

PRODUCT INFORMATION



Elinogrel

Item No. 25976

CAS Registry No.: 936500-94-6
Formal Name: 5-chloro-N-[[[4-[6-fluoro-1,4-dihydro-7-(methylamino)-2,4-dioxo-3(2H)-quinazoliny]phenyl]amino]carbonyl]-2-thiophenesulfonamide

Synonyms: PRT060128, PRT128

MF: C₂₀H₁₅ClFN₅O₅S₂

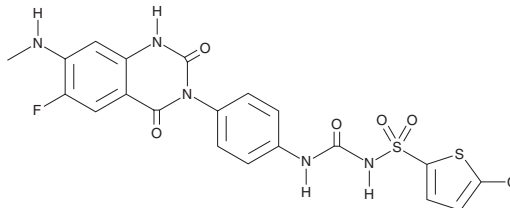
FW: 523.9

Purity: ≥98%

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

Elinogrel is supplied as a solid. A stock solution may be made by dissolving the elinogrel in the solvent of choice, which should be purged with an inert gas. Elinogrel is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

Elinogrel is a reversible antagonist of the platelet purinergic P2Y₁₂ receptor (K_i = 23 nM) that inhibits platelet aggregation in a human platelet-rich plasma (hPRP) assay (IC₅₀ = 2.81 nM).¹ It inhibits thrombosis in a mouse model of vascular injury induced by iron (III) chloride (FeCl₃) when administered at a dose of 60 mg/kg.²

References

1. Zech, G., Hessler, G., Evers, A., *et al.* Identification of high-affinity P2Y₁₂ antagonists based on a phenylpyrazole glutamic acid piperazine backbone. *J. Med. Chem.* **55**(20), 8615-8629 (2012).
2. Haberstoc-Debic, H., Andre, P., Mills, S., *et al.* A clopidogrel-insensitive inducible pool of P2Y(12) receptors contributes to thrombus formation: Inhibition by elinogrel, a direct-acting, reversible P2Y(12) antagonist. *J. Pharmacol. Exp. Ther.* **339**(1), 54-61 (2011).

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, 11/02/2022

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