

# 5'-Nucleotidase (from *Crotalus atrox* venom)

5'-nucleotidase (5'-ribonucleotide phosphohydrolase; EC 3.1.3.5) is partially purified from *Crotalus atrox* (Western Diamondback Rattlesnake) venom and is free of detectable phosphodiesterase activity.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI307-5000	5000KU
----------------	--------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 6 freeze-thaw cycles, but care should be taken that both freezing and thawing steps are done quickly (e.g. liquid N2 or dry ice/EtOH for freezing, rapid finger-rubbing for thawing). Frozen storage of smaller aliquots is recommended if more than 6 freeze-thaws of the entire vial contents (1000µl) are anticipated. Thawed enzyme is stable on ice for at least 2 hrs.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Activity** One U will release 1pmol phosphate per minute from 200µM 5'-AMP, 30°C in a reaction buffer of 10mM Tris-HCl, pH 7.4, 0.2mM MgCl<sub>2</sub>.

**Application Notes** 5'-nucleotidase is useful as the second enzyme in coupled assays of phosphodiesterases (3',5'-cyclic nucleotide phosphodiesterase; EC 3.1.4.17). Phosphodiesterase catalyzed hydrolysis of, for example, cAMP to 5'-AMP is coupled to 5'-nucleotidase cleavage of 5'-AMP to inorganic phosphate and adenosine. Phosphodiesterase activity is then quantified by colorimetric detection of the inorganic phosphate with a reagent such as BIOMOL Green (Prod. No. BML-AK111; see also Cyclic Nucleotide Phosphodiesterase Assay Kit, Prod. No. BML-AK800).

**Formulation** Liquid. In TRIS-HCl, pH 7.4, 150mM sodium chloride, 1mM magnesium chloride and 10% (v/v) glycerol.

**Quantity** 5000kU in 1000µl

**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)