

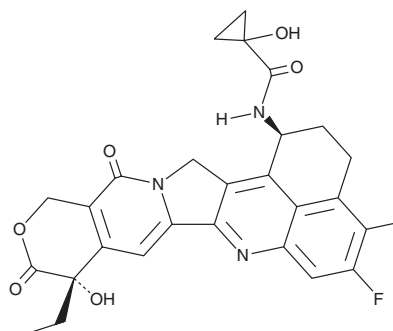
# PRODUCT INFORMATION



## Exatecan-amide-cyclopropanol

Item No. 44470

**CAS Registry No.:** 2414254-34-3  
**Formal Name:** N-[(1S,9S)-9-ethyl-5-fluoro-2,3,9,10,13,15-hexahydro-9-hydroxy-4-methyl-10,13-dioxo-1H,12H-benzo[de]pyrano[3',4':6,7]indolizino[1,2-b]quinolin-1-yl]-1-hydroxy-cyclopropanecarboxamide  
**MF:** C<sub>28</sub>H<sub>26</sub>FN<sub>3</sub>O<sub>6</sub>  
**FW:** 519.5  
**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

Exatecan-amide-cyclopropanol is supplied as a solid. A stock solution may be made by dissolving the exatecan-amide-cyclopropanol in the solvent of choice, which should be purged with an inert gas. Exatecan-amide-cyclopropanol is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

### Description

Exatecan-amide-cyclopropanol is a derivative of the DNA topoisomerase I inhibitor exatecan (Item No. 35452) that has been used as a building block in the synthesis of antibody-drug conjugates (ADCs) with anticancer activity.<sup>1</sup>

### Reference

1. Xu, J., Zhang, Y., Cai, X., *et al.* Ligand-drug conjugate of exatecan analogue, preparation method thereof and application thereof. *Jiangsu Hengrui Medicine Co. and Ltd. Shanghai Hengrui Pharmaceutical Co., Ltd.* US2021/0353764A1 (2021).

#### 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, 11/05/2025

#### 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