

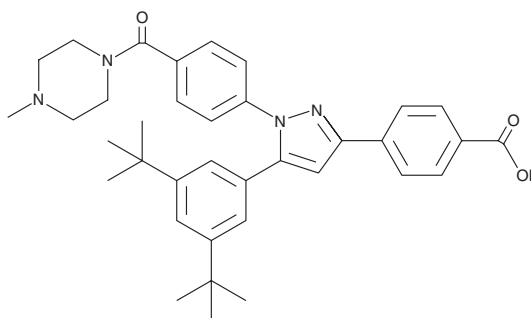
PRODUCT INFORMATION



LY2955303

Item No. 25833

CAS Registry No.: 1433497-19-8
Formal Name: 4-[5-[3,5-bis(1,1-dimethylethyl)phenyl]-1-[4-[(4-methyl-1-piperazinyl)carbonyl]phenyl]-1H-pyrazol-3-yl]-benzoic acid
MF: C₃₆H₄₂N₄O₃
FW: 578.7
Purity: ≥95%
UV/Vis.: λ_{max}: 285 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

LY2955303 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2955303 in the solvent of choice, which should be purged with an inert gas. LY2955303 is soluble in the organic solvent chloroform at a concentration of approximately 25 mg/ml.

Description

LY2955303 is a potent antagonist of retinoic acid receptor γ (RAR γ ; K_i = 1.1 nM).¹ It is selective for RAR γ over RAR α and RAR β (K_is = >1,700 and >2,980 nM, respectively). LY2955303 reduces differential weight bearing between the contralateral and ipsilateral limbs in a rat model of monoiodoacetate-induced osteoarthritis (ED₅₀ = 0.72 mg/kg).

Reference

1. Hughes, N.E., Bleisch, T.J., Jones, S.A., *et al.* Identification of potent and selective retinoic acid receptor gamma (RAR γ) antagonists for the treatment of osteoarthritis pain using structure based drug design. *Bioorg. Med. Chem.* **26(14)**, 3274-3277 (2016).

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, 12/22/2022

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