



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



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 to check the component's completeness. The semblance of the real components may therefore deviate from the drawing.

Description	FLEXSAFE® 2D 50L (TPE)		Design.	
Product P/N	FLS301416 Revision 00		Name	Date
Customer P/N	C-175	Engineered By	Tania Razafimahefa	18/08/2017
Page :	Product	SSB Approved		

N°	Description	PN	Material
1	COUPLING FEMALE 3/8"	CK100771	PC
2	PINCH CLAMP	CL100877	PBT
3	SEALING PLUG	PL100666	PC/SI
4	PRESS IN PLUG 3/8"	PL108501	PP
5	TUBE 3/8"ID X 5/8"OD L40" (1000MM)	TU112108	TPE C-Flex 374®
6	LOT NUMBER LABEL	LB114999	PAPER
7	HOT STAMP FILM (BLACK)	MF106880	PET
8	BAG CHAMBER	SF113810	PE
9	BOAT PORT 3/8" - Plug - 3/8" HB	TC113394	PE



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 to check the component's completeness. The semblance of the real components may therefore deviate from the drawing.

Description	FLEXSAFE® 2D 50L (TPE)		Design.
Product P/N	FLS301416 Revision 00		Name
Customer P/N	C-175	Engineered By	Date
Page :	Product BOM	SSB Approved	Tania Razafimahefa 18/08/2017