



Corning® Plate 3773

round bottom
advanced polymer, not treated

Cat. No.	Amount
CPL-126S	10 pcs.
CPL-126L	50 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a