



Crystal Mount-Base-Vial-Assembly

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

Cat. No.	Amount
HTA-M5-XXLD-BY	20 assemblies

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a