

Hydrogen Peroxide Assay Kit ab102500

82 References **画像数 3**

製品の概要

製品名	Hydrogen Peroxide Assay Kit
サンプルの種類	Serum, Plasma, Other biological fluids, Cell Lysate, Cell culture media
アッセイタイプ	Quantitative
検出感度	> 0.04 μ M
全工程の試験時間	1h 00m
製品の概要	<p>Hydrogen Peroxide Assay Kit ab102500 is a highly sensitive, simple, direct and HTS-ready colorimetric and fluorometric assay for measuring H_2O_2 in biological samples.</p> <p>In the hydrogen peroxide assay protocol, horse radish peroxidase (HRP) reacts with a probe and H_2O_2 to produce a product with color (λ_{max} = 570 nm) and red-fluorescence (Ex/Em=535/587 nm).</p> <p>The detection limit is 2 pmol per assay (or 40 nM concentration) of H_2O_2 in the sensitive fluorometric assay.</p> <p>Hydrogen peroxide assay protocol summary:</p> <ul style="list-style-type: none"> - add standards and samples to wells - add reaction mix and incubate for 10 min - analyze with microplate reader
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K265 Hydrogen Peroxide Colorimetric/Fluorometric Assay Kit. K265-200 is the same size as the 200 test size of ab102500.</p> <p>The kit can perform 200 reactions by fluorometric method or 100 reactions by colorimetric method.</p> <p>Hydrogen Peroxide (H_2O_2) is a reactive oxygen metabolic byproduct that serves as a key regulator for a number of oxidative stress-related states. Functioning through NF-kappaB and other factors, hydroperoxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis, diabetic vasculopathy, osteoporosis, neuro-degenerative diseases, Down's syndrome and immune system diseases.</p>
試験プラットフォーム	Microplate reader

製品の特性

保存方法

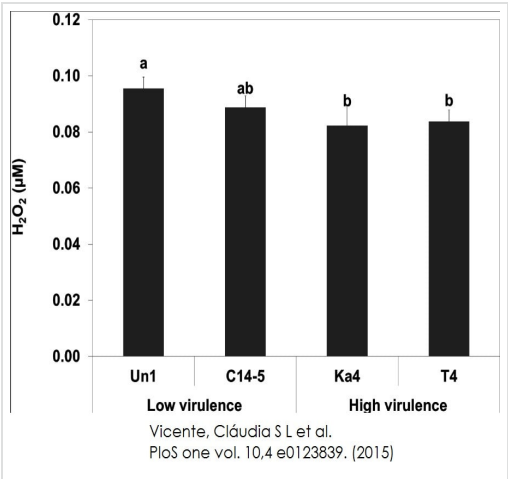
Store at -20°C. Please refer to protocols.

内容	200 tests
Assay Buffer XVIII	1 x 25ml
H2O2 Standard	1 x 0.1ml
Developer Solution V	1 unit
OxiRed Probe	1 x 200µl

関連性

Hydrogen Peroxide (H₂O₂) is a reactive oxygen metabolic byproduct that serves as a key regulator for a number of oxidative stress-related states. Functioning through NF-κB and other factors, hydroperoxide-mediated pathways have been linked to asthma, inflammatory arthritis, atherosclerosis, diabetic vasculopathy, osteoporosis, neuro-degenerative diseases, Down's syndrome and immune system diseases.

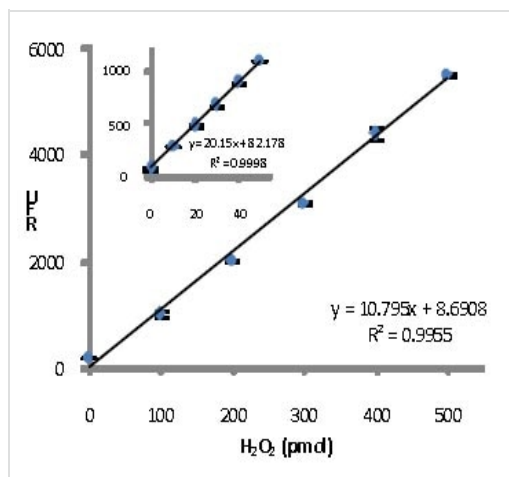
画像



H₂O₂ neutralization of of Bursaphelenchus sp. isolates (high virulence B. xylophilus Ka4 and B. xylophilus T4, and low virulence B. xylophilus C14-5 and B. mucronatus Un1) after 24 hours exposition to H₂O₂. Error bars represent standard deviation. Different letters above the columns indicate significant differences (p < 0.05) between Bursaphelenchus sp. isolates H₂O₂ neutralization, according to post-hoc Duncan's test.

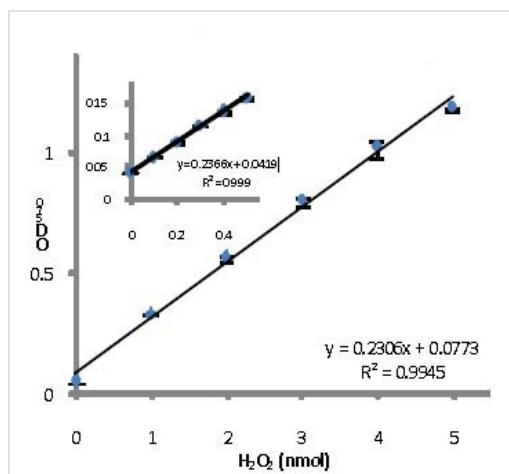
Functional Studies - Hydrogen Peroxide Assay Kit (ab102500)

Vicente, Cláudia S L et al., PloS one vol. 10,4 e0123839., Fig 2, doi:10.1371/journal.pone.0123839



H₂O₂ Fluorimetric Standard Curve

Functional Studies - Hydrogen Peroxide Assay Kit
(ab102500)



H₂O₂ Colorimetric Standard Curve

Functional Studies - Hydrogen Peroxide Assay Kit
(ab102500)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Promise 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