

IM-54

Necrosis inhibitor

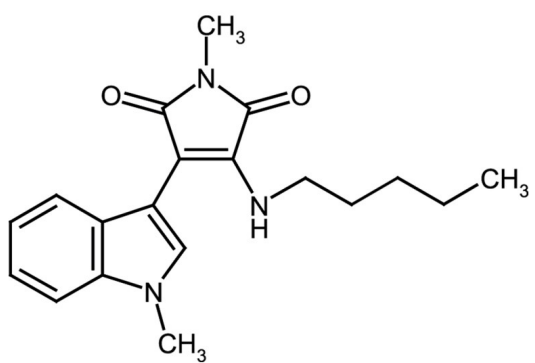
Cell permeable inhibitor of necrosis. Selectively blocks oxidative stress-induced necrotic cell death (~3 μ M IM-54 prevented ~50% cell death in HL60 cells exposed to 100 μ M H₂O₂). Does not protect against etoposide-induced apoptosis or display antioxidant properties. Does not affect S6K1 and PKC kinase activities at concentrations of 50 μ M.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2-(1H-Indol-3-yl)-3-pentylamino-maleimide
Appearance	Yellow powder.
CAS	861891-50-1
Formula	$C_{19}H_{23}N_3O_2$
Identity	Determined by NMR.
MW	325.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (100mg/ml), dichloromethane, 100% ethanol or methanol (10mg/ml).

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-be@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