



Greiner CrystalBridge™ - 45 µl

for 24 well ComboPlates

Cat. No.	Amount
CPB-119	100 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a