

# 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 »](#)

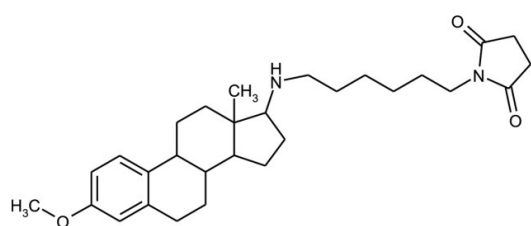
## Ordering Information

[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          | $C_{29}H_{42}N_2O_3$                                                                                                                             |
| 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](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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