

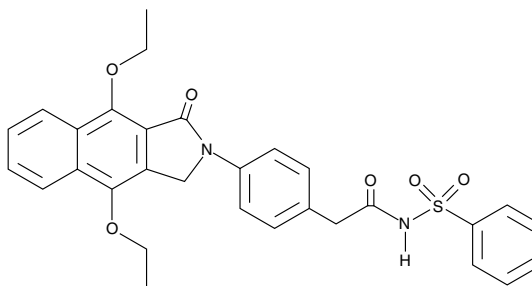
Product Information



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%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 217, 229, 248, 295, 353 nm



Laboratory Procedures

For long term storage, we suggest that GW 627368X be stored as supplied at -20°C. It will be stable for at least two years.

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

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.

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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10009162

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 04/04/2012