



JBScreen Basic 1 - 4

Cat. No.	Amount
CS-125	4 Kits

Shipping: shipped at ambient temperature

Storage Conditions: store at 8-10 °C

Shelf Life: 12 months