



Crystal Mount-Base-Vial-Assembly

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

| Cat. No. | Amount |
|----------------|---------------|
| HTA-M8-XX-BY-R | 20 assemblies |

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a