

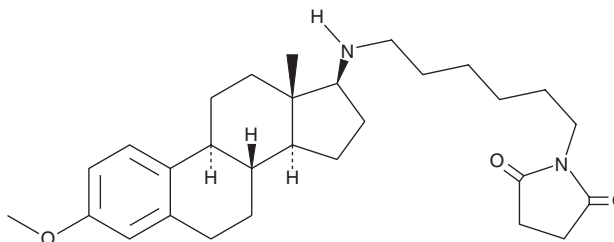
PRODUCT INFORMATION



U-73343

Item No. 17339

CAS Registry No.: 142878-12-4
Formal Name: 1-[6-[[[(17b)-3-methoxyestra-1,3,5(10)-trien-17-yl]amino]hexyl]-2,5-pyrrolidinedione
MF: C₂₉H₄₂N₂O₃
FW: 466.7
Purity: ≥95%
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-73343 is supplied as a crystalline solid. A stock solution may be made by dissolving the U-73343 in the solvent of choice, which should be purged with an inert gas. U-73343 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of U-73343 in these solvents is approximately 0.25, 0.5, and 2 mg/ml, respectively.

Description

U-73122 (Item No. 70740) 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 are 0.6 and 5 μM, respectively.² It also exhibits inhibitory activity against HIV-1 integrase with an IC₅₀ value of 7 μM.⁴ U-73343 is a close structural analog of U-73122 that does not inhibit PLC, making it useful as a negative control for PLC inhibition.¹

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**(2), 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**(2), 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**(5), 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**(21), 4171-4178 (1995).

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, 11/10/2022

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