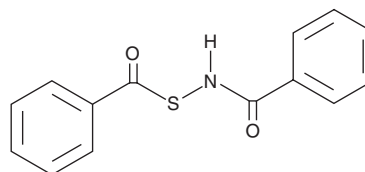


PRODUCT INFORMATION



H₂S Donor 5a Item No. 11238

CAS Registry No.: 134861-13-5
Formal Name: benzenecarbothioic acid, S-(benzoylazanyl) ester
Synonym: N-(Benzoylthio)benzamide
MF: C₁₄H₁₁NO₂S
FW: 257.3
Purity: ≥98%
Supplied as: A 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

H₂S Donor 5a is supplied as a solid. A stock solution may be made by dissolving the H₂S donor 5a in the solvent of choice. H₂S Donor 5a is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

H₂S Donor 5a is a stable, cysteine-activated H₂S donor. In the presence of excess cysteine, the concentration of H₂S released from this compound reaches a maximum value at 18 min.¹ It does not react with potential cellular nucleophiles such as -OH and -NH₂ groups.¹

Reference

1. Zhao, Y., Wang, H., and Xian, M. Cysteine-activated hydrogen sulfide (H₂S) donors. *J. Am. Chem. Soc.* **133**, 15-17 (2011).

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, 11/14/2022

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