






Decision Aid

Select the appropriate Milliflex Oasis® filtration unit for your application


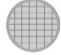




Select the funnel according to the sample type

	Aqueous samples	Antibiotics/solvents	Water for injection
			
Membrane type	0.45 µm MCE	0.45 µm PVDF	0.45 µm MCE
Catalog no. code	MMHA	MMHV	MMHA
Funnel volume	100 mL	100 mL	250 mL

MCE: mixed cellulose ester PVDF: Polyvinylidene fluoride - Durapore®

Summary of available filtration units

PVDF membranes have a yellow label,
0.45 µm MCE membranes have a white label.

	Membrane pore size	Membrane type	Membrane color	Funnel size	Catalog no.
White label	0.45 µm	MCE		100 mL	MMHAWG124
	0.45 µm	MCE		100 mL	MMHABG124
	0.45 µm	MCE		250 mL	MMHAWG224
Yellow label	0.45 µm	PVDF		100 mL	MMHVWP124
	0.45 µm	PVDF		250 mL	MMHVWP224
	0.45 µm	PVDF		100 mL	MMHVMFX24



Units are packed by 8 on 2 detachable trays

Each box contains 24 units: 3 bags of 8 units.