

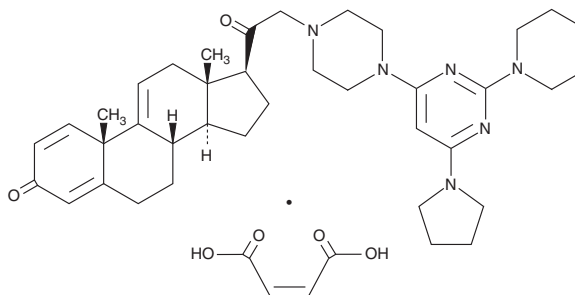
PRODUCT INFORMATION



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: C₃₇H₅₀N₆O₂ • C₄H₄O₄
FW: 726.9
Purity: ≥99%
UV/Vis.: λ_{max}: 237, 287 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

U-74389G is supplied as a crystalline solid. A stock solution may be made by dissolving the U-74389G in the solvent of choice, which should be 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 1.36, 35.3, and 36.3 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 0.5 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.

Description

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).

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, 12/13/2022

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