

**5' ATTO-390 / 3' TAMRA**

Ex: 390 nm / Em: 476 nm

Cat. No.	Amount
FP-420-05	5 to 9 nmol
FP-420-10	10 to 19 nmol
FP-420-20	20 to 29 nmol
FP-420-30	30 to 49 nmol
FP-420-50	50 to 70 nmol

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