

Product datasheet

Purine Nucleoside Phosphorylase Activity Assay Kit (Fluorometric) ab204706

[3 Images](#)

Overview

Product name	Purine Nucleoside Phosphorylase Activity Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Cell Lysate, Purified protein, Tissue Lysate
Assay type	Enzyme activity
Sensitivity	0.005 μ U
Species reactivity	Reacts with: Other species, Mammals
Product overview	Purine Nucleoside Phosphorylase Activity Assay Kit (Fluorometric) (ab204706) is an assay where hypoxanthine formed from the breakdown of inosine is detected via a multi-step reaction, resulting in the generation of an intermediate that reacts with the PNP Probe. The fluorescent product is measured at Ex/Em = 535/587 nm. Limit of quantification is 0.005 μ U recombinant Purine Nucleoside Phosphorylase.

Notes

Purine Nucleoside Phosphorylase (PNP, E.C. 2.4.2.1.) is an enzyme involved in purine metabolism and it 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). PNP 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.

Platform Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Enzyme Mix	1 vial
Hypoxanthine Standard (10 mM)	1 x 100 μ l

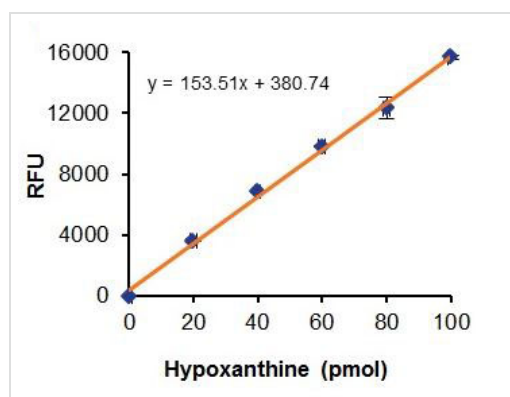
Components	100 tests
Inosine Substrate	1 x 200µl
PNP Assay Buffer (10x)	1 x 10ml
PNP Positive Control	1 vial
PNP Probe (in dry DMSO)	1 x 200µl

Involvement in disease Defects in PNP are the cause of purine nucleoside phosphorylase deficiency (PNP deficiency) [MIM:613179]. It leads to a severe T-cell immunodeficiency with neurologic disorder in children.

Sequence similarities Belongs to the PNP/MTAP phosphorylase family.

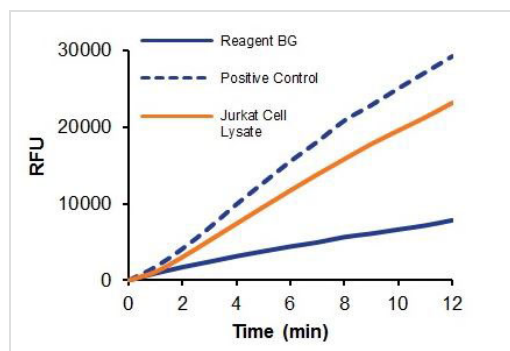
Cellular localization Cytoplasm > cytoskeleton.

Images



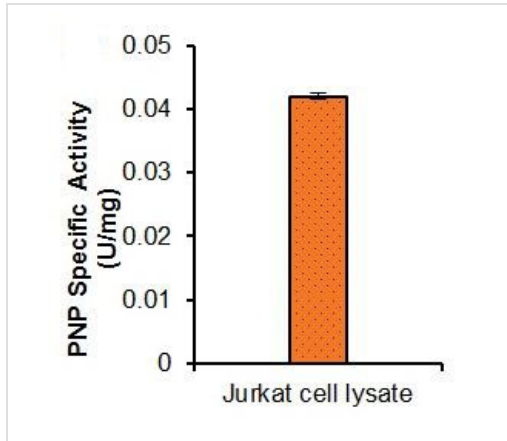
Typical Hypoxanthine Standard calibration curve.

Purine Nucleoside Phosphorylase Activity Assay Kit
(Fluorometric) (ab204706)



Purine Nucleoside Phosphorylase Activity in Jurkat Cell (T-lymphocyte) lysate (315 ng) and Positive Control; BG: Background.

Purine Nucleoside Phosphorylase Activity Assay Kit
(Fluorometric) (ab204706)



PNP specific activity in Jurkat Cell lysate.

Purine Nucleoside Phosphorylase Activity Assay Kit
(Fluorometric) (ab204706)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors