

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

Use/Stability	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 N ₂ 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.
Long Term Storage	-80°C
Shipping	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 ₂ .
-----------------	--

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)