

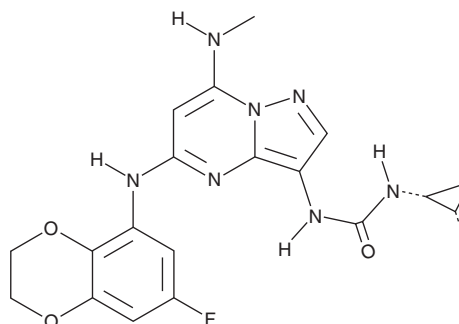
PRODUCT INFORMATION



Cenacitinib

Item No. 44648

CAS Registry No.: 2641636-52-2
Formal Name: N-[(1R,2S)-2-fluorocyclopropyl]-N'-[5-[(7-fluoro-2,3-dihydro-1,4-benzodioxin-5-yl)amino]-7-(methylamino)pyrazolo[1,5-a]pyrimidin-3-yl]-urea
MF: C₁₉H₁₉F₂N₇O₃
FW: 431.4
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

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

Description

Cenacitinib is an inhibitor of tyrosine kinase 2 (TYK2).¹

Reference

1. Bollu, V., Nuss, J., Yuan, S., *et al.* Preparation of a TYK2 inhibitor cross-reference. WO 2024/151992 A1 (2024).

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, 12/04/2025

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