



Corning® plate 3556

1 protein well – round bottom

Cat. No.	Amount
CPL-109S	10 pcs.
CPL-109L	50 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a