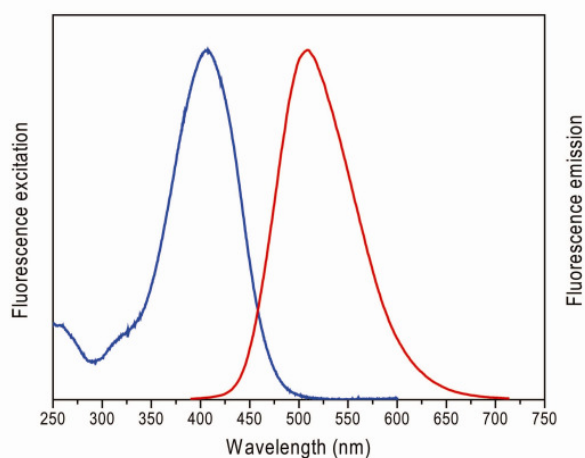


Eterneon™-Azide (394 / 507) green-yellow fluorescent dye

Cat. No.	Amount
CLK-FA014-1	1 mg
CLK-FA014-5	5 mg
CLK-FA014-10	10 mg



Excitation and emission spectrum of Eterneon (394 / 507)

Molecular weight: 538.73 g/mol

Spectroscopic properties:

λ_{abs} 394 nm; λ_{em} 507 nm (in DMSO);
 ϵ 30000 cm⁻¹ M⁻¹ (in DCM)

Quantum yield Φ : 37%

Appearance: yellow-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!