

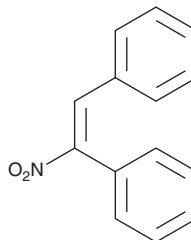
PRODUCT INFORMATION



CID-2818500

Item No. 21919

CAS Registry No.: 1215-07-2
Formal Name: 1,1'-(1-nitro-1,2-ethenediyl)bis-benzene
Synonyms: α -Nitrostilbene, NSC 385
MF: C₁₄H₁₁NO₂
FW: 225.2
Purity: \geq 98%
UV/Vis.: λ_{max} : 228, 316 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

CID-2818500 is supplied as a crystalline solid. A stock solution may be made by dissolving the CID-2818500 in the solvent of choice, which should be purged with an inert gas. CID-2818500 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CID-2818500 in these solvents is approximately 12, 20, and 25 mg/ml, respectively.

CID-2818500 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CID-2818500 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CID-2818500 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CID-2818500 is an inhibitor of protein arginine methyltransferase 1 (PRMT1; IC₅₀ = 11 μ M in a histone H4 methylation assay).¹ It also inhibits histone H4 methylation by PRMT8, but not histone H3.1 methylation by CARM1 or Set7/9, when used at concentrations of 10 and 100 μ M.

Reference

1. Dillon, M.B., Bachovchin, D.A., Brown, S.J., *et al.* Novel inhibitors for PRMT1 discovered by high-throughput screening using activity-based fluorescence polarization. *ACS Chem. Biol.* **7**(7), 1198-1204 (2012).

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