

**5' HEX / 3' BHQ-1®**

Ex: 535 nm / Em: 556 nm

Cat. No.	Amount
FP-124-05	5 to 9 nmol
FP-124-10	10 to 19 nmol
FP-124-20	20 to 29 nmol
FP-124-30	30 to 49 nmol
FP-124-50	50 to 70 nmol

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