

**5' ATTO-425 / 3' ECLIPSE**

Ex: 439 nm / Em: 485 nm

Cat. No.	Amount
FP-415-05	5 to 9 nmol
FP-415-10	10 to 19 nmol
FP-415-20	20 to 29 nmol
FP-415-30	30 to 49 nmol
FP-415-50	50 to 70 nmol

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