

D-Lactate Assay Kit (Colorimetric) ab83429

12 References 画像数 4

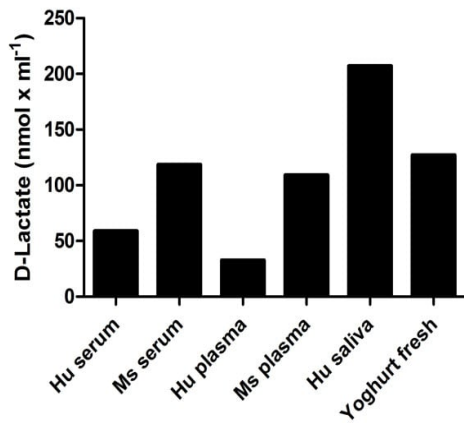
製品の概要

製品名	D-Lactate Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	> 0.01 mM
検出範囲	0.01 mM - 10 mM
全工程の試験時間	0h 40m
製品の概要	<p>D-Lactate Assay Kit (Colorimetric) Kit ab83429 provides a fast, easy way to accurately measure D-lactate in a variety of biological samples.</p> <p>In the D-lactate assay protocol, D-lactate is specifically oxidized by D-lactate dehydrogenase and generates proportional color ($\lambda_{\text{max}} = 450 \text{ nm}$). The kit detects D-Lactate in samples such as serum, plasma, cells, culture and fermentation media.</p> <p>D-lactate assay protocol summary:</p> <ul style="list-style-type: none">- add reaction mix to sample and standard wells- incubate for 30 min- analyze with a microplate reader
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K667 D-Lactate Colorimetric Assay Kit. K667-100 is the same size as the 100 test size of ab83429.</p>
試験プラットフォーム	Microplate reader

製品の特性

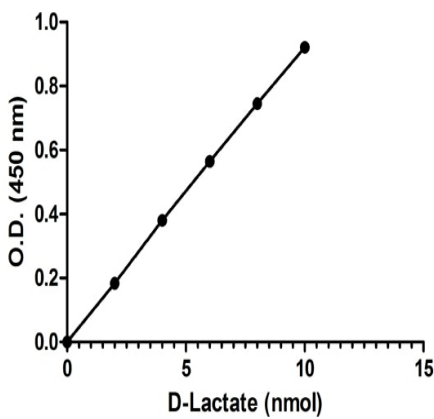
保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Assay Buffer XII	1 x 25ml
Developer Solution III	1 vial



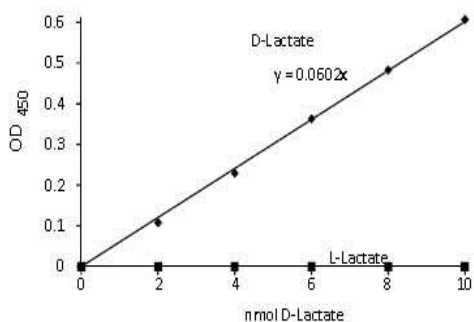
Functional Studies - D-Lactate Assay Kit
(Colorimetric) (ab83429)

D-Lactate measured in human serum, plasma and saliva, mouse serum and plasma and fresh yoghurt showing quantity (nmol) per ml of tested sample. Samples were diluted 3-27 fold.



Colorimetric standard curve: mean of duplicates (+/- SD) with background reads subtracted.

Functional Studies - D-Lactate Assay Kit
(Colorimetric) (ab83429)



D-Lactate Standard Curve. The assay is performed according to the kit instruction. The assay specifically detects D-Lactate in the presence of up to 1000X L-Lactate.

Functional Studies - D-Lactate Colorimetric Assay
Kit (ab83429)

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