



Crystal Mount-Base-Assembly

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

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

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a