



Legend	
Symbol	Description
↑	Fluid path
①	1 cable tie
②	2 cable ties

Description	FLEXSAFE® 2D 20L (TPE/AC)		Design.	
Product P/N	FLS301755 Revision 00		Name	Date
Customer P/N	C-305	Engineered By	Tania Razafimahefa	09/10/2017
Page : 1/4	Product	SSB Approved		

N°	Description	PN	Material
1	ASEPTIQUIK-G 3/8 (HT)	AC114462	PC/PES/PTFE/SI
2	PINCH CLAMP	CL100877	PBT
3	PLUG FEMALE LUER	PR100662	PE
4	CLAVE CONNECTOR	ST107530	ABS/PC/SI
5	LUER FEMALE 1/4"	TC112313	PP
6	TUBE 3/8"ID X 5/8"OD L20" (500MM)	TU111933	TPE C-Flex 374®
7	TUBE 1/4"ID X 7/16"OD L6" (150MM)	TU112236	TPE C-Flex 374®
8	LOT NUMBER LABEL	LB114999	PAPER
9	HOT STAMP FILM (BLACK)	MF106880	PET
10	WELDED HANDLE (ORANGE)	RP114443	PE
11	BAG CHAMBER	SF113810	PE
12	BOAT PORT 3/8" - 1/4" - 3/8" HB	TC113391	PE

		<small>This document is the property of SARTORIUS STEDIM BIOTECH. and must not be disclosed to third party without the written consent of SARTORIUS STEDIM BIOTECH. This drawing represents the components only schematically to illustrate the function and/or assembly of the equipment/device and to enable the customer the check the component's completeness. The semblance of the real components may therefore deviate from the drawing.</small>	
Description	FLEXSAFE® 2D 20L (TPE/AC)		Design.
Product P/N	FLS301755 Revision 00		Name
Customer P/N	C-305	Engineered By	Date
Page : 2/4	Product BOM	SSB Approved	Tania Razafimahefa 09/10/2017