

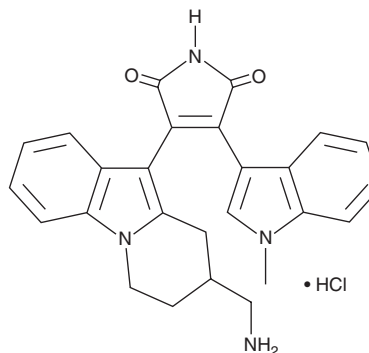
PRODUCT INFORMATION



Bisindolymaleimide X (hydrochloride)

Item No. 17511

CAS Registry No.: 145317-11-9
Formal Name: 3-[8-(aminomethyl)-6,7,8,9-tetrahydropyrido[1,2-a]indol-10-yl]-4-(1-methyl-1H-indol-3-yl)-1H-pyrrole-2,5-dione, monohydrochloride
Synonyms: BIM X, Ro 31-8425
MF: C₂₆H₂₄N₄O₂ • HCl
FW: 461.0
Purity: ≥98%
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

Bisindolymaleimide X (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the bisindolymaleimide X (hydrochloride) in the solvent of choice. Bisindolymaleimide X (hydrochloride) is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

Bisindolymaleimide X is a cell-permeable, reversible, ATP-competitive protein kinase C (PKC) inhibitor ($IC_{50} = 15$ nM, rat brain PKC).¹ It has been shown to inhibit PKC α , β_1 , β_{11} , γ , and ϵ with IC_{50} values of 8, 8, 14, 13, and 39 nM, respectively.¹ It has also been identified as an inhibitor of Cdk2 ($IC_{50} = 200$ nM) and GSK3 α/β .² This compound has been used to activate mesenchymal stem cells, increase the surface expression of homing ligands that bind to intercellular adhesion molecule 1, and target delivery of these cells to sites of inflammation.³

References

1. Wilkinson, S.E., Parker, P.J., and Nixon, J.S. Isoenzyme specificity of bisindolymaleimides, selective inhibitors of protein kinase C. *Biochem. J.* **294**(Pt 2), 335-337 (1993).
2. Brehmer, D., Godl, K., Zech, B., *et al.* Proteome-wide identification of cellular targets affected by bisindolymaleimide-type protein kinase C inhibitors. *Mol. Cell. Proteomics* **3**(5), 490-500 (2004).
3. Levy, O., Mortensen, L.J., Boquet, G., *et al.* A small-molecule screen for enhanced homing of systemically infused cells. *Cell Rep.* **10**(8), 1261-1268 (2015).

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, 12/02/2022

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