

# 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- $\alpha$  and IL-1 $\beta$  production from rheumatoid arthritis synovial primary cell culture and reduces joint swelling in a carageenan/kaolin-injected mouse model.

Citations: 9

[View Online »](#)

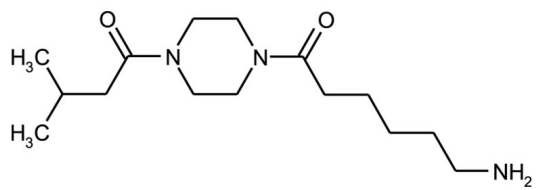
## Ordering Information

[Order Online »](#)

|               |     |
|---------------|-----|
| BML-N110-0005 | 5mg |
|---------------|-----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

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

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-be@enzolifesciences.com](mailto:info-be@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)