

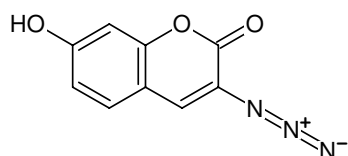


3-Azido-7-hydroxycoumarin

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 research use only!

Shipping: shipped on blue ice

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: > 95 % (HPLC, NMR)

Form: brown solid

Solubility: DMF, DMSO, MeCN, MeOH (Remark: solution in DMSO and DMF are not suitable for longterm storage)

Spectroscopic Properties: λ_{abs} 404 nm (after click reaction), λ_{em} 477 nm (after click reaction)

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.