

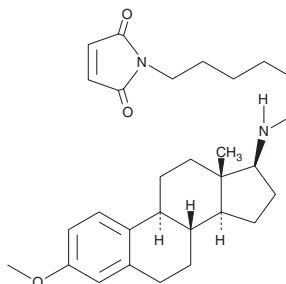
Product Information



U-73122

Catalog No. 70740

CAS Registry No.: 112648-68-7
Formal Name: 1-[6-[[[(17β)-3-methoxyestra-1,3,5(10)-trien-17-yl]amino]hexyl]-1H-pyrrole-2,5-dione
MF: C₂₉H₄₀N₂O₃
FW: 464.6
Purity: ≥99%
Stability: ≥1 year at room temperature
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that U-73122 be stored as supplied at room temperature. It should be stable for at least one year.

U-73122 is supplied as a crystalline solid. A stock solution may be made by dissolving the U-73122 in an organic solvent. U-73122 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of U-73122 in these solvents is approximately 0.5 mg/ml. U-73122 can also be dissolved in ethanol, but the solubility is less than 0.5 mg/ml and it requires 24 hours to go completely into solution.

U-73122 is sparingly soluble in aqueous buffers. Therefore, further dilutions of the organic solvent 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. Store aqueous solutions of U-73122 on ice and use within 12 hours of preparation. Although the aqueous solutions of U-73122 may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

U-73122 is an inhibitor of PLC-dependent processes, however, the mechanism of action remains unclear.¹⁻³ The IC₅₀ values for inhibition of platelet aggregation induced by collagen or thrombin is 0.6 and 5 μM, respectively.² It also exhibits inhibitory activity against HIV-1 integrase with an IC₅₀ value of 7 μM.⁴

References

1. Smith, R.J., Sam, L.M., Justen, J.M., *et al.* Receptor-coupled signal transduction in human polymorphonuclear neutrophils: Effects of a novel inhibitor of phospholipase C-dependent processes on cell responsiveness. *J. Pharmacol. Exp. Ther.* **253**, 688-697 (1990).
2. Bleasdale, J.E., Thakur, N.R., Gremban, R.S., *et al.* Selective inhibition of receptor-coupled phospholipase C-dependent processes in human platelets and polymorphonuclear neutrophils. *J. Pharmacol. Exp. Ther.* **255**, 756-768 (1990).
3. Hildebrandt, J.-P., Plant, T.D., and Meves, H. The effects of bradykinin on K⁺ currents in NG108-15 cells treated with U73122, a phospholipase C inhibitor, or neomycin. *Br. J. Pharmacol.* **120**, 841-850 (1997).
4. Burke, T.R., Jr., Fesen, M.R., Mazumder, A., *et al.* Hydroxylated aromatic inhibitors of HIV-1 integrase. *J. Med. Chem.* **38**, 4171-4178 (1995).

Related Products

D-*myo*-Inositol-1,4,5-triphosphate (potassium salt) - Cat. No. 60960 • D-*myo*-Inositol-1,3,4-triphosphate (potassium salt) - Cat. No. 60970 • D-*myo*-Inositol-1,3,4,5-tetraphosphate (potassium salt) - Cat. No. 60980

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, 01/20/2010

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