

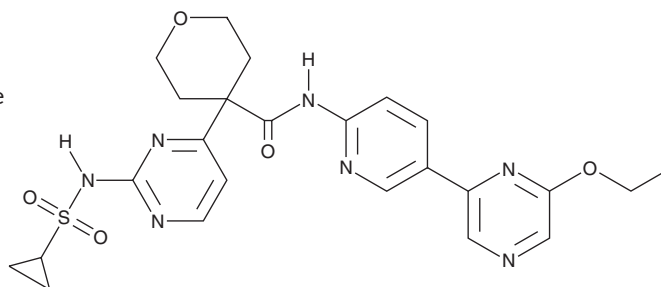
# PRODUCT INFORMATION



## Dencatistat

Item No. 45363

**CAS Registry No.:** 2377000-84-3  
**Formal Name:** 4-[2-[(cyclopropylsulfonyl)amino]-4pyrimidinyl]-N-[5-(6-ethoxy-2-pyrazinyl)-2-pyridinyl]tetrahydro-2H-pyran-4-carboxamide  
**Synonym:** STP938  
**MF:** C<sub>24</sub>H<sub>27</sub>N<sub>7</sub>O<sub>5</sub>S  
**FW:** 525.6  
**Purity:** ≥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

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

### Description

Dencatistat is an inhibitor of CTP synthase 1 (CTPS1; IC<sub>50</sub> = 0.0093 μM), the enzyme that catalyzes the rate-limiting step in the conversion of UTP to CTP.<sup>1</sup> It is selective for CTPS1 over CTPS2 (IC<sub>50</sub> = 12.62 μM). Dencatistat inhibits the proliferation of nine cell lines derived from T cell-based cancers (IC<sub>50</sub>s = 2-183 nM) and 12 cell lines derived from B cell-based cancers (IC<sub>50</sub>s = 3-356 nM). It also inhibits the proliferation of JeKo-1 mantle cell lymphoma and DoHH2 B cell lymphoma cells and induces apoptosis in MOLT-4, Jurkat, and DoHH2 cells. Dencatistat (6-50 mg/kg every other day) reduces tumor growth in a JeKo-1 mouse xenograft model.

### Reference

1. Asnagli, H., Minet, N., Pfeiffer, C., *et al.* CTP synthase 1 is a novel therapeutic target in lymphoma. *Hemasphere* 7(4), e864 (2023).

#### 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, 04/20/2026

#### 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