

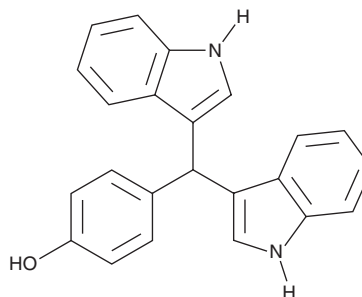
PRODUCT INFORMATION



C-DIM8

Item No. 30694

CAS Registry No.: 151358-47-3
Formal Name: 4-(di-1H-indol-3-ylmethyl)-phenol
Synonyms: DIM-C-*p*PhOH, 1,1-bis-(3'-indolyl)-1-(*p*-hydroxyphenyl)methane
MF: C₂₃H₁₈N₂O
FW: 338.4
Purity: ≥95%
UV/Vis.: λ_{max}: 212, 262 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

C-DIM8 is supplied as a solid. A stock solution may be made by dissolving the C-DIM8 in the solvent of choice, which should be purged with an inert gas. C-DIM8 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of C-DIM8 in ethanol and DMSO is approximately 10 mg/ml and approximately 15 mg/ml in DMF.

C-DIM8 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, C-DIM8 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. C-DIM8 has a solubility of approximately 0.14 mg/ml in a 1:6 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-DIM8 is a *para*-phenyl-substituted diindolylmethane (C-DIM) and an antagonist of the orphan receptor nuclear receptor-related protein 77 (Nur77).¹ It inhibits reporter gene expression induced by the Nur77 agonist C-DIM5 (Item No. 27980) in PANC-28 cells when used at a concentration of 20 μM. C-DIM8 (10, 15, and 20 μM) reduces expression of *BIRC5*, the gene encoding survivin, in PANC-1 pancreatic cancer cells.² *In vivo*, C-DIM8 (30 mg/kg) induces tumor cell apoptosis and reduces tumor volume in an L3.6pl pancreatic carcinoma mouse xenograft model.

References

- Chintharlapalli, S., Burghardt, R., Papineni, S., *et al.* Activation of Nur77 by selected 1,1-Bis(3'-indolyl)-1-(*p*-substituted phenyl)methanes induces apoptosis through nuclear pathways. *J. Biol. Chem.* **280**(26), 24903-24914 (2005).
- Lee, S.-O., Abdelrahim, M., Yoon, K., *et al.* Inactivation of the orphan nuclear receptor TR3/Nur77 inhibits pancreatic cancer cell and tumor growth. *Cancer Res.* **70**(17), 6824-6836 (2010).

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, 12/12/2022

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