



3 Well Crystallization Plate - Polystyrene

Cat. No.	Amount
CPL-150S	10 pcs.
CPL-150L	100 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a