

**5' TET / 3' ECLIPSE**

Ex: 521 nm / Em: 536 nm

| Cat. No.  | Amount        |
|-----------|---------------|
| FP-153-05 | 5 to 9 nmol   |
| FP-153-10 | 10 to 19 nmol |
| FP-153-20 | 20 to 29 nmol |
| FP-153-30 | 30 to 49 nmol |
| FP-153-50 | 50 to 70 nmol |

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