

PRODUCT INFORMATION



3,5-Bis(4-nitrophenoxy)benzoic Acid

Item No. 36553

CAS Registry No.: 173550-33-9

Synonym: Compound W

MF: $C_{19}H_{12}N_2O_8$

FW: 396.3

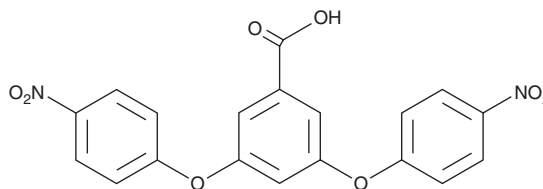
Purity: $\geq 98\%$

UV/Vis.: $\lambda_{\max}$: 299 nm

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

3,5-Bis(4-nitrophenoxy)benzoic acid is supplied as a solid. A stock solution may be made by dissolving the 3,5-bis(4-nitrophenoxy)benzoic acid in the solvent of choice, which should be purged with an inert gas. 3,5-Bis(4-nitrophenoxy)benzoic acid is soluble in organic solvents such as ethanol and DMSO. The solubility of 3,5-bis(4-nitrophenoxy)benzoic acid in these solvents is approximately 100 mM and 5 mM, respectively.

Description

3,5-Bis(4-nitrophenoxy)benzoic acid is a Notch1 signaling modulator.^{1,2} It decreases the production of C-terminally elongated Notch1 amyloid- β -like peptide (N β 25) from the full-length receptor, a marker of Notch1 intracellular domain release and transcription modification, in HEK293 cells.¹ 3,5-Bis(4-nitrophenoxy)benzoic acid (50 μM) also decreases activation of the Notch1 target hes family bHLH transcription factor 1 (HES1) in a luciferase reporter assay using HCT116 cells.²

References

1. Okochi, M., Fukumori, A., Jiang, J., *et al.* Secretion of the Notch-1 A β -like peptide during Notch signaling. *J. Biol. Chem.* **281**(12), 7890-7898 (2006).
2. Sail, V. and Hadden, M.K. Identification of small molecule Hes1 modulators as potential anticancer chemotherapeutics. *Chem. Biol. Drug Des.* **81**(3), 334-342 (2013).

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, 12/20/2022

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