

Product datasheet

NADPH Regeneration Kit ab228547

Overview

Product name NADPH Regeneration Kit

Product overview NADPH Regeneration Kit (ab228547) provides three ready-to-use solutions to regenerate NADPH by a simple mixing. This kit can be used for all NADPH-requiring oxidase assays (cDNA-expressed enzymes and liver fractions). About 300-500 enzyme assays can be performed using this kit. The total number of assays that can be performed depends on the experimental design.

Notes NADPH provides the reducing equivalents for biosynthetic reactions and for oxidation-reduction involved in protection against the toxicity of ROS (reactive oxygen species). NADPH is also used for anabolic pathways, such as lipid synthesis, cholesterol synthesis and fatty acid chain elongation. It is a necessary cofactor in many xenobiotic metabolism reactions. In chloroplasts, NADP is reduced to NADPH by ferredoxin-NADP reductase in the last step of the electron chain in photosynthesis reactions. The NADPH produced is used as reducing power for the biosynthetic reactions in the Calvin cycle of photosynthesis. Many oxidoreductases and all ligases use NADPH as coenzyme. NADPH is required for the measurement of oxidase activity catalyzed by P450s, FMOs, NADPH-P450 reductase, and many other oxidase enzymes.

Properties

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

Components	1000 tests
500X Glucose-6-phosphate dehydrogenase (400 units/mL)	1 x 100µl
Assay Buffer I	1 x 25ml
Assay Buffer II	1 x 25ml

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