

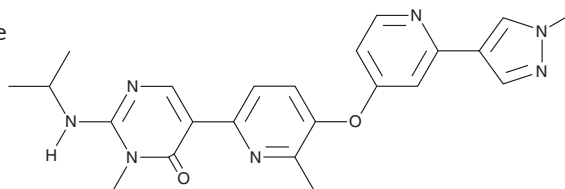
# PRODUCT INFORMATION



## Vimseltinib

Item No. 38783

**CAS Registry No.:** 1628606-05-2  
**Formal Name:** 3-methyl-2-[(1-methylethyl)amino]-5-[6-methyl-5-[[2-(1-methyl-1H-pyrazol-4-yl)-4-pyridinyl]oxy]-2-pyridinyl]-4(3H)-pyrimidinone  
**Synonym:** DCC-3014  
**MF:** C<sub>23</sub>H<sub>25</sub>N<sub>7</sub>O<sub>2</sub>  
**FW:** 431.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 264, 335 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

Vimseltinib is supplied as a solid. A stock solution may be made by dissolving the vimseltinib in the solvent of choice, which should be purged with an inert gas. Vimseltinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of vimseltinib in these solvents is approximately 10 mg/ml.

### Description

Vimseltinib is an inhibitor of colony stimulating factor 1 receptor (CSF1R; IC<sub>50</sub> = 3.7 nM).<sup>1</sup> It is selective for CSF1R over FMS-related tyrosine kinase 3 (FLT3), PDGFRα, PDGFRβ, and KIT (IC<sub>50</sub>s = >3,300, 436, 2,300, and 476 nM, respectively). Vimseltinib inhibits the proliferation of M-NFS-60 leukemia cells (IC<sub>50</sub> = 18 nM). It inhibits CSF1-induced CSF1R autophosphorylation in THP-1 cells (IC<sub>50</sub> = 27 nM) and CSF1- and RANKL-induced osteoclast differentiation (IC<sub>50</sub> = 9.3 nM).

### Reference

1. Caldwell, T.M., Ahn, Y.M., Bulfer, S.L., *et al.* Discovery of vimseltinib (DCC-3014), a highly selective CSF1R switch-control kinase inhibitor, in clinical development for the treatment of Tenosynovial Giant Cell Tumor (TGCT). *Bioorg. Med. Chem. Lett.* **74**, 128928 (2022).

#### 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, 08/09/2023

#### 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