

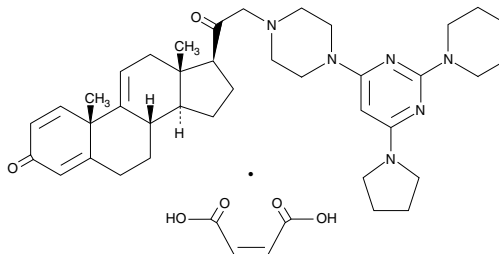
Product Information



U-74389G

Item No. 75860

CAS Registry No.: 153190-29-5
Formal Name: 21-[4-(2,6-di-1-pyrrolidinyl-4-pyrimidinyl)-1-piperazinyl]-pregna-1,4,9(11)-triene-3,20-dione maleate salt
Synonym: Methylated Tirilazad
MF: C₃₈H₅₂N₆O₂ • C₄H₄O₄
FW: 740.9
Purity: ≥99%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 237, 287 nm



Laboratory Procedures

For long term storage, we suggest that U-74389G be stored as supplied at -20°C. It should be stable for at least one year.

U-74389G is supplied as a crystalline solid. A stock solution may be made by dissolving the U-74389G in an organic solvent purged with an inert gas. U-74389G is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of U-74389G in these solvents is approximately 1, 35, and 36 mg/ml, respectively.

U-74389G is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, U-74389G should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. U-74389G has a solubility of approximately 0.05 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

U-74389G, is an antioxidant which prevents iron-dependent lipid peroxidation. It protects against ischemia-reperfusion injury in animal heart, liver, and kidney models.¹⁻⁴

References

1. Ishizaki, N., Zhu, Y., Zhang, S., *et al.* Comparison of various lazaroid compounds for protection against ischemic liver injury. *Transplantation* **63**, 202-208 (1997).
2. Palma-Vargas, J.M., Toledo, A.H., García-Criado, F.J., *et al.* 21-Aminosteroids block neutrophil infiltration and provide liver protection independent of nitrite/nitrate levels. *J. Surg. Res.* **66**, 131-137 (1996).
3. Shoskes, D.A., Xie, Y., Gonzalez-Cadavid, N.F. Nitric oxide synthase activity in renal ischemia-reperfusion injury in the rat. Implications for renal transplantation. *Transplantation* **63**, 495-500 (1997).
4. Perna, A.M., Liguori, P., Bonacchi, M., *et al.* Protection of rat heart from ischaemia-reperfusion injury by the 21-aminosteroid U-74389G. *Pharmacol. Res.* **34**, 25-31 (1996).

Related Products

Celastrol - Item No. 70950 • BHT - Item No. 89910

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 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, 05/02/2011

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