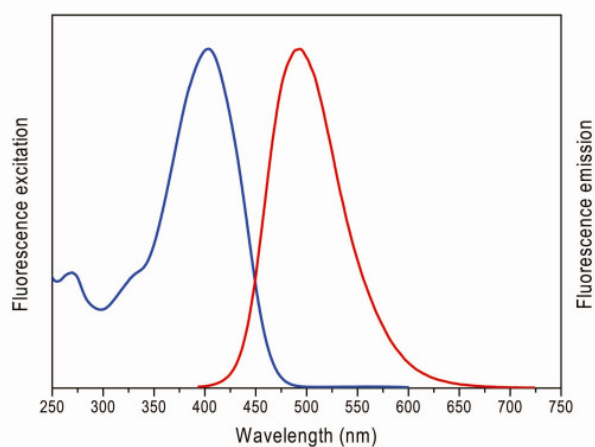


Eterneon™-Azide (384 / 480) green fluorescent dye

Cat. No.	Amount
CLK-FA012-1	1 mg
CLK-FA012-5	5 mg
CLK-FA012-10	10 mg



Excitation and emission spectrum of Eterneon (384 / 480)

Molecular weight: 538.73 g/mol

Spectroscopic properties:

λ_{abs} 384 nm; λ_{em} 480 nm (in DMSO);
 ϵ 18000 cm⁻¹ M⁻¹ (in DCM)

Quantum yield Φ : 42%

Appearance: orange solid

Purity: > 97 %

Solubility: DMF, DCM, DMSO,
Water/Tween® 20 (0.5 %)

Storage conditions: store at 4 °C

Shelf life: 12 months

For research use only!