



Silicone Grease

medium viscosity

Cat. No.	Amount
CGR-101	1 Tube (35 g)

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 12 months