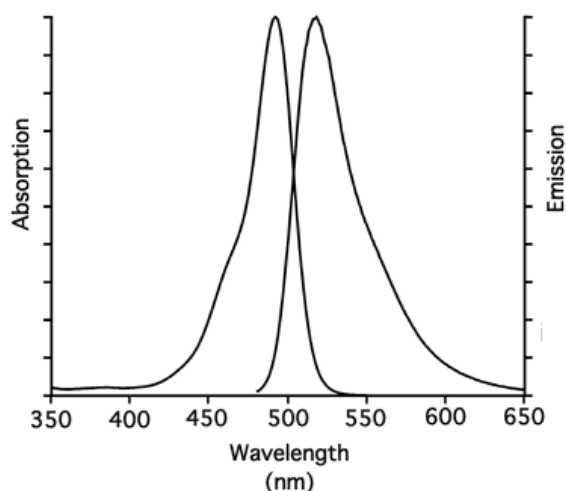
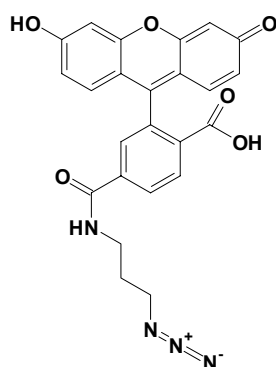


Fluorescein-Azide (6-FAM Azide)

N-(3-Azido-propyl)-2-(6-hydroxy-3-oxo-3H-xanthen-9-yl)-terephthalamic acid

Cat. No.	Amount
CLK-FA001-1	1 mg
CLK-FA001-5	5 mg
CLK-FA001-100	100 mg



Absorption and emission spectrum of 6-FAM

Molecular formula: C₂₄H₁₈N₄O₆

Molecular weight: 458.43 g/mol

Spectroscopic properties:

λ_{abs} 492 nm; λ_{em} 517 nm; ϵ 83000 cm⁻¹M⁻¹

Storage conditions: -20 °C, dry, dark

Purity: > 95 %

Appearance: orange solid

Solubility: DMSO, DMF, MeOH

Shelf life: 12 months

For research use only!

Selected References:

Gramlich *et al.* (2008) Click-Click-Click: Single to Triple Modification of DNA. *Angew. Chem. Int. Ed.* **47**:3442.

Gramlich *et al.* (2008) Postsynthetic DNA Modification through the Copper-Catalyzed Azide-Alkyne Cycloaddition Reaction. *Angew. Chem. Int. Ed.* **47**:8350.

Zhan *et al.* (2005) Synthesis of hemicyanine dyes for 'click' bioconjugation. *Tetrahedron Letters* **46**:1691.