



ELISA PRODUCT INFORMATION & MANUAL

Human PBEF/Visfatin/NAMPT ELISA Kit (Colorimetric) *NBP3-11683*

Enzyme-linked Immunosorbent Assay for quantitative
detection. For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

Table of Contents

1. Intended Use	3
2. Introduction	3
3. General References	4
4. Assay Principle	5
5. Handling & Storage	5
6. Kit Components	5
7. Materials Required but <i>Not</i> Supplied	6
8. General ELISA Protocol	7
8.1. Preparation and Storage of Reagents	7
8.2. Sample Collection, Storage and Dilution	8
8.3. Assay Procedure (Checklist)	9
9. Calculation of Results	10
10. Typical Data	10
11. Performance Characteristics	11-12
12. Technical Hints and Limitations	13
13. Troubleshooting	14
14. Notes	15
Product Specific References	Backcover

1. Intended Use

The Human PBEF/Visfatin/NAMPT ELISA Kit (Colorimetric) is to be used for the *in vitro* quantitative determination of human PBEF/Visfatin/NAMPT in cell lysates or cell-based assays (screening). This ELISA Kit is for research use only.

2. Introduction

Nampt, nicotinamide phosphoribosyl-transferase, is the rate-limiting enzyme of the mammalian NAD, nicotinamide adenine dinucleotide, biosynthesis pathway from nicotinamide to NMN, nicotinamide mononucleotide (1), which is further converted to NAD by nicotinamide/nicotinic acid mononucleotide adenytransferase abbreviated Nmnat. Nampt was originally identified in *Haemophilus ducreyi* (2), and was then found to have a significant homology to the mammalian pre-B cell colony-enhancing factor (PBEF) (3) or visfatin (4), which is also called extracellular Nampt (eNampt) (5) due to the fact that it is found as a circulating form in human serum and secreted from differentiated adipocytes. Although visfatin was originally reported as an insulin-mimicking hormone by its capability of binding to and activating the insulin receptor, subsequent studies do not support this observation (5, 6). However, Revollo et al. observed that haplodeficiency and chemical inhibition of Nampt caused defects in NAD biosynthesis and glucose-stimulated insulin secretion in pancreatic islets *in vivo* and *in vitro* and plasma visfatin and NMN levels were reduced in Nampt heterozygous females (5). This study proposed a model that NMN exists in high amounts in plasma, presumably derived from nicotinamide with help of eNampt. This circulating NMN and nicotinamide are uptaken by beta cells via unknown transport mechanism (s) are converted to NAD by Nmnat and intracellular Nampt (hereinafter abbreviated iNampt), respectively, concluding that Nampt-mediated systemic NAD biosynthesis is critical for insulin secretion presumably via a NAD-dependent histone deacetylase, Sirt1. Since it has been shown that NAD(+) levels in mitochondria remain at physiological levels following genotoxic stress and can maintain cell viability even when nuclear and cytoplasmic pools of NAD(+) are depleted (7), the NAD(+) biosynthetic enzyme Nampt plays a critical role in enhancing life span and protecting against oxidative cell damage. Therefore, measurement of iNampt located at different cell compartments like cytoplasm, mitochondria, and nucleus may be able to give us some biological clue on the intracellular functions of Nampt.

3. General References

- (1) The NAD biosynthesis pathway mediated by nicotinamide phosphoribosyltransferase regulates Sir2 activity in mammalian cells: J.R. Revollo, et al.; J. Biol. Chem. **279**, 50754 (2004)
- (2) Identification of a plasmid-encoded gene from *Haemophilus ducreyi* which confers NAD independence: P.R. Martin, et al.; J. Bacteriol. **183**, 1168 (2001)
- (3) Cloning and characterization of the cDNA encoding a novel human pre-B-cell colony-enhancing factor: B. Samal, et al.; Mol. Cell. Biol. **14**, 1431 (1994)
- (4) Visfatin: a protein secreted by visceral fat that mimics the effects of insulin: A. Fukuhara, et al.; Science **307**, 426 (2005)
- (5) Nampt/PBEF/Visfatin regulates insulin secretion in beta cells as a systemic NAD biosynthetic enzyme: J.R. Revollo, et al.; Cell. Metab. **6**, 363 (2007)
- (6) Molecular characteristics of serum visfatin and differential detection by immunoassays: A. Körner, et al.; J. Clin. Endocrinol. Metab. **92**, 4783 (2007)
- (7) Nutrient-sensitive mitochondrial NAD⁺ levels dictate cell survival: H. Yang, et al.; Cell **130**, 1095 (2007)

4. Assay Principle

This assay is a sandwich Enzyme Linked-Immunosorbent Assay (ELISA) for quantitative determination of human PBEF/Visfatin/NAMPT in cells. A monoclonal antibody specific for PBEF/Visfatin/NAMPT has been precoated onto the 96-well microtiter plate. Standards and samples are pipetted into the wells for binding to the coated antibody. After extensive washing to remove unbound compounds, PBEF/Visfatin/NAMPT is recognized by the addition of a purified polyclonal antibody specific for PBEF/Visfatin/NAMPT (Detection Antibody). After removal of excess polyclonal antibody, HRP conjugated anti-rabbit IgG (HRP) is added. Following a final washing, peroxidase activity is quantified using the substrate 3,3',5,5'-tetramethylbenzidine (TMB). The intensity of the color reaction is measured at 450nm after acidification and is directly proportional to the concentration of PBEF/Visfatin/NAMPT in the samples.

5. Handling & Storage

- Reagent must be stored at 2-8°C when not in use.
- Plate and reagents should be at room temperature before use.
- Do not expose reagents to temperatures greater than 25°C.

6. Kit Components

1 plate coated with human PBEF/Visfatin/NAMPT Antibody	(6 x 16-well strips)	
2 bottles Wash Buffer 10X	(2 x 30 ml)	(Wash Buffer 10X)
2 bottles ELISA Buffer 10X	(2 x 30 ml)	(ELISA Buffer 10X)
1 bottle Lysis Buffer 10X	(12 ml)	(LYSIS Buffer)
1 vial Detection Antibody	(60 µl)	(DET)
1 vial HRP 100X (HRP Conjugated anti-rabbit IgG)	(150 µl)	(HRP 100X)
1 vial human PBEF/Visfatin/NAMPT Standard (lyophilized) (STD)		(32 ng)
1 bottle TMB Substrate Solution	(12 ml)	(TMB)
1 bottle Stop Solution	(12 ml)	(STOP)
2 plate sealers (plastic film)		
2 silica Gel Minibags		

7. Materials Required but *Not* Supplied

- Microtiterplate reader at 450 nm
- Calibrated precision single and multi-channel pipettes. Disposable pipette tips
- Deionized water
- Microtubes or equivalent for preparing dilutions
- Disposable plastic containers for preparing working buffers
- Plate washer: automated or manual
- Glass or plastic tubes for diluting and aliquoting standard
- Phenyl methylsulfonyl fluoride (PMSF)

8. General ELISA Protocol

8.1. Preparation and Storage of Reagents

NOTE: Prepare just the appropriate amount of the buffers necessary for the assay.

- **Wash Buffer 10X** has to be diluted with deionized water 1:10 before use (e.g. 50 ml Wash Buffer 10X + 450 ml water) to obtain Wash Buffer 1X.
- **ELISA Buffer 10X** has to be diluted with deionized water 1:10 before use (e.g. 20 ml ELISA Buffer 10X + 180 ml water) to obtain ELISA Buffer 1X.
- **Lysis Buffer 10X** has to be diluted with deionized water 1:10 before use (e.g. 12 ml Lysis Buffer 10X + 108 ml water) to obtain Lysis Buffer 1X. Add 1 mM PMSF immediately before use.
- **Detection Antibody (DET)** has to be diluted to 1:250 in ELISA Buffer 1X (40 µl DET + 10 ml ELISA Buffer 1X).

NOTE: The diluted Detection Antibody is not stable and cannot be stored!

- **HRP 100X (HRP Conjugated anti-rabbit IgG)** has to be diluted to the working concentration by adding 100 µl in 10 ml of ELISA Buffer 1X (1:100).

NOTE: The diluted Detector is used within one hour of preparation.

- **Human PBEF/Visfatin/NAMPT Standard (STD)** has to be reconstituted with 1 ml of deionized water.
 - This reconstitution produces a stock solution of 32 ng/ml. Mix the standard to ensure complete reconstitution and allow the standard to sit for a minimum of 15 minutes. Mix well prior to making dilutions.

NOTE: The reconstituted standard is aliquoted and stored at -20°C.

- Dilute the standard protein concentrate (STD) (**32 ng/ml**) in ELISA Buffer 1X. A seven-point standard curve using 2-fold serial dilutions in ELISA Buffer 1X is recommended.
- Suggested standard points are:
16 , 8 , 4 , 2 , 1 , 0.5 , 0.25 and 0 ng/ml.

Dilute further for the standard curve:

To obtain	Add	Into
16 ng/ml	300 µl of PBEF/Visfatin/NAMPT (32 ng/ml)	300 µl of ELISA Buffer 1X
8 ng/ml	300 µl of PBEF/Visfatin/NAMPT (16 ng/ml)	300 µl of ELISA Buffer 1X
4 ng/ml	300 µl of PBEF/Visfatin/NAMPT (8 ng/ml)	300 µl of ELISA Buffer 1X
2 ng/ml	300 µl of PBEF/Visfatin/NAMPT (4 ng/ml)	300 µl of ELISA Buffer 1X
1 ng/ml	300 µl of PBEF/Visfatin/NAMPT (2 ng/ml)	300 µl of ELISA Buffer 1X
0.5 ng/ml	300 µl of PBEF/Visfatin/NAMPT (1 ng/ml)	300 µl of ELISA Buffer 1X
0.25 ng/ml	300 µl of PBEF/Visfatin/NAMPT (0.5 ng/ml)	300 µl of ELISA Buffer 1X
0 ng/ml	300 µl of ELISA Buffer 1X	Empty tube

8.2. Sample Collection, Storage and Dilution

Cell Lysates : Grow cell until 90% confluency. Scrap cells off the plate and transfer to an appropriate tube. Keep on ice and microcentrifuge at 1,200 rpm for 5 minutes at 4°C. Remove supernatant, rinse cells once with ice-cold PBS. Remove PBS and add 200 µl ice-cold **Lysis Buffer 1X** supplemented with 1 mM phenyl methylsulfonyl fluoride (PMSF) to ten million cells of interest and incubate on ice for 30 minutes. Microcentrifuge at 12,000 rpm for 5 minutes at 4°C and transfer the supernatant to a new tube. The supernatant is the cell lysate. Use freshly prepared cell lysate samples.

Cell Lysates have to be diluted in ELISA Buffer 1X. Samples containing visible precipitates must be clarified before use.

NOTE: As a starting point, 1/10 to 1/1,000 dilutions of cell lysates are recommended! If sample values fall outside the detection range of the assay, a lower or higher dilution may be required!

8.3. Assay Procedure (Checklist)

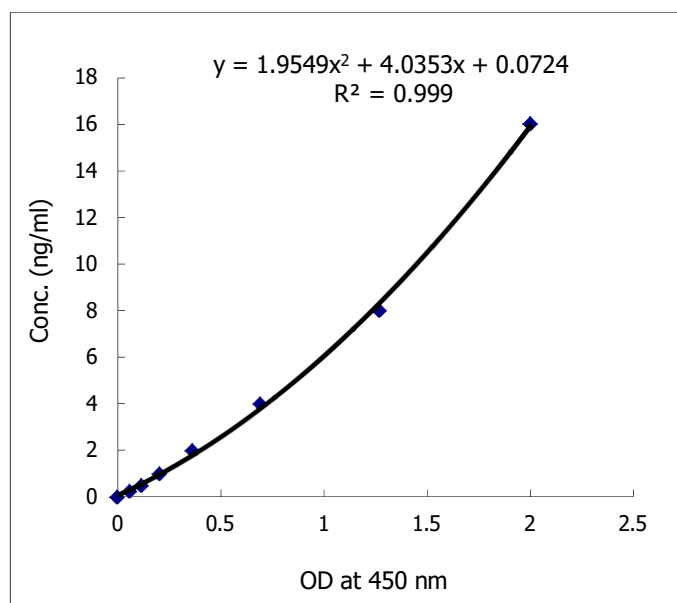
☐	1. Determine the number of 16-well strips needed for the assay and insert them in the frame for current use. The extra strips should be resealed in the foil pouch bag and stored at 4°C. NOTE: Remaining 16-well strips coated with PBEF/Visfatin/NAMPT antibody when opened can be stored at 4°C for up to 1 month.
☐	2. Add 100 µl of the different standards into the appropriate wells in duplicate! At the same time, add 100 µl of diluted cell lysates samples in duplicate to the wells (see 8.1. Preparation and Storage of Reagents and 8.2. Preparation of Samples).
☐	3. Cover the plate with plate sealer and incubate for 1 hour at 37°C.
☐	4. Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of three washes. After the last wash, complete removal of liquid is essential for good performance.
☐	5. Add 100 µl to each well of the Detection Antibody (DET). (see 8.1. Preparation and Storage of Reagents).
☐	6. Cover the plate with plate sealer and incubate for 1 hour at 37°C.
☐	7. Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of three washes. After the last wash, complete removal of liquid is essential for good performance.
☐	8. Add 100 µl to each well of the diluted HRP Conjugated anti-rabbit IgG (HRP) (see 8.1. Preparation and Storage of Reagents).
☐	9. Cover the plate with plate sealer and incubate for 1 hour at 37°C.
☐	10. Aspirate the coated wells and add 300 µl of Wash Buffer 1X using a multichannel pipette or auto-washer. Repeat the process for a total of five washes. After the last wash, complete removal of liquid is essential for good performance.
☐	11. Add 100 µl to each well of TMB Substrate Solution (TMB).
☐	12. Allow the color reaction to develop at room temperature (RT°C) in the dark for 10 minutes.
☐	13. Stop the reaction by adding 100 µl of Stop Solution (STOP). Tap the plate gently to ensure thorough mixing. The substrate reaction yields a blue solution that turns yellow when Stop Solution (STOP) is added.
	! CAUTION: CORROSIVE SOLUTION!
☐	14. Measure the OD at 450 nm in an ELISA reader within 30 minutes.

9. Calculation of Results

- Average the duplicate readings for each standard, control and sample and subtract the average blank value (obtained with the 0 ng/ml point).
- Generate the standard curve by plotting the average absorbance obtained for each standard concentration on the horizontal (X) axis vs. the corresponding PBEF/Visfatin/NAMPT concentration (ng/ml) on the vertical (Y) axis (see **10. TYPICAL DATA**).
- Calculate the PBEF/Visfatin/NAMPT concentrations of samples by interpolation of the regression curve formula as shown above in a form of a quadratic equation
- If the test samples were diluted, multiply the interpolated values by the dilution factor to calculate the concentration of human PBEF/Visfatin/NAMPT in the samples.

10. Typical Data

The following data are obtained using the different concentrations of standard as described in this protocol:



Standard PBEF/Visfatin/NAMPT (ng/ml)	Optical Density (mean)
16	1.953
8	1.124
4	0.612
2	0.306
1	0.167
0.5	0.097
0.25	0.045
0	0

Figure: Standard curve

11. Performance Characteristics

A. Sensitivity (Limit of detection):

The lowest level of PBEF/Visfatin/NAMPT that can be detected by this assay is 30 pg/ml. **NOTE:** *The Limit of detection was measured by adding two standard deviations to the mean value of 50 zero standard.*

B. Assay range: 0.25 ng/ml – 16 ng/ml

C. Specificity:

This ELISA is specific for the measurement of natural and recombinant human PBEF/Visfatin/NAMPT. It does not cross-react with human adiponectin, human resistin, human RELM- β , human leptin, human GPX3, human ANGPTL4, human FABP4, human ANGPTL6, human PAI1.

Mouse PBEF/Visfatin/NAMPT shows weakly 5% cross-reactivity

in this assay. Rat PBEF/Visfatin/NAMPT shows weakly 15%

cross-reactivity in this assay.

D. Intra-assay precision:

Two samples of known concentrations of human PBEF/Visfatin/NAMPT were assayed in replicates 10 times to test precision within an assay.

Samples	Means (ng/ml)	SD	CV (%)	n
1	244.46	6.66	2.73	10
2	248.13	24.21	9.76	10

E. Inter-assay precision:

Two samples of known concentrations of human PBEF/Visfatin/NAMPT were assayed in 6 separate assays to test precision between assays.

Samples	Means (ng/ml)	SD	CV (%)	n
1	209.59	8.42	4.02	6
2	251.04	18.58	7.40	6

F. Recovery:

When samples (cell lysates) are spiked with known concentrations of human PBEF/Visfatin/NAMPT, the recovery averages 98% (range from 90% to 105%).

Samples	Average recovery (%)	Range (%)
1	96.36	95-105
2	102.62	95-105

G. Linearity:

Different human cell lysates samples containing PBEF/Visfatin/NAMPT were diluted several fold (1/100 to 1/3,200) and the measured recoveries ranged from 90% to 105%.

Samples	Sample Dilution	Expected (ng/ml)	Observed (ng/ml)	% of Expected
293 cells	1 : 100	244.04	244.04	100
	1 : 200	122.02	110.55	90.60
	1 : 400	61.01	57.20	93.76
PBEF/Visfat in/NAMPT over-expressed 293 cells	1 : 800	5161.67	5161.67	100
	1 : 1,600	2580.83	2640.78	102.32
	1 : 3,200	1290.42	1207.36	93.56

12. Technical Hints and Limitations

- It is recommended that all standards, controls and samples be run in duplicate.
- Do not combine leftover reagents with those reserved for additional wells.
- Reagents from the kit with a volume less than 100 µl should be centrifuged.
- Residual wash liquid should be drained from the wells after last wash by tapping the plate on absorbent paper.
- Crystals could appear in the 10X solution due to high salt concentration in the stock solutions. Crystals are readily dissolved at room temperature or at 37°C before dilution of the buffer solutions.
- Once reagents have been added to the 16-well strips, DO NOT let the strips DRY at any time during the assay.
- Keep TMB Substrate Solution (TMB) protected from light.
- The Stop Solution (STOP) consists of sulfuric acid. Although diluted, the Stop Solution (STOP) should be handled with gloves, eye protection and protective clothing.

13. Troubleshooting

PROBLEM	POSSIBLE CAUSES	SOLUTIONS
No signal or weak signal	Omission of key reagent	Check that all reagents have been added in the correct order.
	Washes too stringent	Use an automated plate washer if possible.
	Incubation times inadequate	Incubation times should be followed as indicated in the manual.
	Plate reader settings not optimal	Verify the wavelength and filter setting in the plate reader.
	Incorrect assay temperature	Use recommended incubation temperature. Bring substrates to room temperature before use.
High background	Concentration of HRP too high	Use recommended dilution factor.
	Inadequate washing	Ensure all wells are filling wash buffer and are aspirated completely.
Poor standard curve	Wells not completely aspirated	Completely aspirate wells between steps.
	Reagents poorly mixed	Be sure that reagents are thoroughly mixed.
Unexpected results	Omission of reagents	Be sure that reagents were prepared correctly and added in the correct order.
	Dilution error	Check pipetting technique and double-check calculations.

14. Notes