

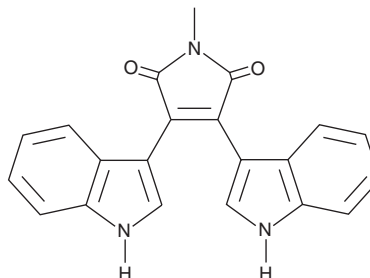
PRODUCT INFORMATION



Bisindolymaleimide V

Item No. 13300

CAS Registry No.: 113963-68-1
Formal Name: 3,4-di-1H-indol-3-yl-1-methyl-1H-pyrrole-2,5-dione
Synonyms: BIM V, Ro 31-6045
MF: C₂₁H₁₅N₃O₂
FW: 341.4
Purity: ≥98%
UV/Vis.: λ_{max}: 278, 375, 471 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 V is supplied as a crystalline solid. A stock solution may be made by dissolving the bisindolymaleimide V in the solvent of choice, which should be purged with an inert gas. Bisindolymaleimide V is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of bisindolymaleimide V in these solvents is approximately 20 mg/ml.

Bisindolymaleimide V is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, bisindolymaleimide V should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Bisindolymaleimide V has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Bisindolymaleimide V (BIM V) is a weak inhibitor of protein kinase C (PKC) demonstrating an IC₅₀ value >100 μM.¹⁻³ While effectively inactive as a PKC inhibitor, BIM V blocks the activation of mitogen-stimulated protein kinase p70s6k/p85s6k (S6K) *in vivo* with an IC₅₀ value of 8 μM.⁴

References

1. Lazareno, S., Popham, A., and Birdsall, N.J.M. Muscarinic interactions of bisindolymaleimide analogues. *Eur. J. Pharmacol.* **360**(2-3), 281-284 (1998).
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. 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).
4. Marmy-Conus, N., Hannan, K.M., and Pearson, R.B. Ro 31-6045, the inactive analogue of the protein kinase C inhibitor Ro 31-8220, blocks *in vivo* activation of p70s6k/p85s6k: Implications for the analysis of S6K signalling. *FEBS Lett.* **519**(1-3), 135-140 (2002).

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, 11/02/2022

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