PRODUCT INFORMATION



U-51605

Item No. 16465

CAS Registry No.:	64192-56-9	
Formal Name:	9α,11α-azoprosta-	
	5Z,13E-dien-1-oic acid	
MF:	C ₂₀ H ₃₂ N ₂ O ₂	
FW:	332.5	N COOH
Purity:	≥98%	
Supplied as:	A solution in methyl acetate	
Storage:	-80°C	
Stability:	≥2 years	

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

Laboratory Procedures

U-51605 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 purged with an inert gas can be used. The solubility of U-51605 in these solvents is approximately 100 mg/ml.

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-51605 is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of U-51605 in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

U-51605 is a stable analog of the endoperoxide prostaglandin H₂ (PGH₂). It is an inhibitor of both prostacyclin (PGI) and thromboxane (TX) synthases with more selectivity towards PGI synthase. U-51605 is also a partial agonist at TP receptors.¹ In human foreskin fibroblasts, U-51605 inhibits PGI synthase at a concentration of 2.8 μ M, whereas, human platelet TX synthase is inhibited at a concentration of 5.6 μ M.^{2,3} At concentrations up to 1 μ M, U-51605 reduced the release of prostacyclin in SHR aorta elicited by the calcium ionophore A-23187 with no effect on TXA₂ production and yet significantly increased PGE₂ and PGF_{2a} release.⁴

References

- 1. Needleman, P., Bryan, B., Wyche, A., et al. Thromboxane synthetase inhibitors as pharmacological tools: Differential biochemical and biological effects on platelet suspensions. Prostaglandins 14, 897-907 (1977).
- 2. Gorman, R.R., Hamilton, R.D., and Hopkins, N.K. Stimulation of human foreskin fibroblast adenosine 3':5'-cyclic monophosphate levels by prostacyclin (prostaglandin I₂). J. Biol. Chem. 254, 1671-1676 (1979).
- Gorman, R.R., Bundy, G.L., Peterson, D.C., et al. Inhibition of human platelet thromboxane synthetase by 9,11-azoprosta-5,13-dienoic acid. Proc. Natl. Acad. Sci. USA 74, 4007-4011 (1977).
- 4. Gluais, P., Paysant, J., Badier-Commander, C., et al. In SHR aorta, calcium ionophore A-23187 releases prostacyclin and thromboxane A₂ as endothelium-derived contracting factors. Am. J. Physiol. Heart Circ. Physiol. 291, H2255-H2264 (2006).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 10/10/2023

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