



Magnetic CryoVial

Cat. No.	Amount
CV-50	50 pcs.
CV-100	100 pcs.
CV-200	200 pcs.

nonsaleable to USA and Japan

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a