



## AxyGem™ Crystallography Plate

Cat. No.	Amount
CPL-135S	10 pcs.
CPL-135L	50 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a