

**5' TAMRA / 3' BHQ-2[®]**

Ex: 564 nm / Em: 579 nm

| Cat. No. | Amount |
|-----------|---------------|
| FP-126-05 | 5 to 9 nmol |
| FP-126-10 | 10 to 19 nmol |
| FP-126-20 | 20 to 29 nmol |
| FP-126-30 | 30 to 49 nmol |
| FP-126-50 | 50 to 70 nmol |

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