

# Visfatin/Nampt monoclonal antibody (P4D5AT)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-922-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for up to 12 months when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	PBEF, Pre-B-Cell colony enhancing factor, Nicotinamide phosphoribosyltransferase
<b>Application</b>	ELISA, WB
<b>Clone</b>	P4D5AT
<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 10% glycerol and 0.02% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human Visfatin (Nampt/PBEF) (aa 1-491).
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Protein G-affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Western Blot (1:1,000 to 1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	P43490

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)