



## 5' Cy5.5 / 3' BBQ®-650

Ex: 683 nm / Em: 705 nm

Cat. No.	Amount
FP-177-05	5 to 9 nmol
FP-177-10	10 to 19 nmol
FP-177-20	20 to 29 nmol
FP-177-30	30 to 49 nmol
FP-177-50	50 to 70 nmol

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

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

**Shelf Life:** 12 months