

PRODUCT INFORMATION

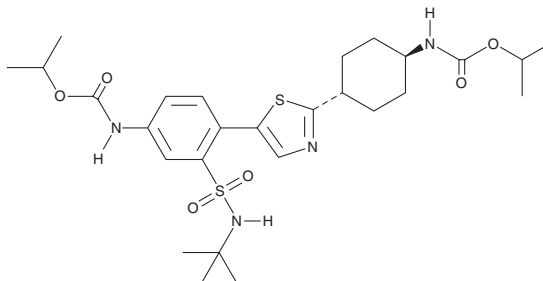


Emzadirib

Item No. 44841

CAS Registry No.: 2301085-04-9
Formal Name: N-[3-[[[(1,1-dimethylethyl)amino]sulfonyl]-4-[2-[trans-4-[[[(1-methylethoxy)carbonyl]amino]cyclohexyl]-5-thiazolyl]phenyl]-, carbamic acid, 1-methylethyl ester

Synonym: CYT0851
MF: C₂₇H₄₀N₄O₆S₂
FW: 580.8
Purity: ≥95%
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

Emzadirib is supplied as a solid. A stock solution may be made by dissolving the emzadirib in the solvent of choice, which should be purged with an inert gas. Emzadirib is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Emzadirib is an inhibitor of RAD51 with anticancer activity *in vitro* and in animal models.¹

Reference

1. Day, M., Maclay, T., Cyr, A., *et al.* Targeting homologous recombination in lymphoid malignancies: Evaluation of four small molecule inhibitors of RAD51. *Blood* **134**(Supp. 1), 2080 (2019).

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, 01/19/2026

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