



Vapor Batch Plate - hydrophilic

Cat. No.	Amount
CPL-129S	10 pcs.
CPL-129L	80 pcs.

For *in vitro* use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a