

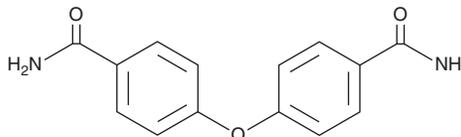
PRODUCT INFORMATION



OUL35

Item No. 28845

CAS Registry No.: 6336-34-1
Formal Name: 4,4'-oxybis-benzamide
Synonym: NSC 39047
MF: C₁₄H₁₂N₂O₃
FW: 256.3
Purity: ≥95%
UV/Vis.: λ_{max}: 258 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

OUL35 is supplied as a crystalline solid. A stock solution may be made by dissolving the OUL35 in the solvent of choice, which should be purged with an inert gas. OUL35 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of OUL35 in these solvents is approximately 30 and 5 mg/ml, respectively.

OUL35 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, OUL35 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. OUL35 has a solubility of approximately 0.20 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

OUL35 is an inhibitor of poly(ADP-ribose) polymerase 10 (PARP10; IC₅₀ = 329 nM).¹ It is selective for PARP10 over a panel of 15 PARP enzymes, including PARP4, PARP14, and PARP16 (IC₅₀s = 22.6, 23.4, and 4.17 μM, respectively). OUL35 (10 μM) reverses inhibition of HeLa cell proliferation induced by PARP10 overexpression. It sensitizes HeLa cells to hydroxyurea (Item No. 23725), increasing hydroxyurea-induced inhibition of cell proliferation when used at a concentration of 5 μM.

Reference

1. Venkannagari, H., Verheugd, P., Koivunen, J., *et al.* Small-molecule chemical probe rescues cells from mono-ADP-ribosyltransferase ARTD10/PARP10-induced apoptosis and sensitizes cancer cells to DNA damage. *Cell Chem. Bio.* **23(10)**, 1251-1260 (2016).

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, 12/08/2022

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