

# Nampt (Visfatin/PBEF) (human), (recombinant)

Visfatin is an adipokine produced by visceral adipose tissue. Visfatin acts as a pro-inflammatory factor for peripheral blood mononuclear cells (PBMCs) and mimics the effects of insulin on adipocytes, monocytes, and hepatocytes.

## Ordering Information

[Order Online »](#)

ENZ-PRT265-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

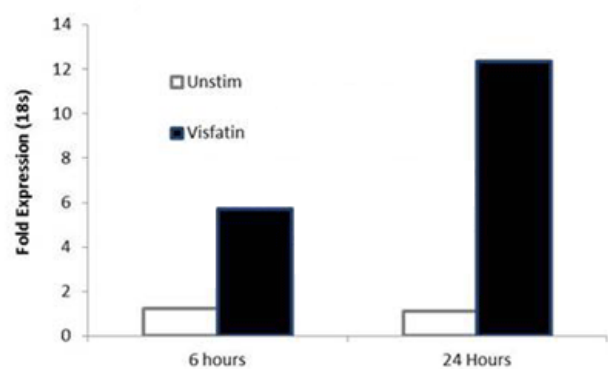


Figure 1: Human Nampt (Visfatin/PBEF) induced expression of IL-6 from human PBMCs.

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Pre-B-cell colony enhancing factor 1, Nicotinamide phosphoribosyltransferase
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% Trifluoroacetic Acid (TFA)
MW	~52.7 kDa
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute in sterile 20 mM HCl at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 466 amino acids.
UniProt ID	P43490



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)