

PRODUCT INFORMATION



CL2A-SN-38 (dichloroacetic acid salt)

Item No. 33941

Formal Name: carbonic acid, [4-[[[(2S)-2-(4-aminobutyl)-35-[4-[[[4-[[[2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)methyl]cyclohexyl]carbonyl]amino]methyl]-1H-1,2,3-triazol-1-yl]-1,4,8-trioxo-6,12,15,18,21,24,27,30,33-nonaoxa-3,9-diazapentatriacont-1-yl]amino]phenyl)methyl (4S)-4,11-diethyl-3,4,12,14-tetrahydro-9-hydroxy-3,14-dioxo-1H-pyrano[3',4':6,7]indolizino[1,2-b]quinolin-4-yl ester, dichloroacetic acid

MF: C₇₃H₉₇N₁₁O₂₂ • XC₂H₂Cl₂O₂

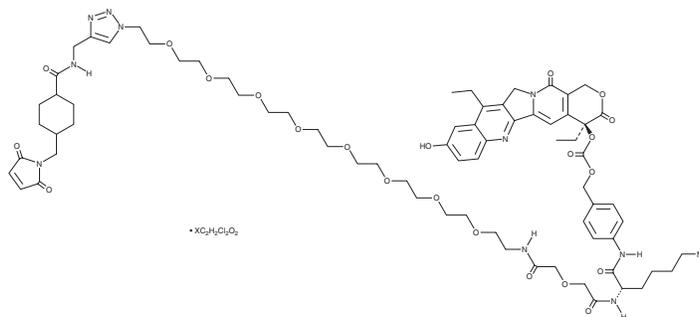
FW: 1,480.6

Purity: ≥90%

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

CL2A-SN-38 (dichloroacetic acid salt) is supplied as a solid. A stock solution may be made by dissolving the CL2A-SN-38 (dichloroacetic acid salt) in the solvent of choice, which should be purged with an inert gas. CL2A-SN-38 (dichloroacetic acid salt) is soluble in the organic solvent methanol at a concentration of approximately 10 mg/ml.

Description

CL2A-SN-38 is a cleavable linker-based antibody-drug conjugate (ADC) containing the topoisomerase I inhibitor SN-38 (Item No. 15632) and a maleimidocaproyl linker.¹ CL2A-SN-38 has been linked to tumor-selective humanized antibodies, such as anti-Trop-2 (hRS7), to target SN-38 to tumor sites. CL2A-SN-38 linked to hRS7 is cytotoxic to Calu-3, COLO 205, Capan-1, PC3, SK-MES-1, and BxPC-3 cancer cells (IC₅₀s = 1.95-23.14 nM). CL2A-SN-38 linked to hRS7 reduces tumor growth in BxPC-3 pancreatic and SK-MES-1 lung cancer mouse xenograft models.

Reference

- Cardillo, T.M., Govindan, S.V., Sharkey, R.M., *et al.* Humanized anti-Trop-2 IgG-SN-38 conjugate for effective treatment of diverse epithelial cancers: Preclinical studies in human cancer xenograft models and monkeys. *Clin. Cancer Res.* **17(10)**, 3157-3169 (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, 10/04/2022

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