

**5' LC<sup>®</sup>Red640 / 3' BBQ<sup>®</sup>-650**

Ex: 625 nm / Em: 640 nm

Cat. No.	Amount
FP-175-05	5 to 9 nmol
FP-175-10	10 to 19 nmol
FP-175-20	20 to 29 nmol
FP-175-30	30 to 49 nmol
FP-175-50	50 to 70 nmol

**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months