

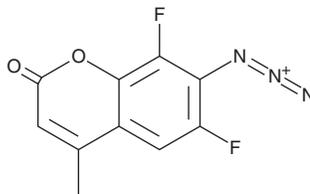
PRODUCT INFORMATION



Difluorinated H₂S Fluorescent Probe 1

Item No. 29835

CAS Registry No.: 2103919-91-9
Formal Name: 7-azido-6,8-difluoro-4-methyl-2H-1-benzopyran-2-one
MF: C₁₀H₅F₂N₃O₂
FW: 237.2
Purity: ≥98%
UV/Vis.: λ_{max}: 296, 326 nm
Ex./Em. Max: 365/450 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Difluorinated H₂S probe 1 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice.

Description

Difluorinated H₂S probe 1 is a fluorescent probe for hydrogen sulfide (H₂S).¹ It selectively fluoresces in the presence of H₂S over Zn²⁺, Fe³⁺, S₂O₃²⁻, ClO⁻, SO₃²⁻, H₂O₂, NO₂⁻, cysteine (Cys), homocysteine (Hcy), and glutathione (GSH) when used at a concentration of 1 μM. Difluorinated H₂S probe 1 displays excitation/emission maxima of 365/450 nm, respectively.

Reference

1. Zhang, J., Gao, Y., Kang, X., *et al.* *o,o*-Difluorination of aromatic azide yields a fast-response fluorescent probe for H₂S detection and for improved bioorthogonal reactions. *Org. Biomol. Chem.* **15**(19), 4212-4217 (2017).

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, 01/30/2024

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