



5' Fluo / 3' BHQ-1®

Ex: 495 nm / Em: 520 nm

Cat. No.	Amount
FP-190-05	5 to 9 nmol
FP-190-10	10 to 19 nmol
FP-190-20	20 to 29 nmol
FP-190-30	30 to 49 nmol
FP-190-50	50 to 70 nmol

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months