

## Specifications for Midi Plus Filter Type LH-7129131

Filter type	Midi Plus filter, hydrophobic, sterile
Filter material	PTFE membrane, PP reinforced
Pore size	0,2 µm
Filter diameter	28,9 mm
Filtration area	2,9 cm <sup>2</sup>
Housing material	Polypropylene, transparent
Bubble Point (visual, water wetted)	≥ 1,4 bar
Burst pressure	> 6,0 bar
Flow rate for air	average ≥ 1,5 l/min at 0,1 bar
Water penetration point	> 4,0 bar
Endotoxin test (LAL)	Acc. to cUSP < 0,25 EU/ml per filter
Sterile filtration capability (Bacterial challenge test)	All filtrates sterile at 4,0 bar with 10 <sup>7</sup> Brevundimonas diminuta CFU/cm <sup>2</sup>
Sterilization and Sterility	Validated EO-sterilization process, monitored by spore strips.

January 28<sup>th</sup>, 2022



Ninna Koivunen  
Head of Quality

Sartorius Biohit Liquid Handling Oy