



5' LC[®]Red640 / 3' BHQ-3[®]

Ex: 625 nm / Em: 640 nm

Cat. No.	Amount
FP-144-05	5 to 9 nmol
FP-144-10	10 to 19 nmol
FP-144-20	20 to 29 nmol
FP-144-30	30 to 49 nmol
FP-144-50	50 to 70 nmol

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months