

PRODUCT INFORMATION



Acebilustat

Item No. 33270

CAS Registry No.: 943764-99-6
Formal Name: 4-[[[(1S,4S)-5-[[4-[4-(2-oxazolyl)phenoxy]phenyl]methyl]-2,5-diazabicyclo[2.2.1]hept-2-yl]methyl]-benzoic acid

Synonym: CTX-4430

MF: C₂₉H₂₇N₃O₄

FW: 481.5

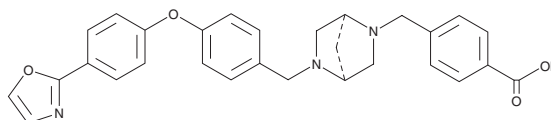
Purity: ≥98%

UV/Vis.: λ_{max}: 274 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

Acebilustat is supplied as a crystalline solid. A stock solution may be made by dissolving the acebilustat in the solvent of choice, which should be purged with an inert gas. Acebilustat is slightly soluble in chloroform, methanol, and DMSO.

Description

Acebilustat is an inhibitor of leukotriene A₄ (LTA₄) hydrolase, the rate-limiting enzyme in LTB₄ synthesis.¹ It reduces the transmigration of polymorphonuclear neutrophils (PMNs) from an epithelial monolayer to airway supernatant derived from sputum isolated from patients with cystic fibrosis when used at a concentration of 1 μM.²

References

1. Elborn, J.S., Horsley, A., MacGregor, G., *et al.* Phase I studies of acebilustat: Biomarker response and safety in patients with cystic fibrosis. *Clin. Transl. Sci.* **10(1)**, 28-34 (2017).
2. Forrest, O.A., Ingersoll, S.A., Preininger, M.K., *et al.* Frontline Science: Pathological conditioning of human neutrophils recruited to the airway milieu in cystic fibrosis. *J. Leukoc. Biol.* **104(4)**, 665-675 (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, 12/08/2022

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