

**Crystal Mount-Base-Assembly**

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

| Cat. No.   | Amount        |
|------------|---------------|
| B-M7-XX-BY | 20 assemblies |

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a