

**5' LC®Red610 / 3' ECLIPSE**

Ex: 590 nm / Em: 610 nm

Cat. No.	Amount
FP-160-05	5 to 9 nmol
FP-160-10	10 to 19 nmol
FP-160-20	20 to 29 nmol
FP-160-30	30 to 49 nmol
FP-160-50	50 to 70 nmol

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