

**5' JOE / 3' DABCYL**

Ex: 522 nm / Em: 548 nm

| Cat. No.  | Amount        |
|-----------|---------------|
| FP-114-05 | 5 to 9 nmol   |
| FP-114-10 | 10 to 19 nmol |
| FP-114-20 | 20 to 29 nmol |
| FP-114-30 | 30 to 49 nmol |
| FP-114-50 | 50 to 70 nmol |

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