



**5' FITC / 3' BHQ-1®**

Ex: 490 nm / Em: 525 nm

Cat. No.	Amount
FP-185-05	5 to 9 nmol
FP-185-10	10 to 19 nmol
FP-185-20	20 to 29 nmol
FP-185-30	30 to 49 nmol
FP-185-50	50 to 70 nmol

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months