

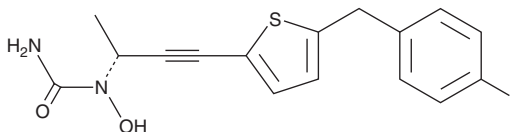
PRODUCT INFORMATION



Atreleuton

Item No. 28736

CAS Registry No.: 154355-76-7
Formal Name: N-[(1R)-3-[5-[(4-fluorophenyl)methyl]-2-thienyl]-1-methyl-2-propyn-1-yl]-N-hydroxy-urea
Synonyms: A-85761.0, ABT-761, VIA-2291
MF: C₁₆H₁₅FN₂O₂S
FW: 318.4
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

Atreleuton is supplied as a solid. A stock solution may be made by dissolving the atreleuton in the solvent of choice, which should be purged with an inert gas. Atreleuton is soluble in the organic solvent methanol.

Description

Atreleuton is an inhibitor of 5-lipoxygenase (5-LO; IC₅₀ = 23 nM in rat basophil leukemia cell lysates).¹ It inhibits leukotriene B₄ (LTB₄; Item No. 20110) formation induced by A23187 (Item Nos. 11016 | 22030) in isolated human whole blood (IC₅₀ = 160 nM). Atreleuton inhibits LTE₄ and LTB₄ biosynthesis in the peritoneal cavity in a rat model of anaphylaxis (ED₅₀s = 1.4 and 0.6 mg/kg, respectively) and bronchoconstriction induced by arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) in anesthetized guinea pigs (ED₅₀ = 3 mg/kg).²

References

1. Brooks, C.D., Stewart, A.O., Basha, A., *et al.* (R)-(+)-N-[3-[5-[(4-fluorophenyl)methyl]-2-thienyl]-1-methyl-2-propynyl]-N-hydroxyurea (ABT-761), a second-generation 5-lipoxygenase inhibitor. *J. Med. Chem.* **38**(24), 4768-4778 (1995).
2. Bell, R.L., Harris, R.R., Malo, P.E., *et al.* ABT-761 attenuates bronchoconstriction and pulmonary inflammation in rodents. *J. Pharmacol. Exp. Ther.* **280**(3), 1366-1373 (1997).

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, 10/13/2022

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