

PRODUCT INFORMATION



LCL-161, phenol

Item No. 40358

CAS Registry No.: 2095244-42-9

Formal Name: N-[(1S)-2-[[[(1S)-1-cyclohexyl-2-[(2S)-2-[4-(3-hydroxybenzoyl)-2-thiazolyl]-1-pyrrolidinyl]-2-oxoethyl]amino]-1-methyl-2-oxoethyl]-N-methyl-carbamic acid, 1,1-dimethylethyl ester

Synonym: cIAP1 Ligand 1

MF: $C_{31}H_{42}N_4O_6S$

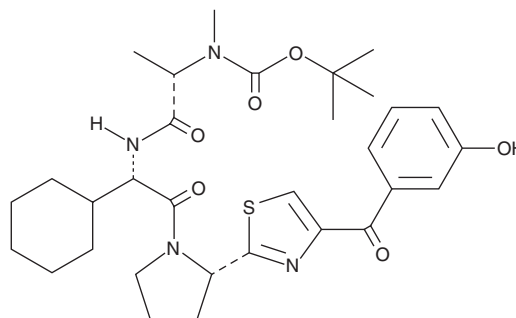
FW: 598.8

Purity: $\geq 98\%$

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

LCL-161, phenol is supplied as a solid. A stock solution may be made by dissolving the LCL-161, phenol in the solvent of choice, which should be purged with an inert gas. LCL-161, phenol is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

LCL-161, phenol is a *tert*-butyloxycarbonyl derivative of the inhibitor of cellular inhibitor of apoptosis (cIAP) family proteins LCL-161 (Item No. 22420) and an intermediate in the synthesis of proteolysis-targeted chimeras (PROTACs).¹ It has been used in the synthesis of PROTACs that drive degradation of androgen receptors.

Reference

1. Shibata, N., Nagai, K., Morita, Y., *et al.* Development of protein degradation inducers of androgen receptor by conjugation of androgen receptor ligands and inhibitor of apoptosis protein ligands. *J. Med. Chem.* **61**(2), 543-575 (2018).

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, 10/01/2024

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