

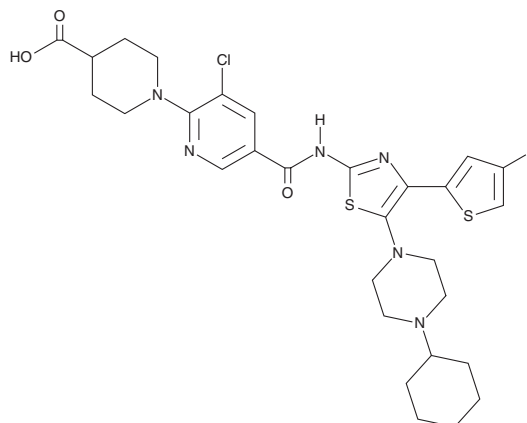
# PRODUCT INFORMATION



## Avatrombopag

Item No. 26535

**CAS Registry No.:** 570406-98-3  
**Formal Name:** 1-[3-chloro-5-[[[4-(4-chloro-2-thienyl)-5-(4-cyclohexyl-1-piperazinyl)-2-thiazolyl]amino]carbonyl]-2-pyridinyl]-4-piperidinecarboxylic acid  
**Synonyms:** AKR 501, AS 1670542, YM477  
**MF:** C<sub>29</sub>H<sub>34</sub>Cl<sub>2</sub>N<sub>6</sub>O<sub>3</sub>S<sub>2</sub>  
**FW:** 649.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 325 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

Avatrombopag is supplied as a crystalline solid. A stock solution may be made by dissolving the avatrombopag in the solvent of choice, which should be purged with an inert gas. Avatrombopag is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

### Description

Avatrombopag is an agonist of the thrombopoietin receptor.<sup>1</sup> It induces proliferation of Ba/F3 cells expressing the thrombopoietin receptor (EC<sub>50</sub> = 3.3 nM) and induces differentiation of hematopoietic progenitor cells into megakaryotes. Avatrombopag is species-specific, inducing STAT5 phosphorylation in human and chimpanzee blood platelets, but not in olive baboon, rhesus monkey, rat, hamster, or several other species blood platelets. It increases the number of human platelets in NOD/SCID mice transplanted with CD34<sup>+</sup> human fetal liver cells when administered at doses of 1 and 3 mg/kg. Formulations containing avatrombopag have been used in the treatment of thrombocytopenia in adults.

### Reference

1. Fukushima-Shintani, M., Suzuki, K., Iwatsuki, Y., *et al.* AKR-501 (YM477) a novel orally-active thrombopoietin receptor agonist. *Eur. J. Haematol.* **82**(4), 247-254 (2009).

#### 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, 12/07/2022

#### 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