

U-74389G

Blocks lipid peroxidation

Lazaroid inhibitor of iron-dependent lipid peroxidation. Antioxidant. Inhibits lipid peroxidation in a variety of experimental systems including cultured brain microvessel endothelial cells, Cu^{2+} treated monocyclic THP1 cells and LLC-PK1 cell layers where it blocks F_2 isoprostane production. Prevented organ damage in intestinal cold storage, preservation and reperfusion injury.

Citations: 15

[View Online »](#)

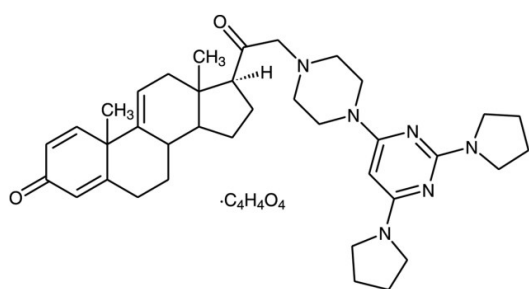
Ordering Information

[Order Online »](#)

BML-FR100-0100	100mg
----------------	-------

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, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	21-(4-(2,6-di-1-Pyrrolidinyl-4-pyrimidinyl)-1-piperazinyl)-pregna-1,4,9(11)-triene-3,20-dione . (Z)-2-butenedioate
Appearance	White solid.
CAS	153190-29-5
Formula	$C_{37}H_{50}N_6O_2 \cdot C_4H_4O_4$
MW	726.9
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (20mg/ml). Insoluble in water.

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