

PRODUCT INFORMATION

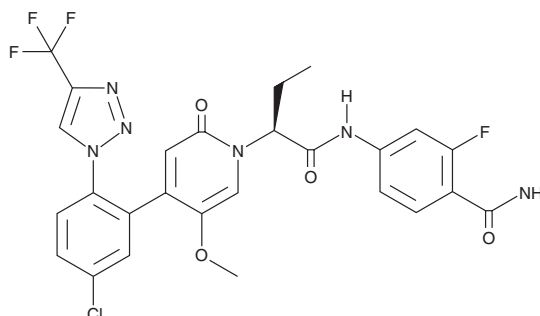


BAY-2433334

Item No. 41101

CAS Registry No.: 2064121-65-7
Formal Name: (αS)-N-[4-(aminocarbonyl)-3-fluorophenyl]-4-[5-chloro-2-[4-(trifluoromethyl)-1H-1,2,3-triazol-1-yl]phenyl]-α-ethyl-5-methoxy-2-oxo-1(2H)-pyridineacetamide

Synonym: Asundexian
MF: C₂₆H₂₁ClF₄N₆O₄
FW: 592.9
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

BAY-2433334 is supplied as a solid. A stock solution may be made by dissolving the BAY-2433334 in the solvent of choice, which should be purged with an inert gas. BAY-2433334 is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

BAY-2433334 is an inhibitor of Factor XIa ($IC_{50} = 0.92$ nM).¹ It is greater than 1,000-fold selective for Factor Xa over FVIIa, FIXa, FXa, thrombin, urokinase, tissue plasminogen activator, and plasmin. BAY-2433334 increases clotting time in isolated human plasma ($EC_{150} = 0.2$ μM). *In vivo*, BAY-2433334 (10 and 30 mg/kg) reduces thrombus weight in a rabbit model of thrombosis induced by ferric chloride injury of the carotid artery.

Reference

1. Roehrig, S., Ackerstaff, J., Jiménez Núñez, E., *et al.* Design and preclinical characterization program toward asundexian (BAY 2433334), an oral factor XIa inhibitor for the prevention and treatment of thromboembolic disorders. *J. Med. Chem.* **66**(17), 12203-12224 (2023).

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, 06/24/2024

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