



MRC Under Oil Crystallization Plate

material UVP

Cat. No.	Amount
CPL-152S	10 pcs.
CPL-152L	100 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a