

U-73343

Phospholipase inhibitor

Inactive analog of U-73122 (Prod. No. BML-ST391) which only slightly inhibits phospholipase C. May be used as a negative control.

Citations: 9

[View Online »](#)

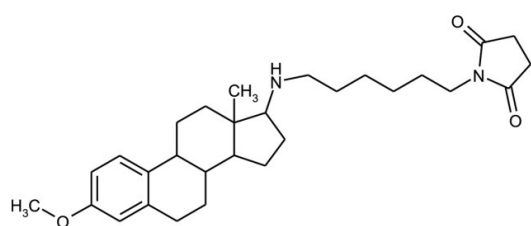
Ordering Information

[Order Online »](#)

BML-ST392-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-[6-[[[(17β)-3-Methoxyestra-1,3,5(10)-trien-17-yl]amino]hexyl]-pyrrolidine-2,5-dione
Appearance	White to off-white solid.
CAS	142878-12-4
Couple Target	Lipoxygenase, Phospholipase, PLC
Couple Type	Inhibitor
Formula	C ₂₉ H ₄₂ N ₂ O ₃
MW	466.7
Purity	≥98% (HPLC)
Solubility	Soluble in methylene chloride or chloroform; slightly soluble in DMSO (2mg/ml) or 100% ethanol (1mg/ml); sparingly soluble in aqueous solutions.

Technical Info / Product Notes Note: Product is not sterile.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com