



## 5' DY-480XL / 3' BHQ-2<sup>®</sup>

Ex: 500 nm / Em: 630 nm

Cat. No.	Amount
FP-194-05	5 to 9 nmol
FP-194-10	10 to 19 nmol
FP-194-20	20 to 29 nmol
FP-194-30	30 to 49 nmol
FP-194-50	50 to 70 nmol

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

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

**Shelf Life:** 12 months