



## Crystal Mount-Base-Assembly

MicroGripper Assembly with specified aperture size (xx) on 18mm/SPINE length rods adhered into specified barcoded bases (By) without cryovials

Cat. No.	Amount
HTB-M7-XX-BY	20 assemblies

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a