

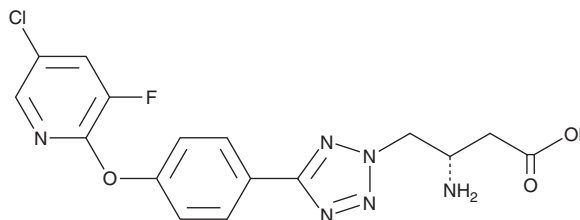
PRODUCT INFORMATION



LYS-006

Item No. 33706

CAS Registry No.: 1799681-85-8
Formal Name: βS-amino-5-[4-[(5-chloro-3-fluoro-2-pyridinyl)oxy]phenyl]-2H-tetrazole-2-butanoic acid
MF: C₁₆H₁₄ClFN₆O₃
FW: 392.8
Purity: ≥98%
UV/Vis.: λ_{max}: 251 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

LYS-006 is supplied as a solid. A stock solution may be made by dissolving the LYS-006 in the solvent of choice, which should be purged with an inert gas. LYS-006 is soluble in the organic solvent DMSO.

Description

LYS-006 is an inhibitor of leukotriene A₄ (LTA₄) hydrolase.^{1,2}

References

1. Loesche, C., Penno, C., Roehn, T., *et al.* LTA₄H inhibitors for the treatment of hidradenitis suppurativa. *Novartis AG. WO2020144604A1* (2020).
2. Folkes, A.S., Hawatmeh, F.Z., Wong, A., *et al.* Emerging drugs for the treatment of hidradenitis suppurativa. *Expert Opin. Emerg. Drugs* **25(3)**, 201-211 (2020).

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, 12/07/2022

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