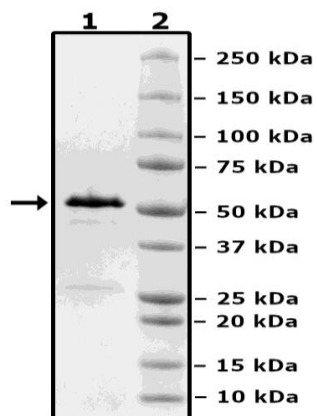


## Product Information

<b>Description:</b>	Human Nicotinamide phosphoribosyltransferase protein (NAMPT), also known as PBEF1 or visfatin (GenBank Accession No. NM_005746), a.a. 2-491 (end), no tag, MW=56 kDa, expressed in an <i>E. coli</i> cell expression system.
<b>Species</b>	Human
<b>Construct:</b>	NAMPT (PBEF1) (2-491(end))
<b>Concentration:</b>	0.86 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
<b>MW:</b>	56 kDa
<b>Genbank Accession:</b>	NM_005746
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Specific Activity:</b>	≥2 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to NAMPT Inhibitor Screening Assay Kit (BPS Bioscience #71276) with NAMPT titrated at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



NAMPT Activity

