



Greiner 381061 – Cap Mat for 1 ml Masterblock

96 well, not pierceable
sterile, individually packed

Cat. No.	Amount
CCM-101	10 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a