

Pyruvate dehydrogenase (PDH) Enzyme Activity Microplate Assay Kit ab109902

★★★★☆ [4 Abreviews](#) [84 References](#) [画像数 4](#)

製品の概要

製品名	Pyruvate dehydrogenase (PDH) Enzyme Activity Microplate Assay Kit
検出方法	Colorimetric
サンプルの種類	Cell culture extracts, Tissue Lysate, Purified mitochondria
アッセイタイプ	Enzyme activity
全工程の試験時間	3h 30m
種交差性	交差種: Mouse, Rat, Cow, Human
製品の概要	<p>Pyruvate dehydrogenase (PDH) Enzyme Activity Microplate Assay Kit (ab109902) can be used to determine the activity of PDH in a human, bovine, mouse, or rat sample. The PDH enzyme is immunocaptured within the wells of the microplate and activity is determined by following the reduction of NAD⁺ to NADH, coupled to the reduction of a reporter dye to yield a colored reaction product with an increase in absorbance at 450 nm at room temperature. Included for performance of the activity assay are buffer, detergent, reagent mix, and a 96-well microplate with monoclonal antibody pre-bound to the wells of the plate, allowing for a stream-lined assay.</p> <p>This assay is optimized for use with whole tissue extract when the amount of total material available for assay is 20-100 µg or more. If using cell extract of cultured cells 1 mg of material is required due to the very low levels of enzyme and reduced levels of mitochondria in the extract.</p> <p>The PDH complex is relatively fragile and sensitive to detergent. Please follow the sample preparation steps provided in the protocol booklet. Other preparation methods may decrease PDH activity. Our Scientific Support team is happy to answer any questions about sample prep.</p> <p>Review our Metabolism Assay Guide to learn about assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also about how to assay metabolic function in live cells using your plate reader.</p>
特記事項	<p>Store 20X Buffer, Detergent, and Microplate at 4°C.</p> <p>Store 5X Stabilizer at -20°C.</p> <p>Store 20X Reagent Mix, 100X Reagent Dye and 100X Coupler at -80°C.</p>
試験プラットフォーム	Microplate reader

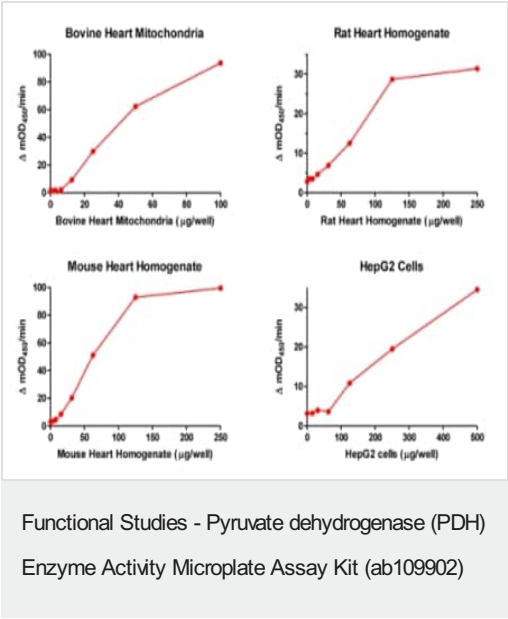
製品の特性

保存方法

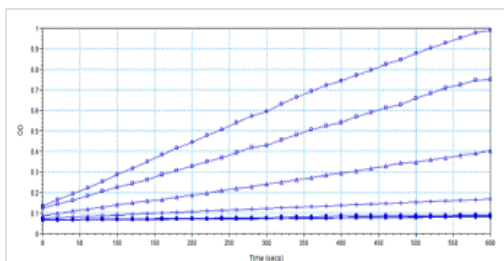
Please refer to protocols.

内容	96 tests
20X Buffer	1 x 15ml
20X Reagent Mix	2 x 600µl
5X Stabilizer	1 x 13ml
96-well Microplate (12 strips)	1 unit
Coupler	1 x 250µl
Detergent	2 x 1ml
Reagent Dye	1 x 250µl

画像

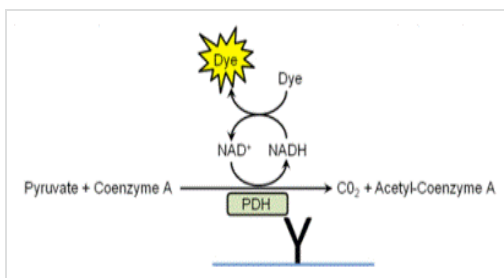


Mitochondria, tissue extracts and whole cultured cell extracts all show linear relationships between signal and sample load at limiting concentrations. The rates shown were determined as change in OD over time, and these are best represented as change in milliOD per minute.



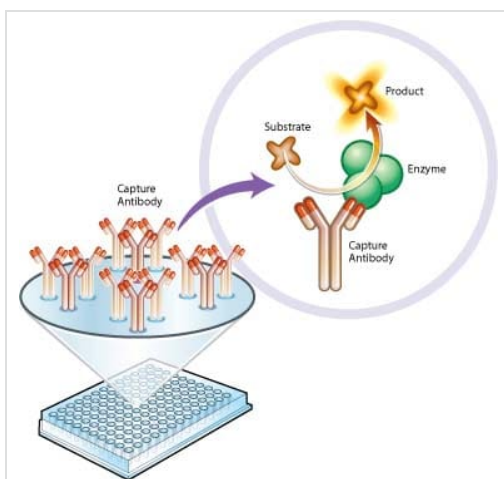
Functional Studies - Pyruvate dehydrogenase (PDH)
Enzyme Activity Microplate Assay Kit (ab109902)

Example of microplate reader recorded data from bovine heart mitochondria (100 µg/well) (top trace) and 2-fold dilutions (stepwise lower traces) using Pyruvate dehydrogenase (PDH) Enzyme Activity Microplate Assay Kit (ab109902).



Pyruvate dehydrogenase (PDH) Enzyme Activity
Microplate Assay Kit (ab109902)

Schematic diagram showing the reaction involved in Pyruvate dehydrogenase (PDH) Enzyme Activity Microplate Assay Kit (ab109902).



Functional Studies - Pyruvate dehydrogenase (PDH)
Enzyme Activity Microplate Assay Kit (ab109902)

Principle of Pyruvate dehydrogenase (PDH) Enzyme Activity
Microplate Assay Kit (ab109902)

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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