



Guide to Recycling and
Responsible Disposal of Lab
Filtration and Chromatography
Products

Simplifying Progress

SARTORIUS

Introduction

Disposable filtration and chromatography consumables are essential tools for scientific research across multiple disciplines. These products can help ensure consistent results, prevent carryover or sample contamination, and improve laboratory safety by eliminating the need for chemical-based sanitization and cleaning processes.

While single-use consumables offer clear benefits, they can also pose challenges for waste management and environmentally friendly disposal. Sartorius is committed to sustainability and offers innovative products that help to maximize throughput and some that can even be reused multiple times to reduce the number of consumables needed in your separation workflows.

In addition, based on a literature review of regional and global standards¹⁻⁵, this guide lists the materials of each component and the packaging of Sartorius lab filtration and chromatography products. It is intended to help you select the most appropriate waste management practices for disposing of our consumables at the end of their useful life.

Many of the materials in Sartorius lab filtration and chromatography products are easy to separate and may therefore be easier to recycle. These are clearly indicated in this guide. For materials that cannot be easily separated (e.g., membranes and housings), Sartorius recommends incineration to recover energy from the waste, preferably in a facility designed to minimize emissions.

Resin Codes

Resin Code in China (GB 16288-2008)	Description
1	PET
2	HDPE
3	PVC
4	LDPE
5	PP
6	PS
7	ABS
14	CA
22	CP
23	CTA
38	MBS
43	PA
58	PC
76	PES
94	PMMA
96	PMP
102 / 103	PP-E / PP-HI
110	PTFE
120	PVDF
125	SAN
126	SB

Resin Code in America (ASTM D7611 / D7611M)	Description
1	PET / PETE
2	PE-HD / HDPE
3	PVC
4	PE-LD / LDPE
5	PP
6	PS
PA	PA
ABS	ABS
7	All other Plastics
20	PAP (corrugated)
21	PAP
40	FE
41	ALU
95	Glass / Plastics

Disclaimers

- The information presented in this document is based on raw material data for Sartorius lab filtration and chromatography products and a literature review of selected SAC^{1,2}, ASTM³ and ISO⁵ standards. It should not be considered definitive, especially in regions where these standards have not been adopted.
- In Europe, resin codes for plastics are not governed by specific European standards but are mostly adopted from the Plastics Industry Association based in the USA.
- In general, recyclability in North America is aligned with European regulations.
- Recycling capabilities vary around the world. You should confirm which materials can be recycled in your region to determine the most appropriate waste management approach for disposal of the products listed in this guide.
- Hazardous waste must be disposed of according to your local guidelines and | or legislation and should not be recycled.

References

1. Standardization Administration of China. (2008). GB 16288-2008: Marking of plastic products. Beijing: SAC.
2. Standardization Administration of China. (2001). GB 18455-2001: Packaging recycling mark. Beijing: SAC.
3. ASTM International. (2019). ASTM D7611 / D7611M-19: Standard practice for coding plastic manufactured articles for resin identification. West Conshohocken, PA: ASTM International. https://doi.org/10.1520/D7611_D7611M-19
4. International Organization for Standardization. (2008). ISO 15270:2008 Plastics --Guidelines for the recovery and recycling of plastics waste. Geneva: ISO. <https://www.iso.org/standard/42845.html>
5. International Organization for Standardization. (2020). ISO/TR 23891:2020 Plastics --Guidance on compounding and converting for recycling of plastics. Geneva: ISO. <https://www.iso.org/standard/76191.html>

Minisart®

Component	Click to Expand
Minisart® NML & NML Plus	
Minisart® High Flow	
Minisart® NY & NY Plus	
Minisart® SRP	
Minisart® RC	



Minisart®



Minisart® NML & NML Plus

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Cellulose Acetate (CA)	14	-	-
Housing	Methyl methacrylate-butadiene-styrene (MBS)	38	ABS	-
Pre-filter	Glass Fiber (GF)	14**	95 (Glass/Plastic)	-
Primary Packaging (K Packs)	Blister: APET Tyvek	-	7	Yes
Primary Packaging (Q Packs)	PA/PE 20/80 mm	43	PA	Yes
Secondary Packaging (K Packs)	Cardboard	FB**	20	Yes
Secondary Packaging (Q Packs)	Non-Corrugated Cardboard	NCFB**	21	Yes

Minisart®



Minisart® High Flow

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Methyl methacrylate-butadiene-styrene (MBS)	38	ABS	-
Primary Packaging (K Packs)	Blister: APET Tyvek	-	7	Yes
Primary Packaging (Q Packs)	PA/PE 20/80 mm	43	PA	Yes
Secondary Packaging (K Packs)	Cardboard	FB**	20	Yes
Secondary Packaging (Q Packs)	Non-Corrugated Cardboard	NCFB**	21	Yes

Minisart®

Minisart® NY & NY Plus

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Nylon (NY)	43	PA	-
Housing	Polypropylene (PP)	5	5	-
Pre-filter	Glass Fiber (GF)	14**	95 (Glass/ Plastic)	-
Primary Packaging (K Packs)	Polyethylene (PE)	2/4	2/4	Yes
Primary Packaging (Q Packs)	PA/PE 20/80 mm	43	PA	Yes
Secondary Packaging (K Packs)	Cardboard	FB**	20	Yes
Secondary Packaging (Q Packs)	Non-Corrugated Cardboard	NCFB**	21	Yes

Minisart®



Minisart® SRP

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polytetrafluoroethylene (PTFE)	110	-	-
Housing	Polypropylene (PP)	5	5	-
Primary Packaging (K Packs)	Blister: APET Tyvek	-	-	Yes
Primary Packaging (non-sterile)	Polyethylene (PE)	2/4	2/4	Yes
Primary Packaging (Q Packs)	PA/PE 20/80 mm	43	PA	Yes
Secondary Packaging (K Packs)	Cardboard	FB**	20	Yes
Secondary Packaging (Q Packs)	Non-Corrugated Cardboard	NCFB**	21	Yes

Minisart®

Minisart® RC

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Regenerated Cellulose (RC)	-	-	-
Housing	Polypropylene (PP)	5	5	-
Primary Packaging (K Packs)	Polyethylene (PE)	2/4	2/4	Yes
Primary Packaging (Q Packs)	PA/PE 20/80 mm	43	PA	Yes
Secondary Packaging (K Packs)	Cardboard	FB**	20	Yes
Secondary Packaging (Q Packs)	Non-Corrugated Cardboard	NCFB**	21	Yes

Claristep®

Component	Click to Expand
Claristep® Single Tray	
Claristep® Filters	



Claristep®

Claristep® Single Tray

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Primary Packaging	Polyethylene (PE)	2/4	2/4	Yes
Secondary Packaging	Corrugated Cardboard	FB*	20	Yes

Claristep®

Claristep® Filters

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Regenerated Cellulose (RC)	-	-	-
Housing	Polypropylene (PP)	5	5	-
Primary Packaging	Polyethylene (PE)	2/4	2/4	Yes
Secondary Packaging (96 Filters)	Cardboard	NCFB**	21	Yes
Secondary Packaging (480 Filters)	Corrugated Cardboard	FB**	20	Yes

Sartolab® P20

Component

Click to Expand

Sartolab® P20



Sartolab® P20 Plus &
Sartolab® P20 Prefilter



Sartolab® P20

Sartolab® P20

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	7	-
Venting Filter	Polytetrafluoroethylene (PTFE)	110	7	-
Filling Bell & Lid	Polycarbonate (PC)	58	7	-
Housing	Polycarbonate (PC)	58	7	-
Primary Packaging	Polyethylene (PE)	2/4	2/4	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes

Sartolab® P20

Sartolab® P20 Plus & Sartolab® P20 Prefilter

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Venting Filter	Polytetrafluoroethylene (PTFE)	110	-	-
Prefilter	Quartz Microfibers	14**	95 (Glass/ Plastic)	-
Filling Bell & Lid	Polycarbonate (PC)	58	-	Yes
Housing	Polycarbonate (PC)	58	-	-
Primary Packaging	Polyethylene (PE)	2/4*	2 /4	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes

Sartolab® RF|BT

Component	Click to Expand
Sartolab® RF BT	
Sartolab® Multistation	



Sartolab® RF|BT

Sartolab® RF|BT

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Funnel, Lid, and Bottle	Polystyrene (PS)	6	6	Yes
Adapter, Tube Adapter, Screw Cap	High Density Polythylene (HDPE)	2	2	Yes
Conical Tube	Polypropylene (PP)	5	5	Yes
Gasket (funnel)	Polyethylene (PE)	2/4	2/4	Yes
Primary Packaging	Polyethylene terephthalate (PET)	1	1	Yes
	Polyethylene (PE)	2/4	2/4	Yes
	Polyamid (PA)	43	PA	Yes
Secondary Packaging	Corrugated Fiberboard	FB**	20	Yes

Sartolab® RF|BT

×

Sartolab® Multistation

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Primary Packaging	Corrugated Fiberboard	FB**	20	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes
General Material	ABS Aluminium Stainless steel	7	41	-

Sartoclear Dynamics® Lab

Component	Click to Expand
Sartoclear Dynamics® Lab P15	+
Sartoclear Dynamics® Lab V50-1000 (filtration devices)	+
Sartoclear Dynamics® Lab DE Filter Aid (filter aid boxes)	+



Sartoclear Dynamics® Lab

Sartoclear Dynamics® Lab P15

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Prefilter	Quartz microfibers	14**	95 (Glass/ Plastic)	-
Housing	Polycarbonate (PC)	58	7	-
Syringe	Polypropylene (PP)	5	5	Yes
Caps	Polyamide (PA)	43	PA	Yes
Filling tube	Polypropylene (PP)	5	5	Yes
Filter Aid	Diatomaceous Earth (DE)	-	-	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes
Syringe Packaging	Polypropylene (PP)	5	5	Yes
Filter Packaging	Polyethylene (PE)	2/4	2/4	Yes

Sartoclear Dynamics® Lab

Sartoclear Dynamics® LabV50-1000



Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Funnel, Lid, and Bottle	Polystyrene (PS)	6	6	Yes
Adapter, Tube Adapter, Screw Cap	High Density Polythylene (HDPE)	2/4	2/4	Yes
Conical Tube	Polypropylene (PP)	5	5	Yes
Gasket (funnel)	Polyethylene (PE)	2/4	2/4	Yes
Primary Packaging	Polyethylene terephtalate (PET)	1	1	Yes
	Polethylene (PE)	2/4	2/4	Yes
	Polyamid (PA)	43	PA	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes

Sartoclear Dynamics® Lab

Sartoclear Dynamics® Lab DE Filter Aid

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Filter Aid	Diatomaceous Earth (DE)	-	-	Yes
Primary Packaging (pouch)	Aluminium (ALU), Polyamid (PA), Polyethylene (PE)	43, 2/4	41, 2/4	Yes
Secondary Packaging	Corrugated Fiberboard	FB**	20	Yes

Vivaspin®

Component	Click to Expand
Vivaspin® 500	+
Vivaspin® 2	+
Vivaspin® 6	+
Vivaspin® 15R Endotest	+
Vivaspin® 20	+
Vivaspin® 100	+
Vivaspin® Filtrate	+
Vivaspin® Turbo 4	+
Vivaspin® Turbo 15	+



Vivaspin®

Vivaspin® 500

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	Yes
Housing	Polycarbonate (PC)	58	7	-
Tube	Polypropylene (PP)	5	5	Yes
Primary Packaging	Polyethylene (PE)	4	4	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® 2



Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
	Cellulose triacetate (CTA)	23	-	-
	Hydrosart® regenerated cellulose (RC)	-	-	-
Gasket	Silicone (SIL)	-	-	-
Housing	Polycarbonate (PC)	58	7	-
Tube	Polycarbonate (PC)	58	7	Yes
Cap	Polycarbonate (PC)	58	7	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® Filtrate

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
	Cellulose triacetate (CTA)	23	-	-
Tube	Polystyrene (PS)	6	6	Yes
Housing	Styrene acrylonitrile (SAN)	125	7	-
Cap	Polyethylene (PE)	2/4	2/4	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes

Vivaspin®

Vivaspin® Turbo 4

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Styrene butadiene copolymer (SBC)	126	-	-
Tube	Polypropylene (PP)	5	5	Yes
Cap	Polypropylene (PP)	5	5	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® 6

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Polycarbonate (PC)	58	7	-
Tube	Polycarbonate (PC)	58	7	Yes
Cap	Polypropylene (PP)	5	5	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® 15R | Endotest

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Hydrosart® regenerated cellulose (RC)	-	-	-
	Cellulose triacetate (CTA)	23	-	-
Housing	Polycarbonate (PC)	58	7	-
Gasket	Silicone (SIL)	-	-	-
Tube	Polycarbonate (PC)	58	7	Yes
Cap	Polypropylene (PP)	5	5	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® Turbo 15

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76/-	-/-	-
	Regenerated cellulose (RC)	-	-	-
Housing	Styrene butadiene copolymer (SBC)	126	7	-
Tube	Polypropylene (PP)	5	5	Yes
Cap	Polypropylene (PP)	5	5	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® 20

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Polycarbonate (PC)	58	7	-
Tube	Polycarbonate (PC)	58	7	-
Cap	Polypropylene (PP)	5	5	-
Pressure Head	Polyoxymethylene (POM)/Aluminium (ALU)	98/-	7/41	-
Diafiltration Cup	Polypropylene (PP)	5	5	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivaspin®

Vivaspin® 100

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Polycarbonate (PC)	58	7	-
Tube	Polycarbonate (PC)	58	7	Yes
Cap	Polypropylene (PP)	5	5	Yes
Pressure Head	Polyoxymethylene (POM)/Aluminium (ALU)	98/-	7/41	-
Primary Packaging	Polyethylene (PE)	4	4	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes

Vivaflow®

Component	Click to Expand
Vivaflow® SU	+
Vivaflow® 50R	+
Vivaflow® 200	+



Vivaflow®

Vivaflow® SU

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Fittings	Polyamide (PA)	43	PA	Yes
	Polypropylene (PP)	5	5	Yes
Gasket	Silicone (SIL)	-	-	-
Housing	Polycarbonate (PC)	58	7	-
Membrane	Hydrosart® regenerated cellulose (RC)	-	-	-
	Polyethersulfone (PES)	76	-	-
Membrane Support	Polyethylene (PE)	4	4	-
Pressure Indicator	Polyamide (PA)	43	PA	Yes
	Polyoxymethylene (POM)	98	7	Yes
	Polypropylene (PP)	5	5	Yes
	Silicone (SIL)	-	-	Yes
	Stainless steel (SS)	-	-	Yes
Tubing	Polyvinyl chloride (PVC)	3	3	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB	21	Yes
	Polyethylene	4	4	Yes
Secondary Packaging	Corrugated fiberboard	FB	20	Yes

Vivaflow®

Vivaflow® 50R

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Fittings	Polyamide (PA)	43	PA	Yes
	Polypropylene (PP)	5	5	Yes
Gasket	Silicone (SIL)	-	-	-
Housing	Polymethyl methacrylate (PMMA)	94	7	-
Membrane	Hydrosart® regenerated cellulose (RC)	-	-	-
Membrane Support	Polypropylene (PP)	5	5	-
Pressure Indicator	Polyamide (PA)	43	PA	Yes
	Polyoxymethylene (POM)	98	7	Yes
	Polypropylene (PP)	5	5	Yes
	Silicone (SIL)	-	-	Yes
	Stainless steel (SS)	-	-	Yes
Tubing	Polyvinyl chloride (PVC)	3	3	Yes
Primary Packaging	Polyethylene	4	4	Yes
Secondary Packaging	Corrugated fiberboard	FB	20	Yes

Vivaflow®

Vivaflow® 200



Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Fittings	Polyamide (PA)	43	PA	Yes
	Polypropylene (PP)	5	5	Yes
Gasket	Silicone (SIL)	-	-	-
Housing	Polymethyl methacrylate (PMMA)	94	7	-
Membrane	Hydrosart® regenerated cellulose (RC)	-	-	-
	Polyethersulfone (PES)	76	-	-
Membrane Support	Polypropylene (PP)	5	5	-
Pressure Indicator	Polyamide (PA)	43	PA	Yes
	Polyoxymethylene (POM)	98	7	Yes
	Polypropylene (PP)	5	5	Yes
	Silicone (SIL)	-	-	Yes
	Stainless steel (SS)	-	-	Yes
Tubing	Polyvinyl chloride (PVC)	3	3	Yes
Primary Packaging	Polyurethane (PU)	113	7	Yes
	Polyethylene	4	4	Yes
Secondary Packaging	Corrugated fiberboard	FB	20	Yes

Vivapore®

Component	Click to Expand
Vivapore®	



Vivapore®

Vivapore®

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Styrene acrylonitrile (SAN)	125	7	-
Primary Packaging	Polyethylene (PE)	4	4	Yes
Secondary Packaging	Corrugated fiberboard	FB	20	Yes
Reservoir	Polypropylene (PP)	5	5	Yes
Stand	Acrylonitrile butadiene styrene (ABS)	7	ABS	Yes

Vivacon®

Component

Click to Expand

Vivacon® 500



Vivacon® 2



Vivacon®

Vivacon® 500

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Hydrosart® regenerated cellulose (RC)	-	-	-
	Cellulose acetate (CA)	14	-	-
Housing	Polycarbonate (PC)	58	7	-
Tube	Polypropylene (PP)	5	5	Yes
Primary Packaging	Polyethylene (PE)	4	4	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	21/20	Yes

Vivacon[®]

Vivacon[®] 2

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Hydrosart [®] regenerated cellulose (RC)	-	-	-
	Cellulose acetate (CA)	14	-	-
Housing	Polycarbonate (PC)	58	7	-
Tube	Polypropylene (PP)	5	5	Yes
Caps	Polypropylene (PP)	5	5	Yes
Primary Packaging	Polyethylene (PE)	4	4	Yes
Secondary Packaging	Corrugated or non-corrugated fiberboard, depending on pack size	NCFB/FB**	20/21	Yes

Depth Filters

Component	Click to Expand
Filter Papers	+
Glass- & Quartz Microfiber Filters (starting with FT-*)	+
Glass- & Quartz Microfiber Filters (starting with 134* or Q3400*)	+



Depth Filters

✕

Filter Papers

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Filter Material	Alpha-Cellulose, Cotton Linters	14	-	-
Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes

Depth Filters

Glass- & Quartz Microfiber Filters (starting with FT-*)

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Filter Material	Glass, Quartz	14**	95	-
Packaging Inlay	Polystyrene (PS)	6	6	-
Packaging Box	Non-Corrugated Fiberboard	NCFB**	21	Yes
Secondary Packaging	Non-Corrugated Fiberboard	NCFB**	21	Yes

Depth Filters

Glass- & Quartz Microfiber Filters (starting with 134* or Q3400*)

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Filter Material	Glass, Quartz	14**	95	-
Packaging Outer Box	Corrugated Fiberboard	FB**	20	Yes
Packaging Box & Blister	Polystyrene (PS)	6	6	-

Sartobind®

Component	Click to Expand
Vivapure® IEX Mini	+
Vivapure® IEX Maxi	+
Vivaclear	+
Sartobind® IEX IDA Lab	+
Sartobind® Rapid A Lab	+



Sartobind®

Vivapure® IEX Mini

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Tube	Polypropylene (PP)	5	5	Yes
Housing	Polycarbonate (PC)	58	7	-
Membrane	Regenerated cellulose (RC)	-	-	-
Primary Packaging	Polyethylene (PE)	2/4	2/4	Yes
Secondary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes

Sartobind®

Vivapure® IEX Maxi

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Tube	Polypropylene (PP)	5	5	Yes
Housing	Polycarbonate (PC)	58	7	-
Membrane	Regenerated cellulose (RC)	-	-	-
Cap	Polypropylene (PP)	5	5	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes

Sartobind®

Vivaclear

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Membrane	Polyethersulfone (PES)	76	-	-
Housing	Polypropylene (PP)	5	5	-
Tube	Polypropylene (PP)	5	5	Yes
Primary Packaging	Polyethylene (PE)	4	4	Yes
Secondary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes

Sartobind®

Sartobind® IEX | IDA Lab

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Housing	Polypropylene (PP)	5	5	-
Membrane	Regenerated cellulose (RC)	-	-	-
Caps	Polycarbonate (PC)	58	7	Yes
Adapters	Polyetheretherketone (PEEK)	68	-	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
Secondary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes

Sartobind®

Sartobind® Rapid A Lab

Component	Material	Resin Code		Easy to Separate
		China	EU/USA	
Housing	Polypropylene (PP)	5	5	-
Membrane	Agarose	-	-	-
Caps	Polycarbonate (PC)	58	7	Yes
Adapters	Polyetheretherketone (PEEK)	68	-	Yes
Primary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes
	Aluminium composite (ALU/OPA/PE)	-/-/4	41/-/4	-
Secondary Packaging	Non-corrugated fiberboard	NCFB**	21	Yes

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0



For further contacts, visit
www.sartorius.com

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906