

# PR39 propeptide

A proline-arginine-rich peptide antibiotic. PR39 inhibits DNA and protein synthesis and the degradation of HIF-1 $\alpha$ . It has been reported to bind to the  $\alpha$ 7 subunit of the 26S proteasome, blocking degradation of the NK- $\kappa$ B inhibitor, I $\kappa$ B $\alpha$  by the ubiquitin-proteasome pathway without affecting overall proteasome activity. More recent studies have demonstrated it to be an efficient inhibitor of all activities of the 20S proteasome.

## Ordering Information

[Order Online »](#)

<b>BML-PW8850-0100</b>	100 $\mu$ g
------------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**MW** 4720.70

**Purity** ≥95%

**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)