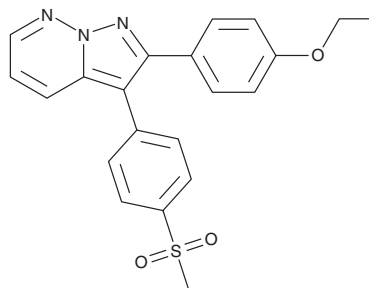


PRODUCT INFORMATION



GW 406381
Item No. 33895

CAS Registry No.: 221148-46-5
Formal Name: 2-(4-ethoxyphenyl)-3-[4-(methylsulfonyl)phenyl]-pyrazolo[1,5-b]pyridazine
Synonym: GW406381X
MF: C₂₁H₁₉N₃O₃S
FW: 393.5
Purity: ≥98%
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

GW 406381 is supplied as a solid. A stock solution may be made by dissolving the GW 406381 in the solvent of choice, which should be purged with an inert gas. GW 406381 is soluble in the organic solvent DMSO (sonicated) at a concentration of approximately 25 mg/ml.

Description

GW 406381 is a non-steroidal anti-inflammatory drug (NSAID) and COX-2 inhibitor (IC₅₀ = 3 nM).¹ It is selective for COX-2 over COX-1 (IC₅₀ = >84.2 μM). GW 406381 (5 mg/kg) reverses chronic constriction injury-induced decreases in paw withdrawal thresholds in a rat model of neuropathic pain.²

References

1. Beswick, P., Bingham, S., Bountra, C., *et al.* Identification of 2,3-diaryl-pyrazolo[1,5-b]pyridazines as potent and selective cyclooxygenase-2 inhibitors. *Bioorg. Med. Chem. Lett.* **14(21)**, 5445-5448 (2004).
2. Zhao, F.-Y., Spanswick, D., Martindale, J.C., *et al.* GW406381, a novel COX-2 inhibitor, attenuates spontaneous ectopic discharge in sural nerves of rats following chronic constriction injury. *Pain* **128(1-2)**, 78-87 (2007).

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, 12/05/2022

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