

**5' Cy3 / 3' ECLIPSE**

Ex: 546 nm / Em: 563 nm

Cat. No.	Amount
FP-162-05	5 to 9 nmol
FP-162-10	10 to 19 nmol
FP-162-20	20 to 29 nmol
FP-162-30	30 to 49 nmol
FP-162-50	50 to 70 nmol

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