



## Individual JBScreen Condition

Indicate screen name, Cat.# and condition # when placing the order

Cat. No.	Amount
CS-IND-10ML	10 ml
CS-IND-100ML	100 ml

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 8-10 °C

**Shelf Life:** 12 months