

PRODUCT INFORMATION

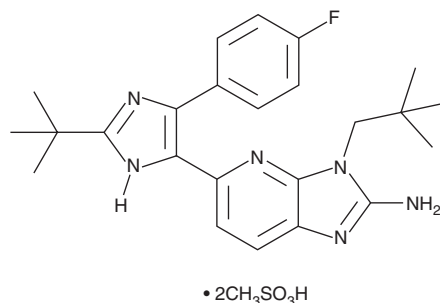


LY2228820 (mesylate)

Item No. 23259

CAS Registry No.: 862507-23-1
Formal Name: 5-[2-(1,1-dimethylethyl)-4-(4-fluorophenyl)-1H-imidazol-5-yl]-3-(2,2-dimethylpropyl)-3H-imidazo[4,5-b]pyridin-2-amine, dimethanesulfonate

Synonym: LSN2322600
MF: C₂₄H₂₉FN₆ • 2CH₃SO₃H
FW: 612.7
Purity: ≥98%
UV/Vis.: λ_{max}: 321 nm
Supplied as: A crystalline 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

LY2228820 (mesylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2228820 (mesylate) in the solvent of choice, which should be purged with an inert gas. LY2228820 (mesylate) is sparingly soluble in ethanol, methanol, and DMSO.

Description

LY2228820 is a potent inhibitor of p38α MAP kinase (IC₅₀ = 7 nM for human recombinant p38α).¹ It inhibits LPS-induced TNF-α production in murine peritoneal macrophages and reduces phosphorylation of MAPKAPK-2 (MK2) in RAW 264.7 cells (IC₅₀s = 5.2 and 34.3 nM, respectively). LY2228820 inhibits TNF-α production in LPS-treated mice and reduces paw swelling and cartilage destruction in a rat model of chronic inflammation with threshold minimum 50% doses (TMED₅₀) values of <1, 1.5, and 1.5 mg/kg, respectively.

Reference

1. Mader, M., de Dios, A., Shih, C., *et al.* Imidazolyl benzimidazoles and imidazo[4,5-b]pyridines as potent p38α MAP kinase inhibitors with excellent in vivo antiinflammatory properties. *Bioorg. Med. Chem. Lett.* **18**(1), 179-183 (2008).

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, 11/22/2022

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