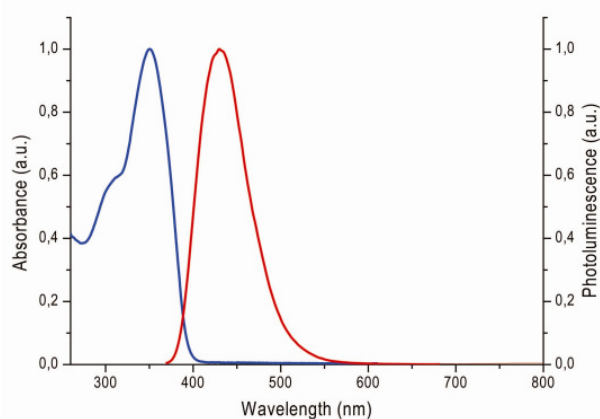


Eterneon™-Azide (350 / 430) blue fluorescent dye

Cat. No.	Amount
CLK-FA010-1	1 mg
CLK-FA010-5	5 mg
CLK-FA010-10	10 mg



Excitation and emission spectrum of Eterneon (350 / 430)

Molecular weight: 456.60 g/mol

Spectroscopic properties:

λ_{abs} 350 nm; λ_{em} 430 nm (in DMSO);
 ϵ 15220 cm⁻¹ M⁻¹ (in DCM)

Quantum yield Φ : 70%

Appearance: yellow 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!