



## 5' 6-FAM / 3' ECLIPSE

Ex: 495 nm / Em: 520 nm

Cat. No.	Amount
FP-182-05	5 to 9 nmol
FP-182-10	10 to 19 nmol
FP-182-20	20 to 29 nmol
FP-182-30	30 to 49 nmol
FP-182-50	50 to 70 nmol

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

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

**Shelf Life:** 12 months