

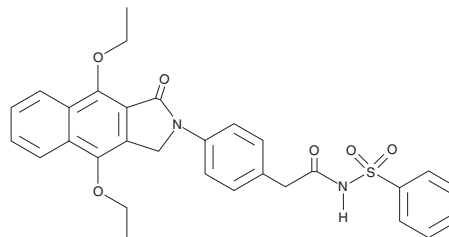
PRODUCT INFORMATION



GW 627368X

Item No. 10009162

CAS Registry No.: 439288-66-1
Formal Name: 4-(4,9-diethoxy-1,3-dihydro-1-oxo-2H-benz[f]isoindol-2-yl)-N-(phenylsulfonyl)benzeneacetamide
MF: C₃₀H₂₈N₂O₆S
FW: 544.6
Purity: ≥96%
UV/Vis.: λ_{max}: 217, 229, 248, 295, 353 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

GW 627368X is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 627368X in the solvent of choice, which should be purged with an inert gas. GW 627368X is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of GW 627368X in these solvents is approximately 10 and 30 mg/ml, respectively.

GW 627368X is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GW 627368X should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GW 627368X has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The effects of prostaglandin E₂ (PGE₂) are transduced by at least four distinct receptors designated EP₁, EP₂, EP₃, and EP₄.¹ GW 627368X is a potent and selective competitive antagonist of the EP₄ receptor with additional human TP receptor affinity. In competition radioligand bioassays, GW 627368X had affinity for human EP₄ and TP receptors with K_i values of 100 nM and 158 nM, respectively.² Affinity for all other human prostanoid receptors is >5.0 μM. In human washed platelets, GW 627368X produced 100% inhibition of U-46619 (EC₁₀₀)-induced aggregation at a concentration of 10 μM.

References

1. Coleman, R.A., Smith, W.L., and Narumiya, S. Classification of prostanoid receptors: Properties, distribution, and structure of the receptors and their subtypes. *Pharmacol. Rev.* **46**, 205-229 (1994).
2. Wilson, R.J., Giblin, G.M.P., Roomans, S., et al. GW627368X ((N-{2-[4-(4,9-diethoxy-1-oxo-1,3-dihydro-2H-benzo[f]isoindol-2-yl)phenyl]acetyl} benzene sulphonamide): A novel, potent and selective prostanoid EP₄ receptor antagonist. *Br. J. Pharmacol.* **148**, 326-339 (2006).

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, 05/02/2024

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