



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
NAMPT Inhibitor Screening Assay Kit
Catalog: 71276-2
Size: 384 reactions

BACKGROUND: NAMPT catalyzes the formation of nicotinamide mononucleotide from nicotinamide and 5-phosphoribosyl-1-pyrophosphate (PRPP). It is the rate limiting component in the mammalian NAD biosynthesis pathway. The protein is thought to be involved in many important biological processes, including metabolism, stress response and aging.

DESCRIPTION: The *NAMPT Inhibitor Screening Assay Kit* is designed to measure NAMPT activity for screening and profiling applications. The NAMPT assay kit comes in a convenient 384-well format, with purified recombinant NAMPT enzyme, NAMPT assay buffer, NAMPT dilution buffer, ATP, Nicotinamide, PRPP, and Ethanol enough for 384 enzyme reactions.

COMPONENTS:

Catalog #	Reagent	Amount	Storage	
91004	NAMPT	60 µg	-80°C	<i>Avoid multiple freeze/thaw cycles!</i>
	NAMPT dilution buffer	3 ml	-80°C	
	4x NAMPT assay buffer	750 µl x3	-80°C	
79686	ATP (400 µM)	250 µl x2	-80°C	
	Nicotinamide (400 µM)	250 µl x2	-80°C	
	PRPP (800 µM)	250 µl x2	-80°C	
	Ethanol (30%)	1 ml	-80°C	
	Low binding black microtiter 384-well plate	1	Room Temp.	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader
Adjustable micropipettor and sterile tips
Aluminum foil

APPLICATIONS: Great for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

STABILITY: Up to 6 months from date of receipt, when stored as recommended.

REFERENCE: Ramsey, K.M., et al. (2009) *Science* **324**: 651-654
Nakahata, Y., et al. (2009) *Science* **324**: 654-657.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

1. Thaw **4x NAMPT assay buffer, ATP, Nicotinamide, Phosphoribosyl pyrophosphate (PRPP), 30% Ethanol.**
2. Thaw **NAMPT** enzyme on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Calculate the amount of **NAMPT** required for the assay and dilute enzyme to 12-25 ng/ μ l with **NAMPT dilution buffer**. Aliquot remaining **NAMPT** enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C . **NAMPT enzyme is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.**
3. Add 6 μ l of diluted **NAMPT** enzyme to the wells designated "Positive Control" and "Test Inhibitor". To the wells designated as "Blank," add 6 μ l **NAMPT dilution buffer**.
4. Add 4 μ l of Inhibitor solution of each well labeled as "Test Inhibitor." For the wells labeled "Positive Control" and "Blank," add 4 μ l of the same solution without inhibitor (Inhibitor buffer). Note: Final DMSO concentration must be $\leq 1\%$. Higher DMSO levels can significantly decrease the enzyme activity. For example, to test an inhibitor dissolved in 100% DMSO at 10 μ M, dilute 1 mM inhibitor with water to make a 50 μ M inhibitor in 5% DMSO (aq). Then, add 4 μ l of the 50 μ M solution into the 20 μ l assay to make a 1% DMSO concentration in the final reaction mixture.
5. Preincubate **NAMPT** with the inhibitors for 30 min at room temperature (slow shaking).
6. During the enzyme preincubation, prepare the master mixture (10 μ l per well): N wells x (5 μ l **4x NAMPT assay buffer** + 1 μ l **ATP** (400 μ M) + 1 μ l **Nicotinamide** (400 μ M) + 1 μ l **Phosphoribosyl pyrophosphate (PRPP)**, 800 μ M) + 1 μ l 30% **Ethanol** + 1 μ l **water**). Add 10 μ l to every well. Store the remaining components at -80°C in single use aliquots.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

	Positive Control	Test Inhibitor	Blank
4x NAMPT assay buffer	5 µl	5 µl	5 µl
ATP (400 µM)	1 µl	1 µl	1 µl
Nicotinamide (400 µM)	1 µl	1 µl	1 µl
PRPP (800 µM)	1 µl	1 µl	1 µl
30% Ethanol	1 µl	1 µl	1 µl
Water	1 µl	1 µl	1 µl
Test Inhibitor	-	4 µl	-
5% DMSO in Water (Inhibitor buffer)	4 µl	-	4 µl
NAMPT dilution buffer	-	-	6 µl
NAMPT (12-25 ng/µl)	6 µl	6 µl	-
Total	20 µl	20 µl	20 µl

7. Initiate the reaction by adding 10 µl of **Master Mixture** to all the wells. Incubate at 30°C for two hours.
8. After the reaction, read sample in a microtiter-plate reading fluorimeter capable of excitation at a wavelength of ~340 nm and detection of emitted light of ~460 nm. "Blank" value is subtracted from all other values.

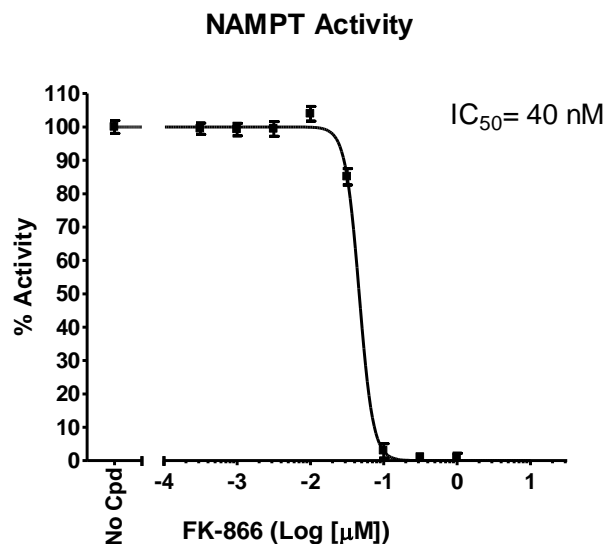
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Example of Assay Results:



NAMPT inhibition by FK866, measured using the NAMPT Inhibitor Screening Assay Kit, BPS Bioscience Cat. #79307. The compound was pre-incubated with the NAMPT enzyme for 30 min at room temperature before the reaction was initiated with the addition of master mix. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
NAMPT (PBEF1), GST tag	91004	50 µg
NAMPT (PBEF1)	71098	50 µg
NMNAT, His-tag	71090	100 µg
NAMPT Inhibitor Screening Assay Kit	71726	96 rxns
CD38, His-Tag (Human), HiP™	71227	100 µg
CD38 Inhibitor Screening Assay Kit	71275	96 rxns
5'-Nucleotidase/CD73, His-tag	71184	50 µg
CD73 Inhibitor Screening Assay Kit	72055	96 rxns
CD73 Inhibitor Screening Assay Kit	72058	384 rxns
Adenosine Deaminase (ADA), His-tag	70016	100 µg
NAD+, Biotin-Labeled	80610	500 µl

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com