

# PRODUCT INFORMATION

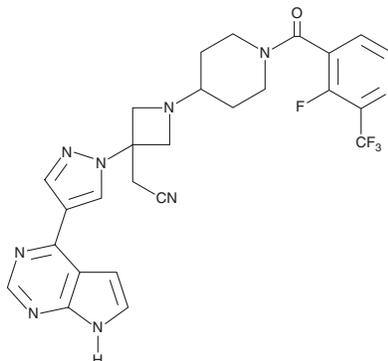


## Itacitinib

Item No. 27597

**CAS Registry No.:** 1334298-90-6  
**Formal Name:** 1-[1-[[3-fluoro-2-(trifluoromethyl)-4-pyridinyl]carbonyl]-4-piperidinyl]-3-[4-(7H-pyrrolo[2,3-d]pyrimidin-4-yl)-1H-pyrazol-1-yl]-3-azetidineaetonitrile

**Synonym:** INCB 039110  
**MF:** C<sub>26</sub>H<sub>23</sub>F<sub>4</sub>N<sub>9</sub>O  
**FW:** 553.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 255, 311 nm  
**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

Itacitinib is supplied as a solid. A stock solution may be made by dissolving the itacitinib in the solvent of choice, which should be purged with an inert gas. Itacitinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of itacitinib in these solvents is approximately 5, 2.5, and 3 mg/ml, respectively.

Itacitinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, itacitinib should first be dissolved in ethanol or DMF and then diluted with the aqueous buffer of choice. Itacitinib has a solubility of approximately 0.02 mg/ml in a 1:40 solution of ethanol:PBS (pH 7.2) or DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Itacitinib is an inhibitor of JAK1.<sup>1</sup> It is greater than 20- and 200-fold selective for JAK1 over JAK2 and JAK3, respectively.

### Reference

1. Norman, P. Selective JAK inhibitors in development for rheumatoid arthritis. *Expert Opin. Investig. Drugs* **23**(8), 1067-1077 (2014).

#### 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, 12/15/2022

#### 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