

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, water wetted)	$\geq 0,9$ bar
Burst pressure	$> 6,0$ bar
Flow rate for air	average $\geq 2,0$ l/min at 0,1 bar
Water penetration point	$> 3,0$ bar

January 26th, 2022



Ninna Koivunen
Head of Quality

Sartorius Biohit Liquid Handling Oy