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 »

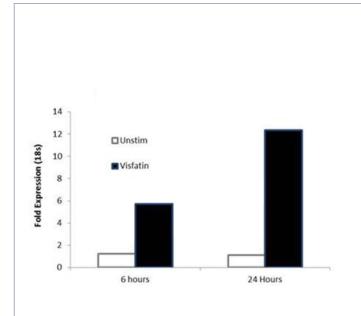


Figure 1: Human Nampt (Visfatin/PBEF) induced expression of IL-6 from human PBMCs.

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 *E. coli*. Contains 466 amino acids.

UniProt ID P43490

Last modified: May 29, 2024

