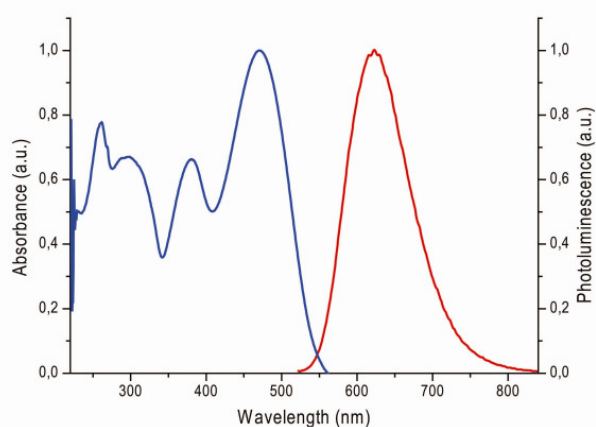


## Eterneon™-Azide (480 / 635) orange-red fluorescent dye

Cat. No.	Amount
CLK-FA015-1	1 mg
CLK-FA015-5	5 mg
CLK-FA015-10	10 mg



Excitation and emission spectrum of Eterneon (480 / 635)

**Molecular weight:** 626.79 g/mol

**Spectroscopic properties:**

$\lambda_{\text{abs}}$  480 nm;  $\lambda_{\text{em}}$  635 nm (in DMSO);  
 $\epsilon$  29900 cm<sup>-1</sup> M<sup>-1</sup> (in DCM)

**Quantum yield  $\Phi$ :** 34%

**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!**