



CrystalClear 'Duo'

Strips with two depressions
Box of 10 contains 10 frames with 12 strips each

Cat. No.	Amount
CPL-133	10 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a