

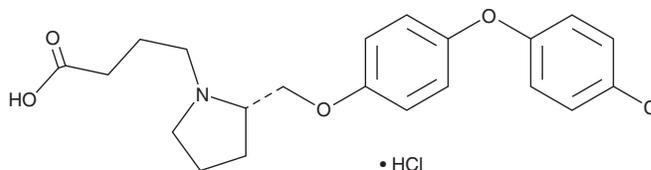
PRODUCT INFORMATION



DG-051

Item No. 21688

CAS Registry No.: 929915-58-2
Formal Name: (2S)-2-[[4-(4-chlorophenoxy)phenoxy]methyl]-1-pyrrolidinebutanoic acid, monohydrochloride
MF: C₂₁H₂₄ClNO₄ • HCl
FW: 426.3
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 279 nm
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

DG-051 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of DG-051 in these solvents is approximately 12.5 and 20 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 DG-051 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of DG-051 in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

DG-051 is a potent, orally bioavailable leukotriene A₄ hydrolase (LTA₄H) inhibitor (K_d = 25 nM; IC₅₀ = 47 nM), the enzyme that catalyzes the rate-determining step in the synthesis of LTB₄ (Item No. 20110).¹ However, it also inhibits the aminopeptidase activity of LTA₄H with IC₅₀ values of 72 and 150 nM when tested using the substrates L-alanine *p*-nitroanaline and prolyl-glycyl-proline peptide (PGP; Item No. 11188), respectively, potentially leading to PGP accumulation *in vivo*.^{1,2} DG-051 does not show any inhibitory activity against a panel of more than 50 additional aminopeptidases, ion channels, or HERG.¹

References

1. Sandanayaka, V., Mamat, B., Mishra, R.K., *et al.* Discovery of 4-[(2S)-2-[[4-(4-chlorophenoxy)phenoxy]methyl]-1-pyrrolidinyl]butanoic acid (DG-051) as a novel leukotriene A₄ hydrolase inhibitor of leukotriene B₄ biosynthesis. *J. Med. Chem.* **53**(2), 573-585 (2010).
2. Low, C.M., Akthar, S., Patel, D.F., *et al.* The development of novel LTA₄H modulators to selectively target LTB₄ generation. *Sci. Rep.* **7**, 44449 (2017).

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, 10/05/2022

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