

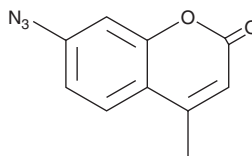
PRODUCT INFORMATION



7-Azido-4-methylcoumarin

Item No. 28307

CAS Registry No.: 95633-27-5
Formal Name: 7-azido-4-methyl-2H-1-benzopyran-2-one
Synonym: AzMC
MF: $C_{10}H_7N_3O_2$
FW: 201.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 297, 324 nm
Ex./Em. Max.: 340/445 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

7-Azido-4-methylcoumarin (AzMC) is supplied as a crystalline solid, which should be purged with an inert gas. A stock solution may be made by dissolving the AzMC in the solvent of choice. AzMC is soluble in the organic solvent DMSO.

Description

AzMC is a coumarin fluorescent probe for hydrogen sulfide (H_2S).¹ AzMC selectively fluoresces in the presence of H_2S over a panel of reactive nitrogen, oxygen, and sulfur species. AzMC displays excitation/emission maxima of 340/445 nm, respectively.

Reference

1. Chen, B., Li, W., Lv, C., *et al.* Fluorescent probe for highly selective and sensitive detection of hydrogen sulfide in living cells and cardiac tissues. *Analyst* **138**(3), 946-951 (2013).

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, 12/02/2022

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