



5' ATTO-Rho13 / 3' BHQ-2®

Ex: 603 nm / Em: 627 nm

| Cat. No. | Amount |
|-----------|---------------|
| FP-196-05 | 5 to 9 nmol |
| FP-196-10 | 10 to 19 nmol |
| FP-196-20 | 20 to 29 nmol |
| FP-196-30 | 30 to 49 nmol |
| FP-196-50 | 50 to 70 nmol |

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months