

**5' 6-FAM / 3' DABCYL**

Ex: 495 nm / Em: 520 nm

Cat. No.	Amount
FP-183-05	5 to 9 nmol
FP-183-10	10 to 19 nmol
FP-183-20	20 to 29 nmol
FP-183-30	30 to 49 nmol
FP-183-50	50 to 70 nmol

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