

PRODUCT INFORMATION

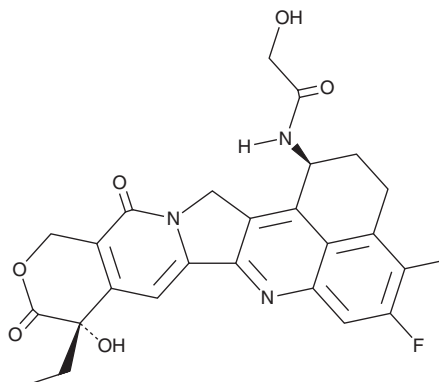


Dxd

Item No. 41699

CAS Registry No.: 1599440-33-1
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]-2-hydroxy-acetamide

MF: $C_{26}H_{24}FN_3O_6$
FW: 493.5
Purity: $\geq 95\%$
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

Dxd is supplied as a solid. A stock solution may be made by dissolving the Dxd in the solvent of choice, which should be purged with an inert gas. Dxd is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Dxd is a derivative of the DNA topoisomerase I inhibitor exatecan (Item No. 35452) and a starting material that has been used in the synthesis of antibody-drug conjugates (ADCs).¹ It has been used in the synthesis of ADCs targeting HER2 with anticancer activity *in vitro* and *in vivo*.

Reference

1. Conilh, L., Fournet, G., Fourmaux, E., *et al.* Exatecan antibody drug conjugates based on a hydrophilic polysarcosine drug-linker platform. *Pharmaceuticals (Basel)* **14**(3), 247 (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, 09/04/2024

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