



## AxyPearl™ Lid

Cat. No.	Amount
CPL-136S	20 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a