

Product datasheet

D-Lactate Assay Kit (Colorimetric) ab83429

4 Images

Overview

Product name	D-Lactate Assay Kit (Colorimetric)
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Quantitative
Sensitivity	> 0.01 mM
Range	0.01 mM - 10 mM
Assay time	0h 40m
Product overview	<p>Abcam's D-Lactate Assay Kit (Colorimetric) Kit provides a fast, easy way to accurately measure D-lactate in a variety of biological samples. In this kit, D-lactate is specifically oxidized by D-lactate dehydrogenase and generates proportional color ($\lambda_{max} = 450 \text{ nm}$). The kit detects D-Lactate in samples such as serum, plasma, cells, culture and fermentation media.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Tested applications	Suitable for: Functional Studies

Properties

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

Components	Identifier	100 tests
D-Lactate Assay Buffer	WM	1 x 25ml
D-Lactate Enzyme Mix (Lyophilised)	Green	1 x 1ng
D-Lactate Standard (100 mM)	Yellow	1 x 100µl
D-Lactate Substrate Mix (Lyophilised)	Red	1 vial

Relevance D-Lactate production in mammals, mainly due to the glyoxalase pathway, is extremely low, with normal serum concentrations in the nano to micromolar range. Typically, elevated D-lactate levels which can rise to millimolar levels, are due to bacterial infection or short bowel syndrome in humans. Abnormally high concentrations of D-lactate are considered indicative of sepsis, ischemia or trauma. Due to slow metabolism and excretion, high D-lactate can cause acidosis and encephalopathy.

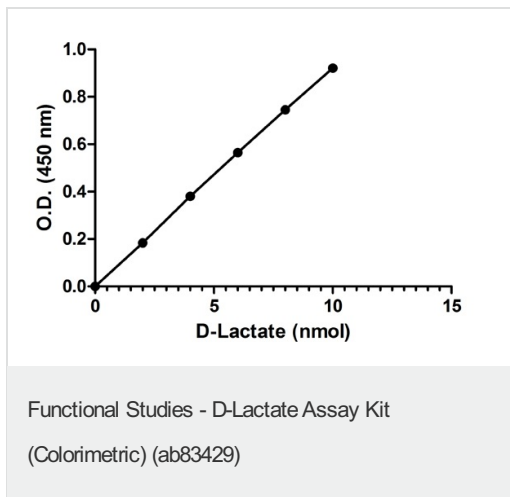
Applications

Our [Abpromise guarantee](#) covers the use of **ab83429** in the following tested applications.

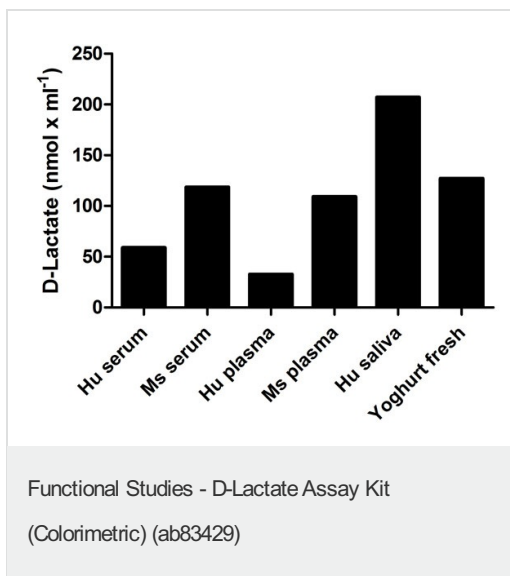
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent dilution.

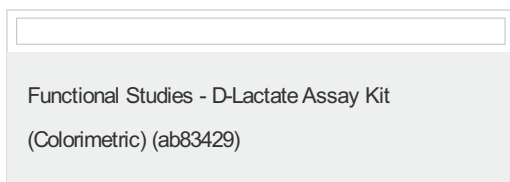
D-Lactate Assay Kit (Colorimetric) images



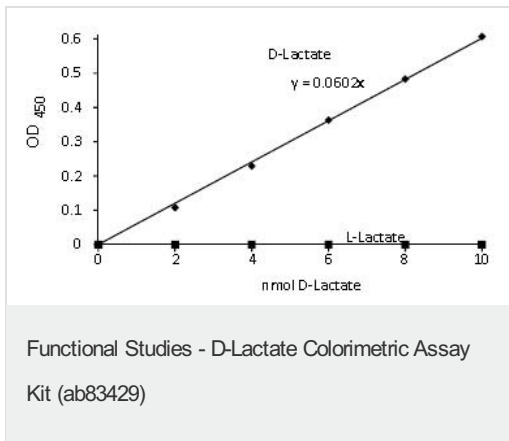
Colorimetric standard curve: mean of duplicates (+/- SD) with background reads subtracted.



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.



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



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.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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