

**5' Cy5.5 / 3' BHQ-3[®]**

Ex: 683 nm / Em: 705 nm

Cat. No.	Amount
FP-142-05	5 to 9 nmol
FP-142-10	10 to 19 nmol
FP-142-20	20 to 29 nmol
FP-142-30	30 to 49 nmol
FP-142-50	50 to 70 nmol

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