



JBScreen Wizard 1 & 2

Cat. No.	Amount
CS-315	2 Kits

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 8-10 °C

Shelf Life: 12 months