

**Apiezon N Cryogenic Vacuum Grease**

Cat. No.	Amount
APZN-1	25 g

nonsaleable to USA and Japan**Shipping:** shipped at ambient temperature**Storage Conditions:** store at ambient temperature**Shelf Life:** 12 months