

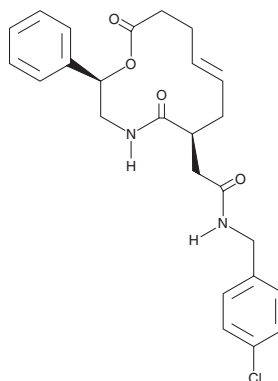
PRODUCT INFORMATION



Robotnikinin

Item No. 13204

CAS Registry No.: 1132653-79-2
Formal Name: N-[(4-chlorophenyl)methyl]-5,12-dioxo-2R-phenyl-1-oxa-4-azacyclododec-8E-ene-6S-acetamide
MF: C₂₅H₂₇ClN₂O₄
FW: 455.0
Purity: ≥95%
Supplied as: A crystalline 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

Robotnikinin is supplied as a crystalline solid. A stock solution may be made by dissolving the robotnikinin in the solvent of choice, which should be purged with an inert gas. Robotnikinin is soluble in ethanol, DMSO, and dimethyl formamide.

Description

Sonic hedgehog (Shh) is the most widely characterized of the three vertebrate Hedgehog homologs, and is essential for proper embryonic development.¹ Shh binds to its receptor, Patched (Ptch1), resulting in the de-repression of Smoothened.^{2,3} This leads to the activation of Gli2, which regulates the transcription of target genes that include Gli1 and Ptch1. Robotnikinin is the first identified small molecule inhibitor of Shh signaling that acts upstream of Smo, binding directly to Shh with a K_d value of 3.1 μM.^{4,5} Robotnikinin represses Gli-mediated transcription in primary human keratinocytes and synthetic human skin in a concentration-dependent manner.⁴

References

1. Heretsch, P., Tzagkaroulaki, L., and Giannis, A. Cyclopamine and hedgehog signaling: Chemistry, biology, medical perspectives. *Angew. Chem. Int. Ed.* **49(20)**, 3418-3427 (2010).
2. Stone, D.M., Hynes, M., Armanini, M., *et al.* The tumour-suppressor gene patched encodes a candidate receptor for sonic hedgehog. *Nature* **384(6605)**, 129-133 (1996).
3. Rohatgi, R., Milenkovic, L., Corcoran, R.B., *et al.* Hedgehog signal transduction by smoothened: Pharmacologic evidence for a 2-step activation process. *Proc. Natl. Acad. Sci. USA* **106(9)**, 3196-3201 (2009).
4. Stanton, B.Z., Peng, L.F., Maloof, N., *et al.* A small molecule that binds hedgehog and blocks its signaling in human cells. *Nat. Chem. Biol.* **5(3)**, 154-156 (2009).
5. Stanton, B.Z. and Peng, L.F. Small-molecule modulators of the Sonic Hedgehog signaling pathway. *Mol. BioSyst.* **6(1)**, 44-54 (2010).

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, 11/03/2022

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