
5497307H2-SS	5497307H3G-SOHB	5498307H2G-SSHB1
5497307H2-SSIT	5497307H3G-SOHB1	5498307H3G-OO
5497307H2G-OO	5497307H3G-SOIT	5498307H3G-OO-1
5497307H2G-OOHB	5497307H3G-SOIT1	5498307H3G-OOHB
5497307H2G-OOIT	5497307H3G-SS	5498307H3G-SO
5497307H2G-SO	5497307H3G-SSHB	5498307H3G-SO-1
5497307H2G-SO-1	5497307H3G-SSHB1	5498307H3G-SOHB
5497307H2G-SOHB	5497307H3G-SSIT	5498307H3G-SOHB1
5497307H2G-SOHB1	5498307H1G-OO	5498307H3G-SS
5497307H2G-SOIT	5498307H1G-OOHB	5498307H3G-SS-1
5497307H2G-SOIT1	5498307H1G-SO	5498307H3G-SSHB

5498307H3G-SSHB1	5601505G9--B	5605304E0-SO-V
5498307H8-SS-A	5602504E1	5605304E0-SS-V
54B2507G2	5602504E2	5605304E7-HH-A
54B2507G3	5602504E3	5605304E7-OO-1
54B2507G3--G	5602504E3--G	5605304E7-OO-A
54B2807G2	5602505G1	5605304E7-ODPA
54B2807G3	5602505G2	5605304E7-ODPB
54B2807G3--G	5602505G2--G	5605304E7-SO-1
5521607P25	5602505G3	5605304E7-SO-A
5521607P25--B	5602505G3--G	5605304E7-SS-A
5521607P32--B	5602506F1	5605304E8-OO-1
5521607P32-DGB	5602506F3	5605304E8-OO-A
5521607P40	5602604E1	5605304E8-SO-1
5521607P40--B	5602605G3	5605304E8-SO-A
5591207P9--GAK	5602704E1	5605304E8-SS-A
5591507P9--GAB	5602704E3	5605304E9-OO-1
5592507P1--GA	5602705G1	5605304E9-OO-A
5592507P2--GA	5602705G2	5605304E9-SO-1
5592507P3--GA	5602706G3	5605304E9-SO-A
5592707P2--GA	5602804E1	5605304E9-SS-A
5592707P3--GA	5602805G1	5605305G0-OO-1
5592807P1--GA	5602805G2	5605305G0-OO-V
5592807P2--GA	5602805G3	5605305G0-SO-1
5592807P3--GA	5605304E0-OO-1	5605305G0-SO-V
5601504E7--B	5605304E0-OO-V	5605305G0-SS-V
5601504E9--B	5605304E0-SO-1	5605305G7-FF-A

5605305G7-OO-1	5607304E2-SS	5608305G1-OO
5605305G7-OO-A	5607304E3-BB	5608305G1-SO
5605305G7-SO-1	5607304E3-OO	5608305G1-SS
5605305G7-SO-A	5607304E3-SB	5608305G2-OO
5605305G7-SS-A	5607304E3-SO	5608305G2-SO
5605305G8-OO-1	5607304E3-SS	5608305G2-SS
5605305G8-OO-A	5607305G1-OO	5608305G3-OO
5605305G8-SO-1	5607305G1-SO	5608305G3-SO
5605305G8-SO-A	5607305G1-SS	5608305G3-SS
5605305G8-SS-A	5607305G2-FF	5621504E9--B
5605305G8-SS-K	5607305G2-OO	5621505G9--B
5605305G9-OO-1	5607305G2-SO	5621506A9--B
5605305G9-OO-A	5607305G2-SS	5621507A9--B
5605305G9-SO-1	5607305G2-SSDP	5622104E1
5605305G9-SO-A	5607305G3-OO	5622106A1
5605305G9-SS-A	5607305G3-SO	5622106A2
5607304E1-BB	5607305G3-SS	5622107A1
5607304E1-OO	5608304E1-OO	5622107A2
5607304E1-SB	5608304E1-SO	5622107A3
5607304E1-SO	5608304E1-SS	5622503E2
5607304E1-SS	5608304E2-OO	5622503E3
5607304E2-BB	5608304E2-SO	5622503E4
5607304E2-FF	5608304E2-SS	5622504E1
5607304E2-OO	5608304E3-OO	5622504E1--H
5607304E2-SB	5608304E3-SO	5622504E2
5607304E2-SO	5608304E3-SS	5622504E3

5622504E3--C	5622704E1	5625304E7-OO-1
5622504E3--SC	5622704E3	5625304E7-OO-A
5622504E4	5622704E3--C	5625304E7-SO-1
5622504E4--SC	5622705G1	5625304E7-SO-A
5622505G1	5622705G2	5625304E7-SS-A
5622505G2	5622705G3	5625304E7-WW-A
5622505G3	5622706A1	5625304E8-FF-A
5622505G3--SC	5622706F1	5625304E8-OO-1
5622506A1	5622706F3	5625304E8-OO-A
5622506A2	5622707A1	5625304E8-SO-1
5622506A3	5622707A2	5625304E8-SO-A
5622506A3--G	5622707A3	5625304E9-OO-1
5622506F1	5622804E1	5625304E9-OO-A
5622506F2	5622804E2	5625304E9-SO-1
5622506F3	5622804E3	5625304E9-SO-A
5622507A1	5622805G1	5625304E9-SS
5622507A2	5622805G2	5625304E9-SS-A
5622507A3	5622805G3	5625304E9-SS-V
5622603E3	5622806A1	5625305G0-OO-1
5622604E1	5625304E0-OO-1	5625305G0-OO-V
5622604E3	5625304E0-OO-V	5625305G0-SO-1
5622604E3--C	5625304E0-SO-1	5625305G0-SO-V
5622605G3	5625304E0-SO-V	5625305G0-SS-V
5622606F3	5625304E0-SS-V	5625305G7-OO-1
5622606F3--C	5625304E0-SSDPV	5625305G7-SO-1
5622703E3	5625304E4-OO-B	5625305G7-SO-A

5625305G8-OO-1	5625306A9-OO-1	5627304E1-BB
5625305G8-OO-A	5625306A9-OO-A	5627304E1-FF
5625305G8-SO-1	5625306A9-SO-1	5627304E1-OO
5625305G8-SO-A	5625306A9-SS-A	5627304E1-SB
5625305G8-SS-A	5625306F0-OO-V	5627304E1-SO
5625305G9-OO-1	5625306F7-OO-A	5627304E1-SS
5625305G9-OO-A	5625306F8-OO-A	5627304E1-SSDP
5625305G9-SO-1	5625306F9-OO-A	5627304E2-BB
5625305G9-SO-A	5625307A0-FF-V	5627304E2-OO
5625305G9-SS-A	5625307A0-OO-1	5627304E2-SB
5625305G9-WW-A	5625307A0-OO-V	5627304E2-SO
5625306A0-OO-1	5625307A0-SO-1	5627304E2-SS
5625306A0-OO-V	5625307A0-SO-V	5627304E3-BB
5625306A0-SO-1	5625307A0-SS-V	5627304E3-FF
5625306A0-SO-V	5625307A7-OO-1	5627304E3-OO
5625306A0-SS-V	5625307A7-OO-A	5627304E3-SB
5625306A7-FF-A	5625307A7-SO-1	5627304E3-SO
5625306A7-HH-A	5625307A7-SS-A	5627304E3-SS
5625306A7-OO-1	5625307A8-OO-1	5627305G1-BB
5625306A7-OO-A	5625307A8-OO-A	5627305G1-FF
5625306A7-SO-1	5625307A8-SO-1	5627305G1-OO
5625306A7-SS-A	5625307A8-SS-A	5627305G1-SB
5625306A8-OO-1	5625307A9-OO-1	5627305G1-SO
5625306A8-OO-A	5625307A9-OO-A	5627305G1-SS
5625306A8-SO-1	5625307A9-SO-1	5627305G2-BB
5625306A8-SS-A	5625307A9-SS-A	5627305G2-FF

5627305G2-OO	5628305G1-OO	5642503A3--AV
5627305G2-SB	5628305G1-SO	5642503A4--AV
5627305G2-SO	5628305G1-SS	5642703A3--AV
5627305G2-SS	5628305G2-OO	5642803A1--AV
5627305G3-BB	5628305G2-SO	5642803A2--AV
5627305G3-FF	5628305G2-SS	5642803A3
5627305G3-OO	5628305G3-OO	5642803A3--AV
5627305G3-SB	5628305G3-SO	5664002J3--KG
5627305G3-SO	5628305G3-SS	5834005J3--AV
5627305G3-SS	5632504A1--AV	5834005J5--AV
5627306A1-BB	5632504A2--AV	5864003J4--WF
5627306A1-SB	5632504A3--AV	BIOSTAT-KIT1
5627306A2-BB	5632704A3--AV	BIOSTAT-KIT2
5627306A2-SB	5632804A1--AV	BIOSTAT-KIT3
5627306A2-SS	5632804A2--AV	CTR-01
5627306A3-BB	5632804A3--AV	CTR-04
5627306A3-SB	5634002J1	FTR-01
5628304E1-OO	5634002J2	IN-61704-117-K01
5628304E1-SO	5634002J3	IN-61704-119-K01
5628304E1-SS	5634002J4	IN-61704-502-K01
5628304E2-OO	5634003J1	IN-61704-515-O03
5628304E2-SO	5634003J2	IN-61704-517-O03
5628304E2-SS	5634003J3	IN-61704-518-K01
5628304E3-OO	5634003J4	IN-61704-525-O03
5628304E3-SO	5642503A1--AV	IT-CA002PS
5628304E3-SS	5642503A2--AV	IT-CA005GF

J1-5152507T3---E	UNIVESSEL-KIT1	5995307G7G-OO--A
J1-5181307T4--OO	UNIVESSEL-KIT2	5995307G7G-SO--A
J1-5181507T9	5445307H7--SSDAA	5995307G8G-FF--A
J1-5182607T2-LGT	5441307H4--SODPB	5995307G8G-FO--A
J1-5182658T2-LGT	5235306D7--SSDPA	5995307G8G-OO--A
J1-5185307T8--SS	5605304E7--SSDPA	5995307G8G-SS--A
J1-5185307T9	5997307G1G-SO	5995307G8G-SO--A
J1-5185307T9-OO	5997307G1G-OO	5995307G9G-FF--A
J1-5231506D9	5997307G2G-SO	5995307G9G-FO--A
J1-5231807H9	5997307G2G-OO	5995307G9G-OO--A
J1-5235307H7--SS	5997307G3G-SO	5995307G9G-SO--A
J1-5235307H9--SS	5997307G3G-OO	5995307G0G-FF--V
J1-5441306G5--OO	5998307G1G-SS	5995307G0G-FO--V
J1-5441307H4--OO	5998307G1G-SO	5995307G0G-OO--V
J1-5445307H7--SS	5998307G1G-OO	5995307G0G-SS--V
J1-5445307H8--OO	5998307G2G-SS	5995307G0G-SO--V
J1-5601504E9	5998307G2G-SO	5992507G1
J1-B65	5998307G2G-OO	5992507G2
T5601304E3--SO	5998307G3G-SS	5992507G3
T5621305G3--SO	5998307G3G-SO	5997307G1G-SS
U1-5441307H1OOAC	5998307G3G-OO	5997307G2G-SS
U1-5441307H4GOOT	5991307G4G-SO--B	5997307G3G-SS
U1-5441307H5GOOT	5991307G4G-OO--B	5991307G4G-SS--B
U1-5441358K1SSAC	5991307G4G-SSVPB	5995307G7G-SS--A
U1-5441358K2SSAC	5995307G7G-FF--A	5995307G9G-SS--A
U1-5441358K3SSAC	5995307G7G-FO--A	5997307G1G-SS--1

5997307G1G-SO--1	5995307G8G-FO--1
5997307G1G-OO--1	5995307G8G-OO--1
5997307G2G-SS--1	5995307G8G-SS--1
5997307G2G-SO--1	5995307G8G-SO--1
5997307G2G-OO--1	5995307G9G-FF--1
5997307G3G-SS--1	5995307G9G-FO--1
5997307G3G-SO--1	5995307G9G-OO--1
5997307G3G-OO--1	5995307G9G-SS--1
5998307G1G-SS--1	5995307G9G-SO--1
5998307G1G-SO--1	5995307G0G-FF--1
5998307G1G-OO--1	5995307G0G-FO--1
5998307G2G-SS--1	5995307G0G-OO--1
5998307G2G-SO--1	5995307G0G-SS--1
5998307G2G-OO--1	5995307G0G-SO--1
5998307G3G-SS--1	5991507G7-----B
5998307G3G-SO--1	5991507G8-----B
5998307G3G-OO--1	5991507G9-----B
5991307G4G-SS--1	
5991307G4G-SO--1	
5991307G4G-OO--1	
5995307G7G-FF--1	
5995307G7G-FO--1	
5995307G7G-OO--1	
5995307G7G-SS--1	
5995307G7G-SO--1	
5995307G8G-FF--1	