

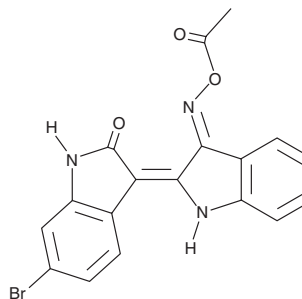
PRODUCT INFORMATION



BIO-Acetoxime

Item No. 16329

CAS Registry No.: 740841-15-0
Formal Name: 3-[3-[(acetyloxy)imino]-1,3-dihydro-2H-indol-2-ylidene]-6-bromo-1,3-dihydro-2H-indol-2-one
Synonyms: 6-Bromoindirubin-3'-acetoxime, GSK3 Inhibitor X
MF: C₁₈H₁₂BrN₃O₃
FW: 398.2
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 248, 290, 325, 523 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BIO-Acetoxime is supplied as a crystalline solid. A stock solution may be made by dissolving the BIO-acetoxime in the solvent of choice. BIO-Acetoxime is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of BIO-acetoxime in these solvents is approximately 10 mg/ml.

Description

BIO-Acetoxime is an analog of the GSK3 inhibitor 6-bromoindirubin-3'-oxime, or BIO (Item No. 13123).^{1,2} It potently inhibits GSK3α/β (IC₅₀ = 10 nM), with selectivity over Cdk5/p25, Cdk2/A, and Cdk1/B (IC₅₀s = 2.4, 4.3, and 63 μM, respectively).^{1,2} BIO-Acetoxime has been used to study the role of GSK3 in the Wnt/β-catenin pathway.³ It also suppresses the cytopathic effects of herpes viruses and reduces viral yields in human oral epithelial cells.⁴

References

1. Meijer, L., Skaltsounis, A.-L., Magiatis, P., *et al.* GSK-3-selective inhibitors derived from tyrian purple indirubins. *Chem. Biol.* **10**, 1255-1266 (2003).
2. Polychronopoulos, P., Magiatis, P., Skaltsounis, A.-L., *et al.* Structural basis for the synthesis of indirubins as potent and selective inhibitors of glycogen synthase kinase-3 and cyclin-dependent kinases. *J. Med. Chem.* **47**, 935-946 (2004).
3. Wang, B., Liu, J., Ma, L.N., *et al.* Chimeric 5/35 adenovirus-mediated Dickkopf-1 overexpression suppressed tumorigenicity of CD44⁺ gastric cancer cells *via* attenuating Wnt signaling. *J. Gastroenterol.* **48**, 798-808 (2013).
4. Hsu, M.J. and Hung, S.L. Antiherpetic potential of 6-bromoindirubin-3'-acetoxime (BIO-acetoxime) in human oral epithelial cells. *Arch. Virol.* **158**, 1287-1296 (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, 06/16/2020

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