

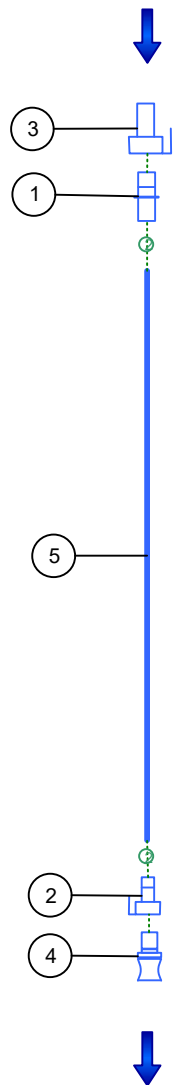


Legend	
Symbol	Description
	Fluid path
	2 cable ties




sartorius stedim
biotech

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	TRANSFER SET MPX/MPX (TPE)		Design.	
Product P/N	FBT123460 Revision 00		Name	Date
Customer P/N	C-032	Engineered By	T.RAZAFIMAHEFA	25/10/2017
Page : 1/4	Product	SSB Approved		

N°	Description	PN	Material
1	COUPLING MALE 1/2"	CK100783	PC/SI
2	COUPLING FEMALE 1/2"	CK108509	PC
3	SEALING CAP	PL100657	PC
4	SEALING PLUG	PL100668	PC / SI
5	TUBE 1/2"ID X 3/4"OD L60" (1500MM)	TU110774	TPE C-Flex 374®

		<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	TRANSFER SET MPX/MPX (TPE)		Design.	
Product P/N	FBT123460 Revision 00		Name	Date
Customer P/N	C-032	Engineered By	T.RAZAFIMAHEFA	25/10/2017
Page : 2/4	Product BOM	SSB Approved		