

PRODUCT INFORMATION



U-46619 Glycine methyl ester

Item No. 10010522

Formal Name: 9,11-dideoxy-9 α ,11 α -methanoepoxy-prosta-5Z,13E-dien-1-oic acid, methyl ester

MF: C₂₄H₃₉NO₅

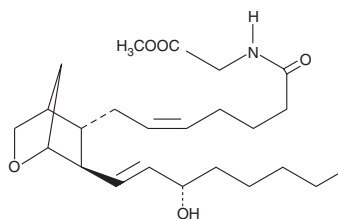
FW: 421.6

Purity: \geq 98%

Supplied as: A solution in methyl acetate

Storage: -20°C

Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

U-46619 Glycine methyl ester is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of U-46619 glycine methyl ester in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of U-46619 glycine methyl ester is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of U-46619 glycine methyl ester in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

U-46619 is a stable analog of the endoperoxide PGH₂, and a TP receptor agonist.¹ It exhibits properties similar to TXA₂, causing platelet shape change, aggregation, and contraction of vascular smooth muscle.^{2,3} Mean EC₅₀ values for shape change in human, rat, and rabbit platelets are 4.8, 6.0, and 7.3 nM respectively, and for aggregation, are 82, 145, and 65 nM, respectively.⁴ U-46619 glycine methyl ester contains a modification at the C-1 position of U-46619 that may uniquely alter its binding properties to the TP receptor or any of the PGH₂-metabolizing enzymes. As a stable PGH₂ analog, it could therefore be a useful tool to explore the inhibition of various enzymes in the arachidonic acid metabolic pathway. U-46619 glycine methyl ester may also serve as a lipophilic prodrug form of U-46619 that will alter its distribution and pharmacokinetic properties. There are no published reports on the biological activity of U-46619 glycine methyl ester.

References

1. Abramovitz, M., Adam, M., Boie, Y., *et al.* The utilization of recombinant prostanoid receptors to determine the affinities and selectivities of prostaglandins and related analogs. *Biochim. Biophys. Acta* **1483**(2), 285-293 (2000).
2. Coleman, R.A., Humphrey, P.P.A., Kennedy, I., *et al.* Comparison of the actions of U-46619, a prostaglandin H₂-analogue, with those of prostaglandin H₂ and thromboxane A₂ on some isolated smooth muscle preparations. *Br. J. Pharmacol.* **73**(3), 773-778 (1981).
3. Liel, N., Mais, D.E., and Halushka, P.V. Binding of a thromboxane A₂/prostaglandin H₂ agonist [³H]U46619 to washed human platelets. *Prostaglandins* **33**(6), 789-797 (1987).
4. Tymkewycz, P.M., Jones, R.L., Wilson, N.H., *et al.* Heterogeneity of thromboxane A₂ (TP-) receptors: Evidence from antagonist but not agonist potency measurements. *Br. J. Pharmacol.* **102**(3), 607-614 (1991).

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, 03/07/2024

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