

Product Information

Purine Nucleoside Phosphorylase Activity Assay Kit (Colorimetric)

Catalog Number **MAK289**

Storage Temperature -20 °C

TECHNICAL BULLETIN

Product Description

Purine Nucleoside Phosphorylase (PNP, E.C. 2.4.2.1.) is an enzyme involved in purine metabolism. PNP catalyzes the cleavage of the glycosidic bond of ribo- or deoxyribonucleosides, in the presence of inorganic phosphate as a second substrate, to generate the purine base, and ribose-1-phosphate or deoxyribose-1-phosphate. It is one of the enzymes of the nucleotide salvage pathways that allows the cell to produce nucleotide monophosphates when the *de novo* synthesis pathway has been interrupted or is non-existent (as is the case in the brain).

Purine Nucleoside Phosphorylase is a cytosolic enzyme. PNP deficiency disease leads to toxic buildup of deoxyguanosine in T-cells leading to T-lymphocytopenia with no apparent effects on B-lymphocyte function. Inhibition of PNP is an important target for chemotherapeutic applications and treatment of T-cell mediated autoimmune diseases. PNP deficiency is also associated with neurological problems.

In this Purine Nucleoside Phosphorylase Activity Assay, hypoxanthine formed by the breakdown of inosine is further converted to uric acid using a developer. The uric acid is measured at a wavelength of 293 nm. Limit of quantification is 0.1 µU recombinant Purine Nucleoside Phosphorylase.

Unit Definition: One unit of Purine Nucleoside Phosphorylase Activity is the amount of enzyme that hydrolyzes inosine to yield 1.0 µmole of hypoxanthine per minute at room temperature.

Components

The kit is sufficient for 100 assays in 96 well plates.

PNP Assay Buffer (10×) Catalog Number MAK289A	10 mL
Developer Catalog Number MAK289B	1 vial
Inosine Substrate Catalog Number MAK289C	200 µL
Hypoxanthine Standard (10 mM) Catalog Number MAK289D	100 µL
PNP Positive Control Catalog Number MAK289E	1 vial
UV Transparent Plate (96 well) Catalog Number MAK289F	1 each

Reagents and Equipment Required but Not Provided.

- 96 well flat-bottom plate – Use provided plate for this UV assay.
- Spectrophotometric multiwell plate reader
- Protease Inhibitor Cocktail
- Dounce Homogenizer

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Briefly centrifuge vials before opening. Use ultrapure water for the preparation of reagents. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

PNP Assay Buffer (10×) – Make 1× buffer by adding one volume of 10× Assay Buffer to nine volume of water. Store at -20°C or 4°C . Bring to room temperature before use.

Developer – Reconstitute with 210 μL of 1× PNP Assay Buffer and mix gently by pipetting. Briefly centrifuge to collect the contents in the bottom of the tube. Aliquot and store at -20°C . Avoid repeated freeze/thaw cycles.

Inosine Substrate – Aliquot and store at -20°C . Avoid repeated freeze/thaw cycles.

PNP Positive Control – Reconstitute with 22 μL of 1× PNP Assay Buffer and mix gently by pipetting. Briefly centrifuge to collect the contents in the bottom of the tube. Aliquot and store at -20°C . Avoid repeated freeze/thaw cycles.

Storage/Stability

The kit is shipped on wet ice and storage at -20°C , protected from light, is recommended.

Procedure

All samples and standards should be run in duplicate.

Sample Preparation

Rinse tissue and transfer ~100 mg of fresh or frozen tissue (stored at -80°C) to a pre-chilled homogenizer. Add 300 μL of cold 1× PNP Assay Buffer containing protease inhibitor cocktail (not provided) and thoroughly homogenize tissue on ice. Transfer the tissue homogenate to a cold microfuge tube.

To prepare cell extract, add 150–300 μL of cold 1× PNP Assay Buffer containing protease inhibitor cocktail (not provided) to $1-5 \times 10^6$ fresh or frozen cells and pipette several times to disrupt the cells. Transfer cell homogenate including cell debris to a cold microfuge tube and agitate on a rotary shaker at 4°C for at least 15 minutes.

Centrifuge the tissue or cell homogenate at $10,000 \times g$ at 4°C for 15 minutes. Transfer the clarified supernatant to a fresh pre-chilled tube and store on ice. Use lysates immediately to assay PNP activity.

Note: Lysates can be aliquoted and snap frozen in liquid nitrogen before storing at -80°C . Avoid freeze/thaw cycles.

Hypoxanthine Standard

Dilute Hypoxanthine Standard to 1 mM by adding 10 μL of 10 mM Hypoxanthine Standard to 90 μL of 1× PNP Assay Buffer. Add 0, 2, 4, 6, 8, and 10 μL of prepared 1 mM Hypoxanthine Standard into a series of wells in the provided 96 well plate to generate 0, 2, 4, 6, 8, and 10 nmole/well Hypoxanthine Standard. Adjust the volume to 50 μL /well with 1× PNP Assay Buffer.

Purine Nucleoside Phosphorylase Activity Assay

Add 2–50 μL of sample into desired well(s) in the provided 96 well plate. For Positive Control, add 2 μL of Positive Control for the assay. Make up the volume of samples and Positive Control to 50 μL /well with 1× PNP Assay Buffer. Add 50 μL of 1× PNP Assay Buffer to one well as reagent Background Control.

Notes: For unknown samples, performing a pilot experiment and testing several sample concentrations to ensure the readings are within the Standard Curve range is suggested.

Small molecules such as xanthine and hypoxanthine in the samples will contribute to the background. If the background level is too high, remove these molecules by passing the sample through a desalting column or by buffer exchange using a 10 kDa spin column. Use this treated sample for the assay.

Optional: Prepare a parallel sample well as sample background control to ensure that the small molecules are removed by either using a desalting or spin column.

Sample Mixes

Prepare enough reagents for the number of assays to be performed. Make 50 μL of the appropriate mix for each well, see Table 1.

Table 1.

Preparation of Sample Mixes

Reagent	Reaction Mix	Background Control Mix
1× PNP Assay Buffer	46 μL	48 μL
Developer	2 μL	2 μL
Inosine Substrate	2 μL	–

Add 50 μL of Reaction Mix into each sample, reagent background control, and Positive Control wells, and 50 μL of Background Control mix to Standards and sample background control well(s). Mix well.

Measurement

Measure absorbance (A_{293}) in kinetic mode for at least 30 minutes at room temperature. Choose two time points (T_1 and T_2) in linear range (can be as short as 2 minutes) of plot and obtain corresponding absorbance for sample (A_{S1} and A_{S2}) and reagent background control (A_{BG1} and A_{BG2}). Read the Hypoxanthine Standard Curve along with the samples.

ResultsCalculations

Subtract 0 Standard reading from all Standard Readings. Plot the Hypoxanthine Standard Curve. Subtract reagent background control reading from sample reading. Compare the ΔA_{293} [$(A_{S2} - A_{BG2}) - (A_{S1} - A_{BG1})$] to the Standard Curve to obtain B (nmole) of Hypoxanthine generated by the sample during the reaction time ($\Delta T = T_2 - T_1$).

$$\text{PNP activity} = \frac{B}{(\Delta T \times \mu\text{g})} \times \text{DF}$$

$$(\text{nmole}/\text{min}/\mu\text{g} = \text{mU}/\mu\text{g})$$

B = Hypoxanthine amount from Standard Curve (pmole)

ΔT = the reaction time (minutes)

μg = the amount of protein/well (μg)

DF = is the dilution factor of the sample

Sample PNP Activity can also be expressed as U/mg ($\mu\text{moles}/\text{min}$ hypoxanthine generated per mg) of protein.

Unit Definition: One unit of Purine Nucleoside Phosphorylase Activity is the amount of enzyme that hydrolyzes inosine to yield 1.0 μmole of hypoxanthine per minute at room temperature.

Troubleshooting Guide

Problem	Possible Cause	Suggested Solution
Assay Not Working	Cold assay buffer	Assay Buffer must be at room temperature
	Omission of step in procedure	Refer and follow Technical Bulletin precisely
	Plate reader at incorrect wavelength	Check filter settings of instrument
	Type of 96 well plate used	Use provided plate for this UV assay.
Samples with erratic readings	Samples prepared in different buffer	Use the Assay Buffer provided or refer to Technical Bulletin for instructions
	Cell/Tissue culture samples were incompletely homogenized	Repeat the sample homogenization, increasing the length and extent of homogenization step.
	Samples used after multiple freeze-thaw cycles	Aliquot and freeze samples if samples will be used multiple times
	Presence of interfering substance in the sample	If possible, dilute sample further
	Use of old or inappropriately stored samples	Use fresh samples and store correctly until use
Lower/higher readings in samples and standards	Improperly thawed components	Thaw all components completely and mix gently before use
	Use of expired kit or improperly stored reagents	Check the expiration date and store the components appropriately
	Allowing the reagents to sit for extended times on ice	Prepare fresh Reaction Mix before each use
	Incorrect incubation times or temperatures	Refer to Technical Bulletin and verify correct incubation times and temperatures
	Incorrect volumes used	Use calibrated pipettes and aliquot correctly
Non-linear standard curve	Use of partially thawed components	Thaw and resuspend all components before preparing the reaction mix
	Pipetting errors in preparation of standards	Avoid pipetting small volumes
	Pipetting errors in the Reaction Mix	Prepare a Reaction Mix whenever possible
	Air bubbles formed in well	Pipette gently against the wall of the plate well
	Standard stock is at incorrect concentration	Refer to the standard dilution instructions in the Technical Bulletin
	Calculation errors	Recheck calculations after referring to Technical Bulletin
	Substituting reagents from older kits/lots	Use fresh components from the same kit
Unanticipated results	Samples measured at incorrect wavelength	Check the equipment and filter settings
	Samples contain interfering substances	If possible, dilute sample further
	Sample readings above/below the linear range	Concentrate or dilute samples so readings are in the linear range

SJ,MAM 04/16-1