



## Silicone Grease

high viscosity

Cat. No.	Amount
CGR-103	1 Tube (35 g)

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** 12 months