



5' FITC / 3' TAMRA

Ex: 490 nm / Em: 525 nm

Cat. No.	Amount
FP-184-05	5 to 9 nmol
FP-184-10	10 to 19 nmol
FP-184-20	20 to 29 nmol
FP-184-30	30 to 49 nmol
FP-184-50	50 to 70 nmol

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months