

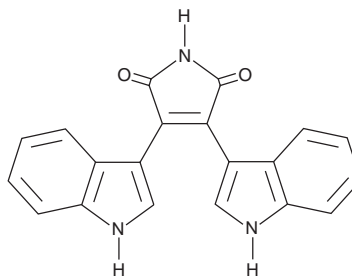
PRODUCT INFORMATION



Bisindolymaleimide IV

Item No. 13299

CAS Registry No.: 119139-23-0
Formal Name: 3,4-di-1H-indol-3-yl-1H-pyrrole-2,5-dione
Synonyms: Arcyriarubin A, BIM IV
MF: C₂₀H₁₃N₃O₂
FW: 327.3
Purity: ≥98%
UV/Vis.: λ_{max}: 276, 373, 461 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

Bisindolymaleimide IV is supplied as a crystalline solid. A stock solution may be made by dissolving the bisindolymaleimide IV in the solvent of choice, which should be purged with an inert gas. Bisindolymaleimide IV is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Bisindolymaleimide IV in these solvents is approximately 20 mg/ml.

Bisindolymaleimide IV is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, bisindolymaleimide IV should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Bisindolymaleimide IV has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Bisindolymaleimide IV (BIM IV) is a cell permeable inhibitor of protein kinase C (PKC) with IC₅₀ values reported to range from 0.10 to 0.55 μM.¹⁻⁴ BIM IV was designed to be more discriminative than its parent compound, the non-selective PKC inhibitor, staurosporine (Item No. 81590). However, BIM IV also inhibits protein kinase A with IC₅₀ values ranging from 2 to 11.8 μM.^{1,2}

References

1. Davis, P.D., Hill, C.H., Lawton, G., *et al.* Inhibitors of protein kinase C. 1.1 2,3-bisarylmalimides. *J. Med. Chem.* **35**(1), 177-184 (1992).
2. Toullec, D., Pianetti, P., Coste, H., *et al.* The bisindolymaleimide GF 109203X is a potent and selective inhibitor of protein kinase C. *J. Biol. Chem.* **266**(24), 15771-15781 (1991).
3. Fabre, S. and Prudhomme, M. Protein kinase C inhibitors; structure-activity relationships in K252c-related compounds. *Bioorg. Med. Chem.* **1**(3), 193-196 (1993).
4. Lazareno, S., Popham, A., and Birdsall, N.J.M. Muscarinic interactions of bisindolymaleimide analogues. *Eur. J. Pharmacol.* **360**(2-3), 281-284 (1998).

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, 04/24/2023

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