



## 5' ATTO-Rho13 / 3' BHQ-3®

Ex: 603 nm / Em: 627 nm

Cat. No.	Amount
FP-197-05	5 to 9 nmol
FP-197-10	10 to 19 nmol
FP-197-20	20 to 29 nmol
FP-197-30	30 to 49 nmol
FP-197-50	50 to 70 nmol

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

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

**Shelf Life:** 12 months