



## Crystal Mount-Base-Vial-Assembly

MicroMesh Assembly with specified aperture size (xx) on 18mm/SPINE length rods adhered into specified barcoded reusable bases (By) with cryovials

Cat. No.	Amount
HTA-M3-XX-BY-R	20 assemblies

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a