

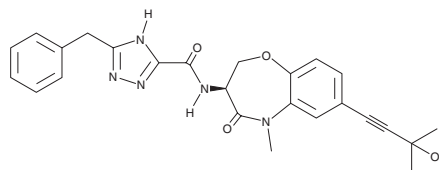
PRODUCT INFORMATION



Ocadusertib

Item No. 46372

CAS Registry No.: 2382811-41-6
Formal Name: 3-(phenylmethyl)-N-[(3S)-2,3,4,5-tetrahydro-7-(3-hydroxy-3-methyl-1-butyne-1-yl)-5-methyl-4-oxo-1,5-benzoxazepin-3-yl]-1H-1,2,4-triazole-5-carboxamide
Synonyms: LY3871801, R552
MF: C₂₅H₂₅N₅O₄
FW: 459.5
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

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

Description

Ocadusertib is an inhibitor of receptor-interacting serine/threonine kinase 1 (RIPK1; IC₅₀ = 0.038 μM).¹ It is selective for RIPK1 over RIPK2 and RIPK3 (IC₅₀s = >20 μM for both) and a panel of 105 kinases at 10 μM. Ocadusertib inhibits necroptosis induced by TNF and zVAD in U937 cells (EC₅₀ = 3 nM) and apoptosis induced by TNF and (5Z)-7-oxozeaenol (Item No. 17459) in primary human T cells (EC₅₀ = 0.4 nM). It inhibits hypothermia and decreases serum cytokine levels in a mouse model of sterile sepsis at 5 mg/kg. Ocadusertib also decreases disease severity in the Sharpincpdm mouse model of chronic proliferative dermatitis.

Reference

1. Shaw, S.J., Taylor, V.C., Sims, J.T., *et al.* Early development of ocadusertib, a selective receptor-interacting serine/threonine-protein kinase 1 inhibitor. *Clin. Transl. Sci.* **19**(3), (2026).

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, 08/26/2026

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