



## 5' Cy3 / 3' BBQ®-650

Ex: 546 nm / Em: 563 nm

| Cat. No.  | Amount        |
|-----------|---------------|
| FP-170-05 | 5 to 9 nmol   |
| FP-170-10 | 10 to 19 nmol |
| FP-170-20 | 20 to 29 nmol |
| FP-170-30 | 30 to 49 nmol |
| FP-170-50 | 50 to 70 nmol |

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months