

# PRODUCT INFORMATION

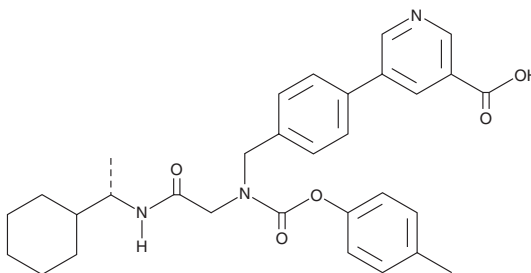


## Tie2 Inhibitor 7

Item No. 19535

**CAS Registry No.:** 1020412-97-8  
**Formal Name:** 5-[4-[[[2-[[[(1S)-1-cyclohexylethyl]amino]-2-oxoethyl] [(4-methylphenoxy)carbonyl]amino]methyl]phenyl]-3-pyridinecarboxylic acid

**MF:** C<sub>31</sub>H<sub>35</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 529.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 257 nm  
**Supplied as:** A crystalline 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

Tie2 Inhibitor 7 is supplied as a crystalline solid. A stock solution may be made by dissolving the Tie2 inhibitor 7 in the solvent of choice, which should be purged with an inert gas. Tie2 Inhibitor 7 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Tie2 inhibitor 7 in these solvents is approximately 1, 20, and 14 mg/ml, respectively.

Tie2 Inhibitor 7 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Tie2 inhibitor 7 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Tie2 Inhibitor 7 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Tie2 Inhibitor 7 blocks Tie2 kinase activity with a K<sub>i</sub> value of 1.3 μM.<sup>1</sup> It has been shown to inhibit angiopoietin 1-induced Tie2 autophosphorylation and downstream signaling with an IC<sub>50</sub> value of 0.3 μM.<sup>1</sup> This compound can prevent endothelial cell tube formation and aberrant vessel growth in a rat model of Matrigel-induced choroidal neovascularization.<sup>1</sup>

### Reference

1. Liu, J., Lin, T.H., Cole, A.G., *et al.* Identification and characterization of small-molecule inhibitors of Tie2 kinase. *FEBS Lett.* **582**(5), 785-791 (2008).

#### 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](http://WWW.CAYMANCHEM.COM)