



Greiner 780271 – Masterblock 2 ml

96 well, V-bottom
sterile, individually packed

Cat. No.	Amount
CMB-102	10 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a