



## 3 Well Crystallization Plate Low Profile - UVP

Cat. No.	Amount
CPL-164S	10 pcs.
CPL-164L	100 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a