

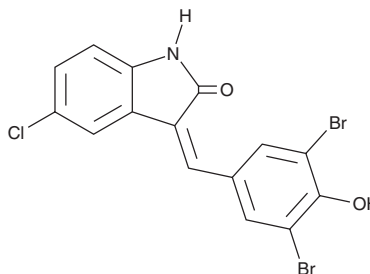
PRODUCT INFORMATION



CID-25014542

Item No. 29837

CAS Registry No.: 220904-99-4
Formal Name: 5-chloro-3-[(3,5-dibromo-4-hydroxyphenyl)methylene]-1,3-dihydro-2H-indol-2-one
MF: C₁₅H₈Br₂ClNO₂
FW: 429.5
Purity: ≥95%
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

CID-25014542 is supplied as a solid. A stock solution may be made by dissolving the CID-25014542 in the solvent of choice, which should be purged with an inert gas. CID-25014542 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

CID-25014542 is an inhibitor of leucine-rich repeat kinase 2 (LRRK2; IC₅₀ = 15 nM).¹ It is selective for LRRK2 over a panel of 44 other kinases but does inhibit MAPK4 and Pim-2 at 1 μM. CID-25014542 (10 μM) also inhibits FMS-related tyrosine kinase 3 (FLT3). It decreases the production of amyloid-β (1-42) (Aβ42) in a reporter assay using H4 cells overexpressing Aβ42 (IC₅₀ = 15 μM). CID-25014542 (20 μM) inhibits EGF-induced activation of MAPK in serum-starved cells.² It increases the survival of primary rat granule neurons in low potassium media when used at concentrations of 1, 5, or 25 μM.³

References

1. Göring, S., Taymans, J.M., Baekelandt, V., *et al.* Indolinone based LRRK2 kinase inhibitors with a key hydrogen bond. *Bioorg. Med. Chem. Lett.* **24(19)**, 4630-4637 (2014).
2. Wood, E., Crosby, R.M., Dickerson, S., *et al.* A prodrug approach to the design of cRaf1 kinase inhibitors with improved cellular activity. *Anticancer Drug Des.* **16(1)**, 1-6 (2001).
3. Balderamos, M., Ankati, H., Akubathini, S.K., *et al.* Synthesis and structure-activity relationship studies of 3-substituted indolin-2-ones as effective neuroprotective agents. *Exp. Biol. Med. (Maywood)* **233(11)**, 1395-1402 (2008).

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, 05/21/2025

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