



HD Clear™ Sealing Tape - 3 inches wide

Cat. No.	Amount
CTP-101M	1 roll (50 m)

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a