



Crystal Mount-Base-Assembly

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

Cat. No.	Amount
B-M8-XX-BY	20 assemblies

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a