



## Corning® plate 3557

5 protein wells – conical flat-crystal basin

Cat. No.	Amount
CPL-115S	10 pcs.
CPL-115L	50 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a