

NADPH Fluorometric Kit ab287848

[3 Images](#)

Overview

Product name	NADPH Fluorometric Kit
Detection method	Fluorescent
Sample type	Tissue Extracts, Cell culture media, Purified mitochondria
Assay type	Quantitative
Sensitivity	< 40 nM
Assay duration	Multiple steps standard assay
Product overview	NADP (nicotinamide dinucleotide) is a major electron donor for anabolic reactions such as lipid and nucleic acid synthesis. The reduced form, NADPH is a cofactor for many redox enzymes. It is of increasing interest to be able to measure low level of intracellular NADPH. The New NADPH Quantitation Assay kit ab287848 (Previously known as PicoProbe NADPH Quantitation Fluorometric Assay Kit) provides a highly sensitive method to detect intracellular NADPH. NADH is not detected as the NADPH Recycling Enzyme Mix specifically recognizes NADPH in the enzyme recycling reaction. The assay is rapid, simple and does not require purifying NADPH from samples. The kit can measure less than 40 nM NADPH in a variety of samples.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called K349 PicoProbe™ NADPH Quantitation Fluorometric Assay Kit. K349-100 is the same size as the 100 test size of ab287848.
Platform	Microplate (12 x 8 well strips)

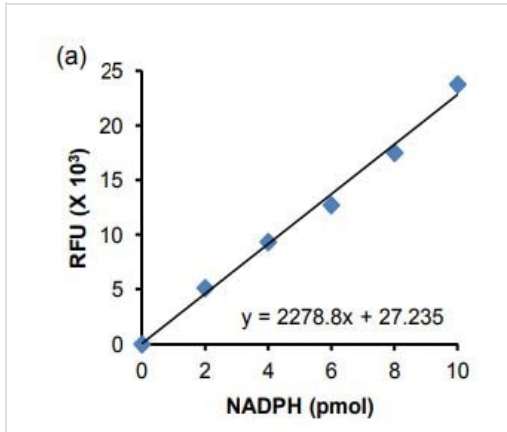
Properties

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

Components	100 tests
NADPH Cycling buffer	1 x 15ml
NADPH Cycling Enzyme Mix	1 x 0.2ml
NADPH Extraction buffer	1 x 50ml
NADPH Standard (Lyophilized)	1 vial

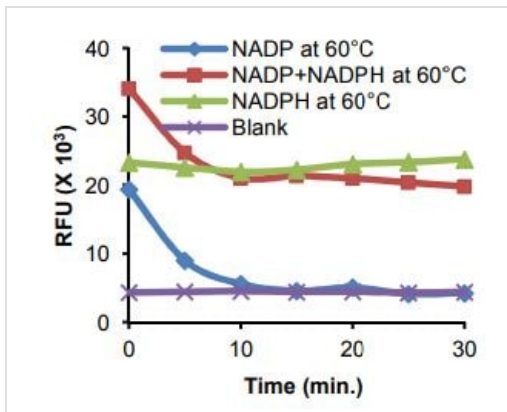
Components	100 tests
PicoProbe	1 x 0.4ml

Images



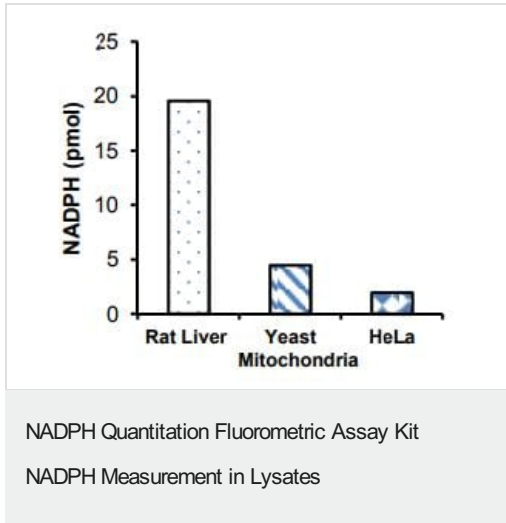
NADPH Standard Curve

NADPH Quantitation Fluorometric Assay Kit
Standard Curve



Decomposition of NADP, but not NADPH at 60°C

NADPH Quantitation Fluorometric Assay Kit
NADPH Decomposition



Measurement of NADPH in rat liver lysate (150 µg), yeast mitochondrial lysate (150 µg) and HeLa lysate (150 µg).

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