

PRODUCT INFORMATION



Lusutrombopag

Item No. 26533

CAS Registry No.: 1110766-97-6
Formal Name: (2E)-3-[2,6-dichloro-4-[[[4-[3-[(1S)-1-(hexyloxy)ethyl]-2-methoxyphenyl]-2-thiazolyl]amino]carbonyl]phenyl]-2-methyl-2-propenoic acid

MF: C₂₉H₃₂Cl₂N₂O₅S
FW: 591.5

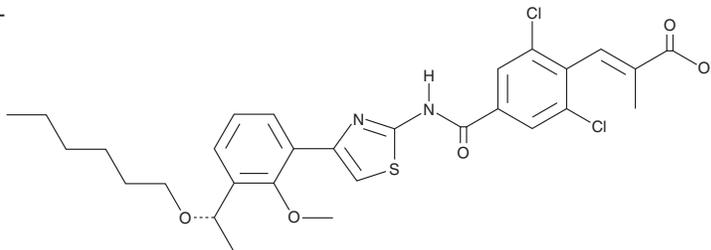
Purity: ≥98%

UV/Vis.: λ_{max}: 218, 260, 322 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

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

Description

Lusutrombopag is an agonist of the thrombopoietin receptor.¹ It induces proliferation of Ba/F3 cells expressing human *MPL*, the gene for the thrombopoietin receptor, with an EC₅₀ value of 84 nM but not of those expressing mouse *Mpl*.² Lusutrombopag (0.3-10 mg/kg per day) increases the number of circulating platelets in a transgenic mouse model expressing human *MPL* and increases the number of megakaryocytes in the bone marrow when administered at a dose of 10 mg/kg per day. Formulations containing lusutrombopag have been used in the treatment of thrombocytopenia in adult patients with chronic liver disease scheduled to undergo invasive procedures.

References

1. Takayama, M. and Kurose, N. Pharmaceutical composition containing optically active compound having thrombopoietin receptor agonist activity, and intermediate therefor. *Shionogi & Co., Ltd.*, **US00888.9722B2** (2014).
2. Yoshida, H., Yamada, H., Nogami, W., *et al.* Development of a new knock-in mouse model and evaluation of pharmacological activities of lusutrombopag, a novel, nonpeptidyl small-molecule agonist of the human thrombopoietin receptor c-Mpl. *Exp. Hematol.* **59**, 30-39 (2018).

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