

PRODUCT INFORMATION

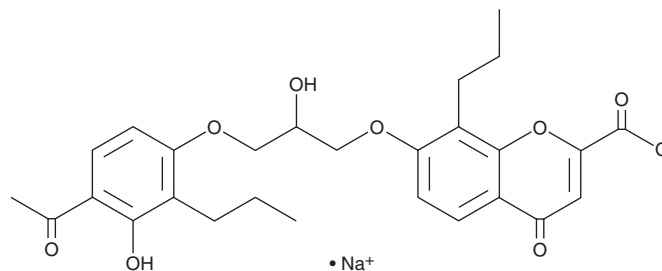


FPL 55712 (sodium salt)

Item No. 35433

CAS Registry No.: 40786-08-1
Formal Name: 7-[3-(4-acetyl-3-hydroxy-2-propylphenoxy)-2-hydroxypropoxy]-4-oxo-8-propyl-4H-1-benzopyran-2-carboxylic acid, monosodium salt

MF: $C_{27}H_{29}O_9 \cdot Na$
FW: 520.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 214, 286 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

FPL 55712 (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the FPL 55712 (sodium salt) in the solvent of choice, which should be purged with an inert gas. FPL 55712 (sodium salt) is soluble in DMSO.

Description

FPL 55712 (sodium salt) is a leukotriene (LT) receptor antagonist.¹ It inhibits the binding of LTD₄ (Item No. 20310) to guinea pig lung homogenates ($IC_{50} = 4 \mu M$) and reduces LTD₄-induced contraction of isolated guinea pig tracheal smooth muscle ($pA_2 = 1 \mu M$). FPL 55712 (sodium salt) inhibits bronchoconstriction induced by LTD₄, LTC₄ (Item No. 20210), or LTE₄ (Item No. 20410) in guinea pigs ($ID_{50}S = 2, 6, \text{ and } 1 \text{ mg/kg}$, respectively).

Reference

1. Cohen, N., Weber, G., Banner, B.L., *et al.* 3,4-Dihydro-2H-1-benzopyran-2-carboxylic acids and related compounds as leukotriene antagonists. *J. Med. Chem.* **32(8)**, 1842-1860 (1989).

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, 12/12/2022

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