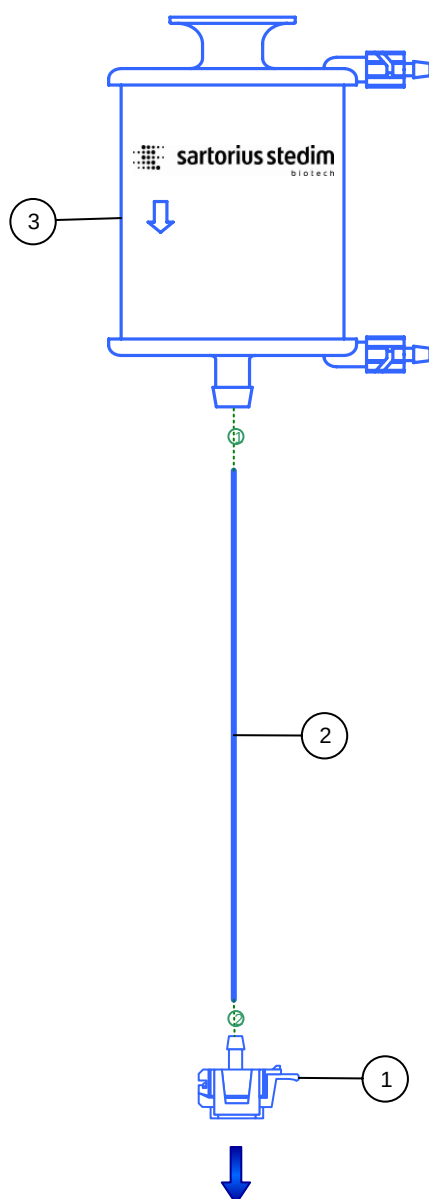


Legend	
Symbol	Description
↑	Fluid path
①	1 cable tie
②	2 cable ties



Description	TRANSFER SET SARTOPORE®/AC (TPE)		Design.	
Product P/N	FBT302624 Revision 00		Name	Date
Customer P/N	C-286	Engineered By	Tania Razafimahefa	31/01/2018
Page : 1/4	Product	SSB Approved		

