

# PRODUCT INFORMATION

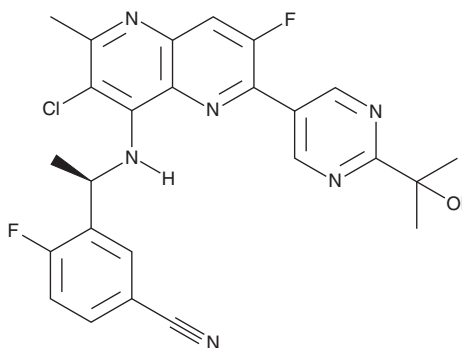


## TNF- $\alpha$ -IN-2

Item No. 45397

**CAS Registry No.:** 2074702-04-6  
**Formal Name:** 3-[(1R)-1-[[3-chloro-7-fluoro-6-[2-(1-hydroxy-1-methyl ethyl)-5-pyrimidinyl]-2-methyl-1,5-naphthyridin-4-yl]amino]ethyl]-4-fluoro-benzonitrile

**Synonym:** TNF-Inhibitor 2  
**MF:** C<sub>25</sub>H<sub>21</sub>ClF<sub>2</sub>N<sub>6</sub>O  
**FW:** 494.9  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

TNF- $\alpha$ -IN-2 is supplied as a solid. A stock solution may be made by dissolving the TNF- $\alpha$ -IN-2 in the solvent of choice, which should be purged with an inert gas. TNF- $\alpha$ -IN-2 is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

TNF- $\alpha$ -IN-2 is an inhibitor of TNF- $\alpha$ .<sup>1</sup> It inhibits TNF- $\alpha$ -induced secretion of a secreted placental alkaline phosphatase (SEAP) reporter in HEK-Blue cells (IC<sub>50</sub> = 27 nM). TNF- $\alpha$ -IN-2 (10 mg/kg) reduces paw TNF- $\alpha$ , IL-1 $\beta$ , and IL-6 levels and swelling in a mouse model of  $\alpha$ -collagen antibody-induced arthritis.

### Reference

1. Xiao, H. Y., Li, N., Duan, J. J., et al. Biologic-like in vivo efficacy with small molecule inhibitors of TNF $\alpha$  identified using scaffold hopping and structure-based drug design approaches. *J. Med. Chem.* **63**(23), 15050-15071 (2020).

#### 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, 05/04/2026

#### 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