

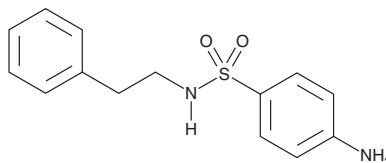
PRODUCT INFORMATION



C-7280948

Item No. 21601

CAS Registry No.: 587850-67-7
Formal Name: 4-amino-N-(2-phenylethyl)-benzenesulfonamide
MF: C₁₄H₁₆N₂O₂S
FW: 276.4
Purity: ≥98%
UV/Vis.: λ_{max}: 265 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

C-7280948 is supplied as a crystalline solid. A stock solution may be made by dissolving the C-7280948 in the solvent of choice, which should be purged with an inert gas. C-7280948 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of C-7280948 in these solvents is approximately 1, 20, and 25 mg/ml, respectively.

C-7280948 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, C-7280948 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. C-7280948 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

C-7280948 is a protein arginine methyltransferase 1 (PRMT1) inhibitor (IC₅₀ = 12.8 μM, *in vitro*).^{1,2} PRMT1 enzymes transfer a methyl group from S-adenosylmethionine to arginine residues on histones, which leads to transcription.² PRMT1 is essential for mixed lineage leukemia oncogenic transformation and thus its inhibition is a potential treatment for these aggressive leukemias.³

References

1. Luo, M. Current chemical biology approaches to interrogate protein methyltransferases. *ACS Chem Biol.* **7**(3), 443-463 (2012).
2. Heinke, R., Spannhoff, A., Meier, R., *et al.* Virtual screening and biological characterization of novel histone arginine methyltransferase PRMT1 inhibitors. *Chem. Med. Chem.* **4**(1), 69-77 (2009).
3. Muntean, A.G. and Hess, J.L. The pathogenesis of mixed-lineage leukemia. *Annu. Rev. Pathol.* **7**, 283-301 (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