

**5' ATTO-495 / 3' ECLIPSE**

Ex: 498 nm / Em: 526 nm

Cat. No.	Amount
FP-419-05	5 to 9 nmol
FP-419-10	10 to 19 nmol
FP-419-20	20 to 29 nmol
FP-419-30	30 to 49 nmol
FP-419-50	50 to 70 nmol

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