

Product datasheet

NADPH Assay Kit (Colorimetric) ab186031

1 Image

Overview

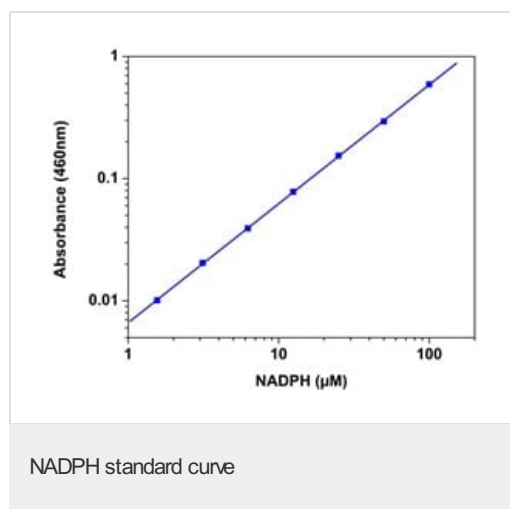
Product name	NADPH Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Tissue, Cell Lysate
Assay type	Quantitative
Sensitivity	> 3 µM
Range	3.13 µM - 100 µM
Assay time	2h 00m
Product overview	<p>Nicotinamide adenine dinucleotide (NAD⁺) and nicotinamide adenine dinucleotide phosphate (NADP⁺) are two important cofactors found in cells. NADH is the reduced form of NAD⁺. NAD forms NADP with the addition of a phosphate group to the 2' position of the adenyl nucleotide through an ester linkage. The traditional NAD/NADH and NADP/NADPH assays are based on monitoring the changes in NADH or NADPH absorption at 340 nm. The short UV wavelength of NAD/NADH and NADP/NADPH assays makes the traditional methods suffer low sensitivity and high interference.</p> <p>Abcam's Colorimetric NADPH Assay Kit (ab186031) provides a convenient method for detecting NADPH. The NADPH probe is a chromogenic sensor that has its maximum absorbance at 460 nm upon NADPH reduction. The absorbance maximum increases to ~ 635 nm if the enhancer is added to the assay system. The absorption of the NADPH probe is directly proportional to the concentration of NADPH in the solution. The Colorimetric NADPH Assay Kit provides a sensitive assay to detect as little as 3 µM NADPH in a 100 µL assay volume.</p>
Notes	<p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.</p> <p>It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
Platform	Microplate reader

Properties

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

Components	400 tests
Assay Buffer	1 x 16ml
Lysis Buffer	1 x 10ml
NADPH Probe	1 x 4ml
NADPH Standard	1 x 167µg

Images



NADPH dose response was measured with the Colorimetric NADPH Assay Kit in a 96-well white/clear bottom plate using a microplate reader. As low as 3 µM of NADPH can be detected with 30min incubation (n=3) with absorbance measurement at 460nm.

The absorbance in blank wells (with the PBS buffer only) is used as a control, and is subtracted from the values for those wells with the NADPH reactions.

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