

PRODUCT INFORMATION



MN-001

Item No. 11948

CAS Registry No.: 125961-82-2
Formal Name: 4-[6-acetyl-3-[3-[(4-acetyl-3-hydroxy-2-propylphenyl)thio]propoxy]-2-propylphenoxy]-butanoic acid

Synonyms: KCA 757, Tipelukast

MF: C₂₉H₃₈O₇S

FW: 530.7

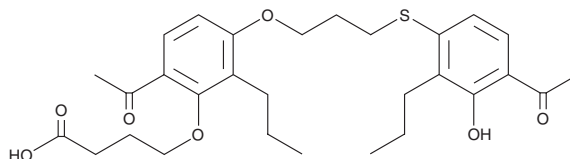
Purity: ≥98%

UV/Vis.: λ_{max}: 216, 252, 311 nm

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

MN-001 is supplied as a solid. A stock solution may be made by dissolving the MN-001 in the solvent of choice, which should be purged with an inert gas. MN-001 is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml. MN-001 is also slightly soluble in chloroform and ethyl acetate.

Description

MN-001 is an anti-inflammatory agent.¹ *In vivo*, MN-001 (10, 30, and 50 mg/kg) reduces the frequency of bladder voids, increases the intercontractile interval, and decreases the number of non-voiding contractions in a rat model of bladder hyperactivity induced by ovalbumin-induced mast cell stimulation.

Reference

1. Rajasekaran, M., Locke, K.W., and Parsons, C.L. MN-001, a novel oral anti-inflammatory agent, suppresses bladder hyperactivity in a rat model. *BJU Int.* **98(2)**, 430-434 (2006).

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, 11/11/2022

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