

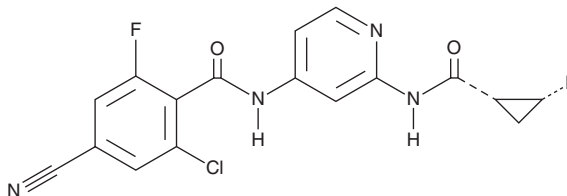
PRODUCT INFORMATION



TC JL 37

Item No. 43110

CAS Registry No.: 1258294-34-6
Formal Name: 2-chloro-4-cyano-6-fluoro-N-[2-
[[[(1R,2R)-2-fluorocyclopropyl]carbonyl]
amino]-4-pyridinyl]-benzamide
MF: C₁₇H₁₁ClF₂N₄O₂
FW: 376.7
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

TC JL 37 is supplied as a solid. A stock solution may be made by dissolving the TC JL 37 in the solvent of choice, which should be purged with an inert gas. TC JL 37 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

A TYK2 inhibitor ($K_i = 1.6$ nM); inhibits IL-12-induced phosphorylation of STAT4 in cells and IL-12- and IL-18-induced IFN- γ production in isolated human PBMCs (EC_{50} s = 224 and 168 nM, respectively); inhibits IL-12- and IL-18-induced increases in the serum levels of IFN- γ in mice at 10 to 100 mg/kg

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, 04/10/2025

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