

PRODUCT INFORMATION

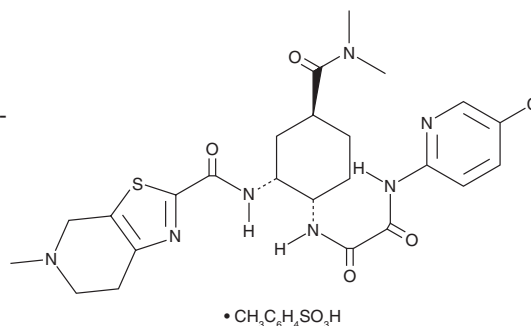


Edoxaban (tosylate)

Item No. 24030

CAS Registry No.: 480449-71-6
Formal Name: N¹-(5-chloro-2-pyridinyl)-N²-[(1S,2R,4S)-4-[[dimethylamino]carbonyl]-2-[[[(4,5,6,7-tetrahydro-5-methylthiazolo[5,4-c]pyridin-2-yl)carbonyl]amino]cyclohexyl]-4-methylbenzenesulfonate, ethanediamide

Synonym: DU-176b
MF: C₂₄H₃₀ClN₇O₄S • C₇H₈O₃S
FW: 720.3
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

Edoxaban (tosylate) is supplied as a solid. A stock solution may be made by dissolving the edoxaban (tosylate) in the solvent of choice. Edoxaban (tosylate) is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

Edoxaban is an orally bioavailable and potent inhibitor of factor Xa (K_i s = 0.561 and 0.903 nM for free and complexed human factor Xa, respectively).¹ It is >10,000-fold selective for factor Xa over thrombin, FVIIa/sTF, FXIa, tPA, aPC, trypsin, plasmin, and chymotrypsin. Edoxaban prolongs prothrombin, activated partial thromboplastin, and thrombin clotting times in a concentration-dependent manner *ex vivo* in human plasma. Oral administration of edoxaban (0.5-12.5 mg/kg) reduces thrombus formation and prolongs prothrombin time in a dose-dependent manner in rat and rabbit models of venous stasis thrombosis. It also reduces thrombus formation in a rat model of platinum wire-induced venous thrombosis. Formulations containing edoxaban have been used to prevent stroke in patients with atrial fibrillation.

Reference

1. Furugohri, T., Isobe, K., Honda, Y., *et al.* DU-176b, a potent and orally active factor Xa inhibitor: *In vitro* and *in vivo* pharmacological profiles. *J. Thromb. Haemost.* **6(9)**, 1542-1549 (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, 12/22/2022

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