

# PRODUCT INFORMATION



## Tolebrutinib

Item No. 37723

**CAS Registry No.:** 1971920-73-6  
**Formal Name:** 4-amino-1,3-dihydro-1-[(3R)-1-(1-oxo-2-propen-1-yl)-3-piperidinyl]-3-(4-phenoxyphenyl)-2H-imidazo[4,5-c]pyridin-2-one

**MF:** C<sub>26</sub>H<sub>25</sub>N<sub>5</sub>O<sub>3</sub>

**FW:** 455.5

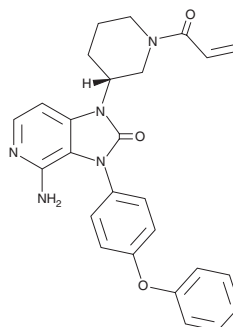
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 222 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

Tolebrutinib is supplied as a solid. A stock solution may be made by dissolving the tolebrutinib in the solvent of choice, which should be purged with an inert gas. Tolebrutinib is soluble in DMSO.

### Description

Tolebrutinib is an irreversible inhibitor of Bruton's tyrosine kinase (BTK; IC<sub>50</sub> = 1 nM).<sup>1</sup> It decreases the number of CD69<sup>+</sup> B cells in isolated mouse whole blood (IC<sub>50</sub> = 200 nM).

### Reference

1. Hopkins, B.T., Bame, E., Bajrami, B., *et al.* Discovery and preclinical characterization of BIIB091, a reversible, selective BTK inhibitor for the treatment of multiple sclerosis. *J. Med. Chem.* **65**(2), 1206-1224 (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, 06/12/2023

#### 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