

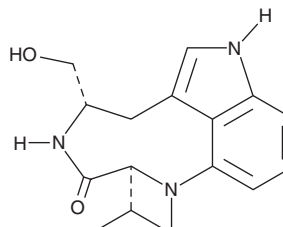
PRODUCT INFORMATION



(-)-Indolactam V

Item No. 14647

CAS Registry No.: 90365-57-4
Formal Name: (2S,5S)-1,2,4,5,6,8-hexahydro-5-(hydroxymethyl)-1-methyl-2-(1-methylethyl)-3H-pyrrolo[4,3,2-gh]-1,4-benzodiazonin-3-one
MF: C₁₇H₂₃N₃O₂
FW: 301.4
Purity: ≥95%
UV/Vis.: λ_{max}: 203, 229, 301 nm
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

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

Description

(-)-Indolactam V is a tumor-promoting teleocidin compound that activates protein kinase C (PKC). It binds to PKC regulatory domains of mouse skin PKC η and rat brain PKC γ with K_i values of 3.4 nM and 1 μ M, respectively.^{1,2} At 1 μ M, (-)-indolactam V induces differentiation of human embryonic stem cells into pancreatic progenitors through activation of PKC signaling.³

References

1. Nakagawa, Y., Irie, K., Nakamura, Y., *et al.* Synthesis and biological activities of indolactone-V, the lactone analogue of the tumor promoter (-)-indolactam-V. *Biosci. Biotechnol. Biochem.* **61(8)**, 1415-1417 (1997).
2. Masuda, A., Irie, K., Nakagawa, Y., *et al.* Binding selectivity of conformationally restricted analogues of (-)-indolactam-V to the C1 domains of protein kinase C isozymes. *Biosci. Biotechnol. Biochem.* **66(7)**, 1615-1617 (2002).
3. Chen, S., Borowiak, M., Fox, J.L., *et al.* A small molecule that directs differentiation of human ESCs into the pancreatic lineage. *Nat. Chem. Biol.* **5(4)**, 258-265 (2009).

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, 11/23/2022

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