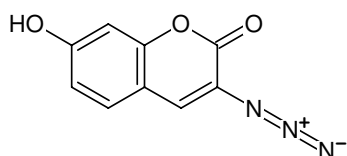


**3-Azido-7-hydroxycoumarin**

Abs/Em = 404/477 nm

3-Azido-7-hydroxy-chromen-2-one

Cat. No.	Amount
CLK-FA047-1	1 mg
CLK-FA047-5	5 mg



Structural formula of 3-Azido-7-hydroxycoumarin

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₉H₅N₃O₃**Molecular Weight:** 203.15 g/mol**Exact Mass:** 203.03 g/mol**Purity:** ≥ 90 % (HPLC)**Form:** solid**Color:** brown**Solubility:** DMF, DMSO, MeCN, MeOH (Remark: solution in DMSO and DMF are not suitable for longterm storage)**Spectroscopic Properties:** λ_{abs} 348 nm (dye before click reaction in 100 mM Tris-HCl pH 7.5), λ_{abs} 404 nm (after click reaction with L-Homopropargylglycine in 100 mM potassium phosphate buffer pH 7.0), λ_{em} 477 nm (after click reaction with L-Homopropargylglycine in 100 mM potassium phosphate buffer pH 7.0)**Selected References:**Laughlin *et al.* (2009) Imaging the glycome. *Proc Natl Acad Sci USA*. **106** (1):12.Li *et al.* (2010) Fluorogenic click reaction for labeling and detection of DNA in proliferating cells. *Biotechniques*. **49** (1):525.Sivakumar *et al.* (2004) A fluorogenic 1,3-dipolar cycloaddition reaction of 3-azidocoumarins and acetylenes. *Org Lett*. **6** (24):4603.