

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 µU
Species reactivity	Reacts with: Mammals, Other species
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 µU recombinant Purine Nucleoside Phosphorylase.

Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K767 Purine Nucleoside Phosphorylase Activity Assay Kit (Fluorometric). K767-100 is the same size as the 100 test size of ab204706.</p> <p>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.</p>
--------------	---

Platform	Microplate reader
-----------------	-------------------

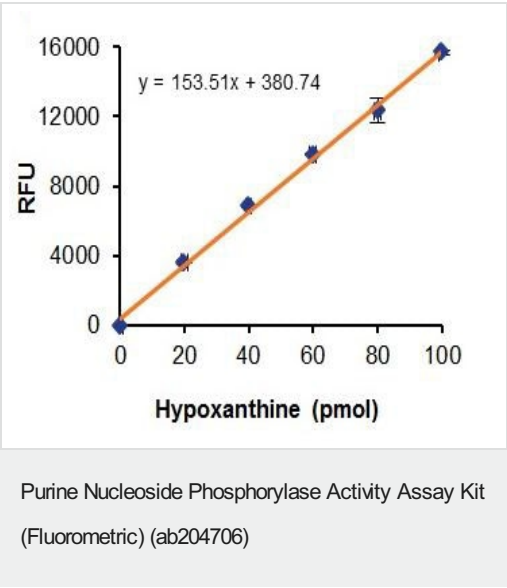
Properties

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

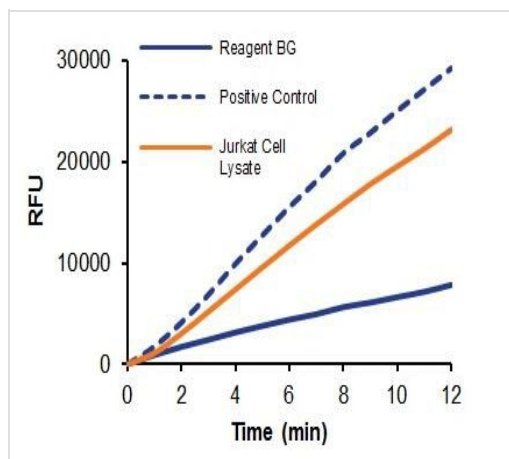
Components	100 tests
10X PNP Assay Buffer	1 x 10ml
Hypoxanthine Standard	1 x 100µl
Inosine Substrate	1 x 200µl
OxiRed Probe	1 x 200µl
PNP Enzyme Mix	1 vial
PNP Positive Control	1 vial

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

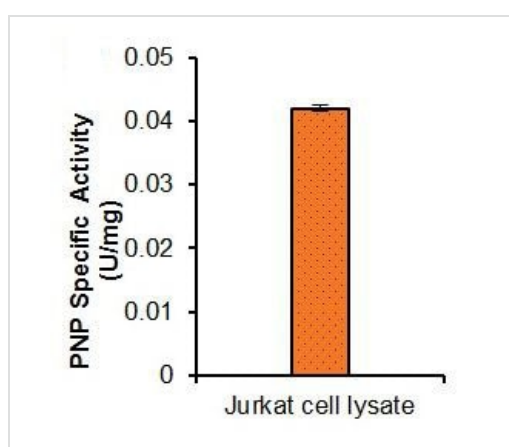


Typical Hypoxanthine Standard calibration curve.



Purine Nucleoside Phosphorylase Activity 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