

U-73122

Inhibitor of PLC activation

Inhibits agonist induced phospholipase C (PLC) activation ($IC_{50}=1-2.1 \mu M$) in human platelets. It also inhibits agonist induced down-regulation of muscarinic receptors in SK-N-SH neuroblastoma cells. It is a useful tool to investigate receptor-mediated PI turnover in signal transduction. It is a potent inhibitor of human neutrophil adhesion to biological surfaces as well as adhesion-dependent granule exocytosis and oxidative burst. Inhibits 5-lipoxygenase *in vivo* and *in vitro* ($IC_{50} = 30nM$ for recombinant enzyme and $2.4\mu M$ for PMNL homogenates).

Citations: 28

[View Online »](#)

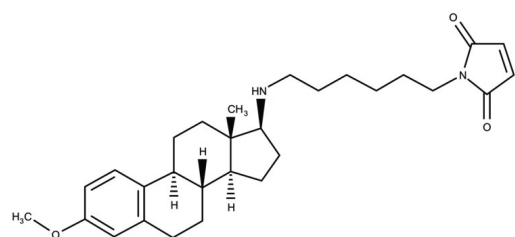
Ordering Information

[Order Online »](#)

BML-ST391-0005	5mg
BML-ST391-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability

As indicated on product label or CoA when stored as recommended. U73122 is only sparingly soluble in aqueous media. Therefore the compounds are delivered to cells either after dissolving in organic solvents or complexing with serum proteins. Serum albumin (1-10mg/ml). Dissolve a weighed amount of U-73122 in methylene chloride or chloroform. Dispense in aliquots sufficient for one experiment into suitable vessels. Evaporate solvent with nitrogen. Stock solutions in DMSO or ethanol are stable for 2 months when stored at -20°C. Preferably dissolve just before use. Use caution in reusing stored DMSO solutions. Discard solutions that have turned to a pink color, which indicates a loss of inhibitor activity. Dried aliquots prepared from chloroform solutions are stable for up to 1 month when stored at -20°C. Useful range of concentrations for U-73122 is 0.1 to 10µM. The appropriate concentration will depend upon the cell type, cell density, medium concentration (the maleimide moiety of U-73122 may react with thiol containing compounds) and the particular cell parameter that is being measured.

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)-1H-pyrrole-2,5-dione

Appearance

white to off-white solid

CAS

112648-68-7

Couple Target

Lipoxygenase, Phospholipase, PLC

Couple Type

Inhibitor

Formula

$C_{29}H_{40}N_2O_3$

MW

464.7

Purity

≥98% (HPLC)



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

Soluble in methylene chloride or chloroform. Slightly soluble in DMSO (2mg/ml) or 100% ethanol (1mg/ml), sparingly soluble in aqueous solutions.

ENZO LIFE SCIENCES
& Luxembourg
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

ENZO LIFE SCIENCES
& Luxembourg
Phone: +33 472 440 655
info-fr@enzolifesciences.com

ENZO LIFE SCIENCES
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

ENZO LIFE SCIENCES
Phone: +44 1392 825900
info-uk@enzolifesciences.com

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