

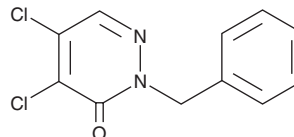
PRODUCT INFORMATION



EcDsbB-IN-9

Item No. 44749

CAS Registry No.: 41933-33-9
Synonyms: EcDsbB Inhibitor 9, *E. coli* DsbB-IN-9,
E. coli DsbB Inhibitor 9
Formal Name: 4,5-dichloro-2-(phenylmethyl)-3(2H)-pyridazinone
MF: C₁₁H₈Cl₂N₂O
FW: 255.1
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

EcDsbB-IN-9 is supplied as a solid. A stock solution may be made by dissolving the EcDsbB-IN-9 in the solvent of choice, which should be purged with an inert gas. EcDsbB-IN-9 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

EcDsbB-IN-9 is an inhibitor of the disulfide bond formation protein DsbB (IC₅₀ = 1.7 μM for the *E. coli* enzyme).¹ It is active against *E. coli* (MIC = 5.7 μM). EcDsbB-IN-9 is also active against the nematodes *B. xylophilus*, *D. destructor*, and *A. besseyi* when used at a concentration of 100 μg/ml.²

References

1. Landeta, C., Blazyk, J.L., Hatahet, F., *et al.* Compounds targeting disulfide bond forming enzyme DsbB of Gram-negative bacteria. *Nat. Chem. Biol.* **11(4)**, 292-298 (2015).
2. Lu, S., Zhu, Z., Du, T., *et al.* Nematicidal activity and preliminary mechanism of pyridazine compounds. *Phys. Mol. Plant Pathol.* **139**, 102836 (2025).

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, 01/26/2026

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