

Specifications for Midi Plus Filter Type LH-7129120

Filter type	Midi Plus filter, hydrophobic, non-sterile
Filter material	PTFE membrane, PP reinforced
Pore size	0,45 µm
Filter diameter	28,9 mm
Filtration area	2,9 cm ²
Housing material	Polypropylene, transparent
Bubble Point (visual, ethanol wetted)	≥ 0,9 bar
Burst pressure	> 6,0 bar
Flow rate for air	average ≥ 2,0 L/min at 0,1 bar
Water penetration point	> 3,0 bar

December 12th, 2016



Seppo Riikonen
Director, Quality & Process Development

Sartorius Biohit Liquid Handling Oy