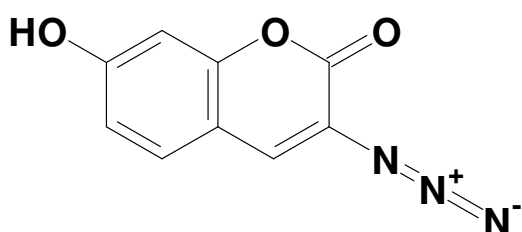


## 3-Azido-7-hydroxycoumarin

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

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



**Molecular Formula:** C<sub>9</sub>H<sub>5</sub>N<sub>3</sub>O<sub>3</sub>

**Molecular Weight:** 203.15 g/mol

**Storage conditions:** store at -20°C and dark

**Purity:** > 95 % (HPLC & NMR)

**Appearance:** brown solid

**Solubility:** DMSO, DMF, MeOH, MeCN

Remark: solution in DMSO and DMF are not suitable for longterm storage

**Shelf life:** 12 months

**Spectroscopic properties:**

after click reaction  $\lambda_{\text{obs}}$  404 nm and  $\lambda_{\text{em}}$  477 nm

**For research use only!**

**This product is not available in USA.**

#### 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.