

ENMD-1068

Proteinase-activated receptor ligand

Selective PAR-2 antagonist. Dose dependently inhibits calcium signaling in Lewis lung carcinoma cells stimulated with a soluble peptide PAR-2 agonist or trypsin, but not with a soluble PAR-1 agonist. Does not inhibit platelet aggregation stimulated by thrombin suggesting that it does not inhibit PAR-3 or PAR-4 signaling. Reduces TNF- α and IL-1 β production from rheumatoid arthritis synovial primary cell culture and reduces joint swelling in a carageenan/kaolin-injected mouse model.

Citations: 9

[View Online »](#)

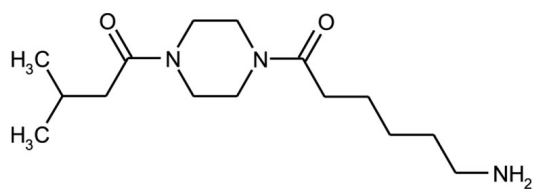
Ordering Information

[Order Online »](#)

BML-N110-0005	5mg
BML-N110-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White solid.
Couple Target	Proteinase-activated receptor
Couple Type	Ligand
Formula	$C_{15}H_{29}N_3O_2 \cdot HCl \cdot H_2O$
MW	283.3 . 36.5 . 18.0
Purity	≥98% (TLC)
Solubility	Soluble in water (5mg/ml), DMSO (2.5mg/ml) or 100% ethanol (5mg/ml).



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