

**5' ATTO-Rho14 / 3' BHQ-2®**

Ex: 625 nm / Em: 646 nm

Cat. No.	Amount
FP-199-05	5 to 9 nmol
FP-199-10	10 to 19 nmol
FP-199-20	20 to 29 nmol
FP-199-30	30 to 49 nmol
FP-199-50	50 to 70 nmol

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