

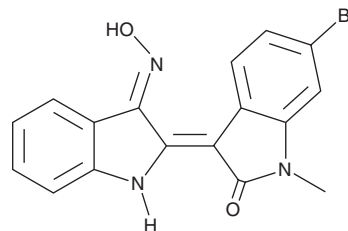
PRODUCT INFORMATION



MeBIO

Item No. 19573

CAS Registry No.: 667463-95-8
Formal Name: (3Z)-6-bromo-3-[(3E)-1,3-dihydro-3-(hydroxyimino)-2H-indol-2-ylidene]-1,3-dihydro-1-methyl-2H-indol-2-one
Synonyms: Me6BIO, 1-Methyl-6-bromo-indirubin-3'-oxime
MF: C₁₇H₁₂BrN₃O₂
FW: 370.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Special Conditions: Keep cool and dry. Protect from light and moisture.

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

Laboratory Procedures

MeBIO is supplied as a solid. A stock solution may be made by dissolving the MeBIO in the solvent of choice. MeBIO is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas.

Description

MeBIO is a control analog of the GSK3 α/β inhibitor, BIO (Item No. 13123). It displays minimal activity against Cdk1/Cyclin B, GSK3 α/β , and Cdk5/p25 (IC₅₀s = 92.0, 44-100, and >100 μ M, respectively).^{1,2} MeBIO has been identified as an aryl hydrocarbon receptor (AhR) agonist (EC₅₀s = 20 and 93 nM for yeast and a hepatoma cell line, respectively) that causes AhR to translocate to the nucleus.³

References

1. 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(4)**, 935-946 (2004).
2. Meijer, L., Skaltsounis, A.L., Magiatis, P., *et al.* GSK-3-selective inhibitors derived from tyrian purple indirubins. *Chem. Biol.* **10(12)**, 1255-1266 (2003).
3. Knockaert, M., Blondel, M., Bach, S., *et al.* Independent actions on cyclin-dependent kinases and aryl hydrocarbon receptor mediate the antiproliferative effects of indirubins. *Oncogene* **23(25)**, 4400-4412 (2004).

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, 12/22/2022

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