

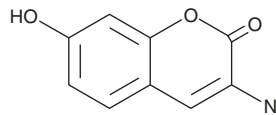
# PRODUCT INFORMATION



## 3-azido-7-hydroxy Coumarin

Item No. 10596

**CAS Registry No.:** 817638-68-9  
**Formal Name:** 3-azido-7-hydroxy-2H-1-benzopyran-2-one  
**Synonyms:** 3-azido-7-Hydroxycoumarin, AHC, Azidohydroxycoumarin  
**MF:** C<sub>9</sub>H<sub>5</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 203.2  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Ex./Em. Max:** 404/477 nm



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

3-azido-7-hydroxy Coumarin is supplied as a solid. A stock solution may be made by dissolving the 3-azido-7-hydroxy coumarin in the solvent of choice, which should be purged with an inert gas. 3-azido-7-hydroxy Coumarin is soluble in methanol, acetonitrile and DMSO.

### Description

3-azido-7-hydroxy Coumarin is a click-reactive fluorescent dye.<sup>1,2</sup> Upon copper(I)-catalyzed azide/alkyne cycloaddition (CuAAC) in a biological system, a fluorescent product is produced that displays excitation/emission maxima of 404 and 477 nm, respectively. 3-azido-7-hydroxy Coumarin has been used for spatially restricted metabolite labeling and the detection of foodborne pathogens.

### References

1. You, H., Wang, M., Wang, S., *et al.* Ultrasensitive and specific phage@DNAzyme probe-triggered fluorescent click chemistry for on-site detection of foodborne pathogens using a smartphone. *Anal. Chem.* **95(30)**, 11211-11218 (2023).
2. Mishra, P.K., Sharma, N., Kim, H., *et al.* Gen-Click: Genetically encodable click reactions for spatially restricted metabolite labeling. *ACS Cent. Sci.* **9(8)**, 1650-1657 (2023).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/01/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)