

**5' ATTO-Rho6G / 3' BHQ-2®**

Ex: 533 nm / Em: 557 nm

| Cat. No. | Amount |
|-----------|---------------|
| FP-405-05 | 5 to 9 nmol |
| FP-405-10 | 10 to 19 nmol |
| FP-405-20 | 20 to 29 nmol |
| FP-405-30 | 30 to 49 nmol |
| FP-405-50 | 50 to 70 nmol |

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