

**Low MWCO Filtration Plate**

MWCO: 100 KDa

Cat. No.	Amount
MP-MWCO-5-L	5 pcs.

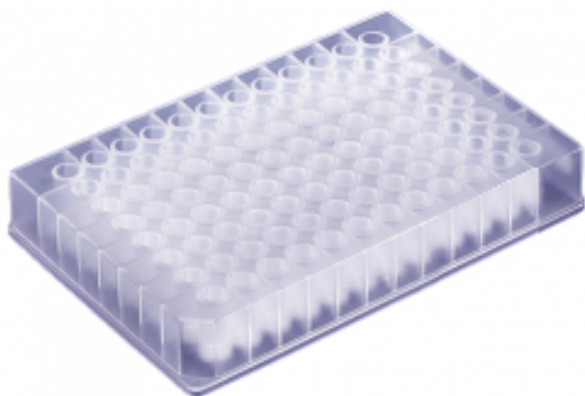
**Description:**

Plate Material: Polypropylene, hydrophobic

Number of Wells: 96

Well Capacity: 400 µl

Filter Membrane: Polyethersulfone (PES)

Molecular Weight Cutoff: 100 KDa

The plates are used to separate proteins/protein detergent complexes by their sizes (diameters). Polyethersulfone membranes have very low non-specific binding to proteins and thus optimize yields and purity.

For *in vitro* use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** n/a