

PRODUCT INFORMATION



EMD 534085

Item No. 35719

CAS Registry No.: 858668-07-2

Formal Name: N-[2-(dimethylamino)ethyl]-N'-
[[[(2R,4aS,5R,10bS)-3,4,4a,5,6,10b-
hexahydro-5-phenyl-9-
(trifluoromethyl)-2H-pyrano[3,2-c]
quinolin-2-yl]methyl]-urea

MF: C₂₅H₃₁F₃N₄O₂

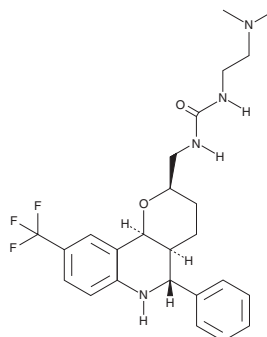
FW: 476.5

Purity: ≥99%

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

EMD 534085 is supplied as a solid. A stock solution may be made by dissolving the EMD 534085 in the solvent of choice, which should be purged with an inert gas. EMD 534085 is soluble in DMSO.

Description

EMD 534085 is an inhibitor of the kinesin spindle protein Eg5 (IC₅₀ = 8 nM).¹ It inhibits proliferation of HCT116 colon cancer cells (IC₅₀ = 30 nM), as well as induces mitotic arrest and apoptosis in HL-60 leukemia cells.^{1,2} EMD 534085 (15 and 30 mg/kg) reduces tumor growth in a COLO 205 colon cancer mouse xenograft model.¹

References

1. Schiemann, K., Finsinger, D., Zenke, F., *et al.* The discovery and optimization of hexahydro-2H-pyrano[3,2-c]quinolines (HHPQs) as potent and selective inhibitors of the mitotic kinesin-5. *Bioorg. Med. Chem. Lett.* **20**(5), 1491-1495 (2010).
2. Tang, Y., Orth, J.D., Xie, T., *et al.* Rapid induction of apoptosis during Kinesin-5 inhibitor-induced mitotic arrest in HL60 cells. *Cancer Lett.* **310**(1), 15-24 (2011).

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, 12/13/2022

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