

NAMPT Inhibitor Screening Assay Kit

Description

The NAMPT Inhibitor Screening Assay Kit is a 96-well fluorescent based assay designed to measure the activity of NAMPT (nicotinamide phosphoribosyltransferase) for screening and profiling applications. The kit contains enough purified NAMPT enzyme, NAMPT Assay Buffer, NAMPT Dilution Buffer, ATP, Nicotinamide, PRPP and ethanol for 100 enzyme reactions.

Background

Nicotinamide phosphoribosyltransferase (NAMPT) catalyzes the formation of nicotinamide mononucleotide from nicotinamide and 5-phosphoribosyl-1-pyrophosphate (PRPP). It is the rate limiting enzyme in the mammalian NAD (nicotinamide adenine dinucleotide) biosynthesis pathway. NAMPT is thought to be involved in many important biological processes, including metabolism, stress response and aging. NAMPT is often overexpressed in cancer cells, an adaptation required to sustain the higher energetic need of the cells. In addition, it also plays roles in proliferation, invasion, stemness, angiogenesis, immune regulation, and drug resistance of tumors, making it an attractive target candidate for cancer therapy.

Applications

Study enzyme kinetics and screen small molecule inhibitors for drug discovery and High Throughput Screen (HTS) applications.

Supplied Materials

Catalog #	Name	Amount	Storage
91004	Human NAMPT (PBEF1), GST-Tag*	50 µg	-80°C
	NAMPT Dilution Buffer	3 ml	-80°C
79946	4x NAMPT Assay Buffer	2 x 750 µl	-80°C
79686	400 µM ATP	250 µl	-80°C
	400 µM Nicotinamide	250 µl	-80°C
	800 µM PRPP	250 µl	-80°C
	30% Ethanol	1 ml	-80°C
79685	Low binding black microtiter 96-well plate	1	Room Temp.

* The concentration of protein is lot-specific and will be indicated on the tube containing the protein.

Materials Required but Not Supplied

- Fluorescent microplate reader capable of reading $\lambda_{exc}/\lambda_{em}$ = 340 nm/460 nm
- Adjustable micropipettor and sterile tips

Stability



This assay kit will perform optimally for up to **6 months** from date of receipt when the materials are stored as directed.

Safety

This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous and is harmful by inhalation, in contact with skin, eyes, clothing, and if swallowed. If contact occurs, wash thoroughly.

Contraindications

The final concentration of DMSO in the assay should not exceed 1%.

Assay Protocol

- All samples and controls should be tested in duplicate.
 - The assay should include “Blank”, “Positive Control” and “Test Inhibitor” conditions.
 - We recommend maintaining the diluted protein on ice during use.
 - For detailed information on protein handling please refer to Protein FAQs (bpsbioscience.com).
 - We recommend using FK-866 as internal control. If not running a dose response curve for the control inhibitor, we recommend running the control inhibitor at 0.1X, 1X and 10X the IC₅₀ value shown in the validation data below.
1. Thaw **4x NAMPT Assay Buffer, 400 μM ATP, 400 μM Nicotinamide, 800 μM PRPP, NAMPT Dilution Buffer** and **30% Ethanol** on ice.
 2. Thaw **NAMPT** enzyme on ice. Briefly spin the tube to recover the full content of the tube.
 3. Dilute NAMPT with NAMPT Dilution Buffer to 20-50 ng/μl (10 μl/well).
 4. Add 10 μl of diluted NAMPT to the “Positive Control” and “Test Inhibitor” wells.
 5. Add 10 μl of NAMPT Dilution Buffer to the “Blank” wells.
 6. Prepare Test Inhibitor (5 μl/well): for a titration prepare serial dilutions at concentrations 10-fold higher than the desired final concentrations. The final volume of the reaction is 50 μl.
 - 6.1. If the Test Inhibitor is soluble in water, make a dilution in distilled water at a concentration 10-fold higher than the final desired concentration.

For the positive and negative controls, use distilled water (Diluent Solution).

OR

6.2. If the Test Inhibitor is soluble in DMSO: Prepare the test inhibitor at 100-fold the highest desired concentration in 100% DMSO, then dilute the inhibitor 10-fold in distilled water to prepare the highest concentration of the 10-fold intermediate dilutions. The concentration of DMSO is now 10%.

Prepare serial dilutions of the Test Inhibitor at 10-fold the desired final concentrations using 10% DMSO in distilled water to keep the concentration of DMSO constant.

For positive and negative controls, prepare 10% DMSO in water so that all wells contain the same amount of DMSO (Diluent Solution).

Note: The final concentration of DMSO in the assay should not exceed 1%.

7. Add 5 µl of Test Inhibitor to the “Test Inhibitor” wells.
8. Add 5 µl of Diluent Solution to the “Positive Control” and “Blank” wells.
9. Pre-incubate NAMPT with the inhibitors for 30 minutes at Room Temperature (RT), with gentle agitation.
10. Prepare a Master Mix (35 µl/well): N wells x (12.5 µl of 4x NAMPT Assay Buffer + 2.5 µl of 400 µM ATP + 2.5 µl of 400 µM Nicotinamide + 2.5 µl of 800 µM PRPP + 2.5 µl of 30% Ethanol + 12.5 µl of distilled water).
11. Start the reaction by adding 35 µl of Master Mix to all wells.
12. Incubate at 30°C for 2 hours.

Component	Blank	Test Inhibitor	Positive Control
Master Mix	35 µl	35 µl	35 µl
Test Inhibitor	-	5 µl	-
Diluent Solution	5 µl	-	5 µl
NAMPT Dilution Buffer	10 µl	-	-
NAMPT (20-50 ng/µl)	-	10 µl	10 µl
Total	50 µl	50 µl	50 µl

13. Measure the fluorescence intensity in a microtiter plate-reader fluorimeter capable of excitation at $\lambda=340$ nm and emission at $\lambda=460$ nm.
14. The “Blank” should be subtracted from all other values (background value).

Example Results

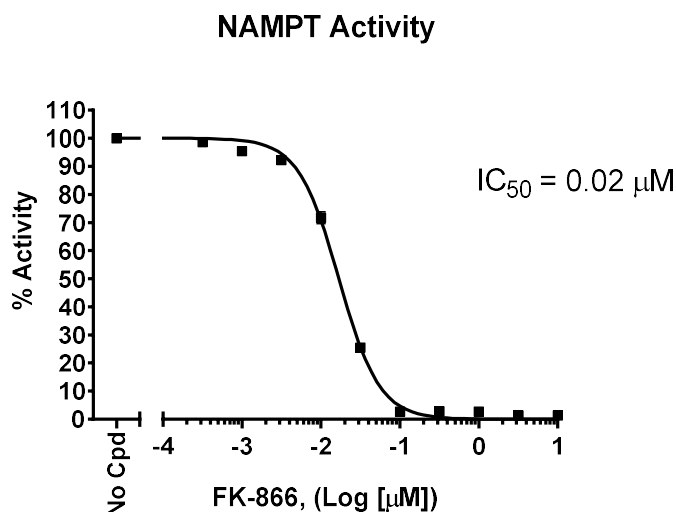


Figure 1: NAMPT inhibition by FK-866.

NAMPT activity was measured in the presence of increasing concentrations of FK-866. Results are expressed as percent of control activity (measured in the absence of inhibitor and set at 100%).

Data shown is representative. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com

Troubleshooting Guide

Visit bpsbioscience.com/assay-kits-faq for detailed troubleshooting instructions. For all further questions, please email support@bpsbioscience.com

References

Ramsey K., *et al.*, 2009, *Science* 324 (5927): 651-4.
 Nakahata Y., *et al.*, 2009, *Science* 324 (5927) : 654-7.
 Gasparini M., *et al.*, 2022, *Int J Biochem Cell Biol.* 145 : 106189.

Related Products

Products	Catalog #	Size
NAMPT (PBEF1)	71098	20 μ g
NMNAT1, His-Tag	71090	100 μ g
CHS-828	27333	25 mg
CD38, His-Tag (Human), HiP™	71277	100 μ g
CD38 Inhibitor Screening Assay Kit (Cyclase Activity)	71275	96 reactions
CD73, His-Tag (Human)	71184	50 μ g/500 μ g

Version 020724