

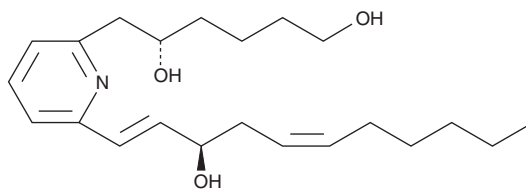
PRODUCT INFORMATION



U-75302

Item No. 70705

CAS Registry No.: 119477-85-9
Formal Name: 6-(6-(3R-hydroxy-1E,5Z-undecadien-1-yl)-2-pyridinyl)-1,5S-hexanediol
MF: C₂₂H₃₅NO₃
FW: 361.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

U-75302 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of U-75302 in these solvents is approximately 50, 10, and 7.1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of U-75302 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of U-75302 in PBS (pH 7.2) is approximately 80 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

U-75302 is an LTB₄ receptor antagonist with a K_i of 159 nM on guinea pig lung membranes.^{1,2} This activity is specific for the BLT₁ receptor; U-75302 does not antagonize the binding of [³H]-LTB₄ to the human BLT₂ receptor.³

References

1. Tebo, J., Der, S., Frevel, M., *et al.* Heterogeneity in control of mRNA stability by AU-rich elements. *J. Biol. Chem.* **278**(14), 12085-12093 (2003).
2. Lin, A.H., Morris, J., Wishka, D.G., *et al.* Novel molecules that antagonize leukotriene B₄ binding to neutrophils. *Ann. N.Y. Acad. Sci.* **524**, 196-200 (1988).
3. Wang, S., Gustafson, E., Pang, L., *et al.* A novel hepatointestinal leukotriene B₄ receptor. *J. Biol. Chem.* **275**(52), 40686-40694 (2000).

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, 02/01/2024

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