

Fumarate Assay Kit ab102516

★★★★★ [1 Abreviews](#) [18 References](#) [5 Images](#)

Overview

Product name	Fumarate Assay Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Quantitative
Sensitivity	> 1 nmol/well
Assay time	0h 40m
Product overview	Fumarate Assay Kit ab102516 provides a convenient tool for sensitive detection of Fumarate in biological samples.

In the fumarate assay protocol, the fumarate enzyme mix recognizes fumarate as a specific substrate leading to proportional color development. The amount of fumarate can therefore be easily quantified using a colorimetric assay ($\lambda = 450 \text{ nm}$). It can detect as low as 1 nmol of fumarate per well (20 μM).

Fumarate assay protocol summary:

- add samples and standards to wells
- add reaction mix and incubate for 1 hr
- analyze with a microplate reader

Notes

This product is manufactured by BioVision, an Abcam company and was previously called K633 Fumarate Colorimetric Assay Kit. K633-100 is the same size as the 100 test size of ab102516.

Fumarate ($\text{HO}_2\text{CCH}=\text{CHCO}_2\text{H}^-$) is an intermediate in the Krebs cycle used by cells to metabolize food to form ATP. In the mammalian liver, Fumarate is also a product of the Urea cycle where its release in the cytosol leads to its conversion into malate and subsequently oxaloacetate while generating NADH in the cytosol.

Platform Microplate reader

Properties

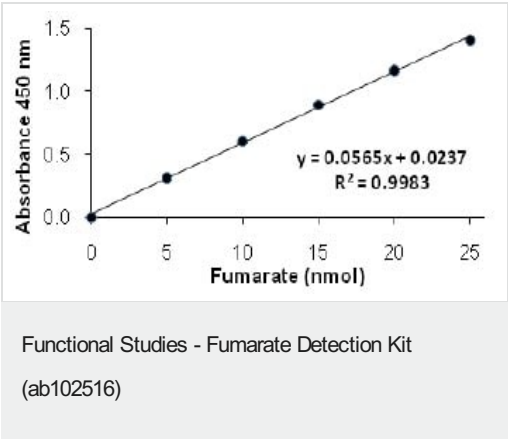
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer LXIV	1 x 25ml
Developer Solution III	1 vial
Fumarate Enzyme Mix	1 vial
Fumarate Solution	1 x 0.2ml

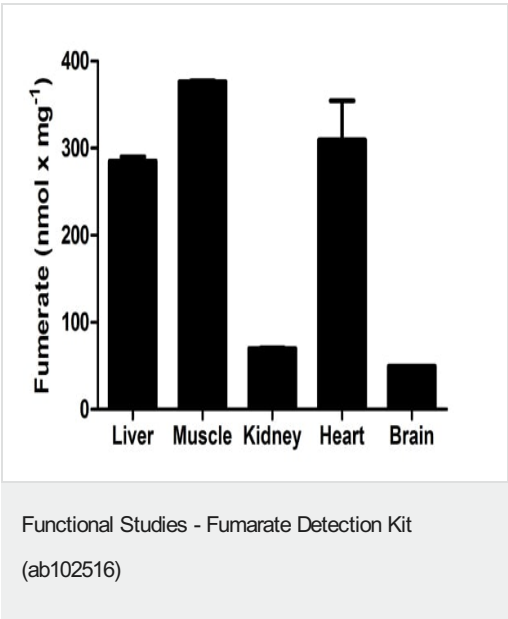
Relevance

Fumarate ($\text{HO}_2\text{CCH}=\text{CHCO}_2\text{H}^-$) is an intermediate in the Krebs's cycle used by cells to metabolize food to form ATP. In the mammalian liver, Fumarate is also a product of the Urea cycle where its release in the cytosol leads to its conversion into malate and subsequently oxaloacetate while generating NADH in the cytosol. The human skin naturally produces fumaric acid when exposed to sunlight. In fact, fumaric acid esters have been used to treat psoriasis, possibly due to an impaired production of fumaric acid in the skin. Fumaric acid has also been used in beverages, baking powders and candy.

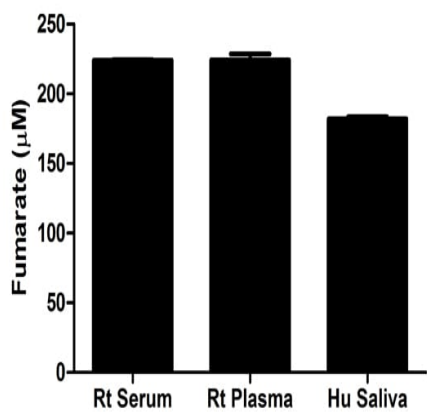
Images



Fumurate Standard Curve

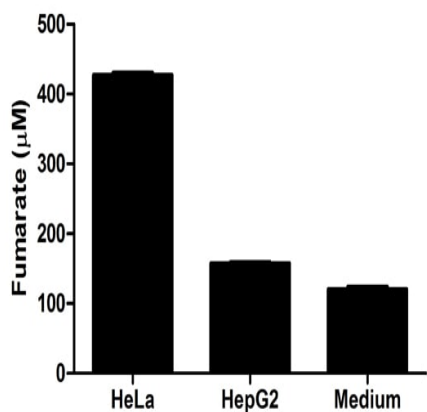


Fumarate measured in mouse tissue lysates (mg of extracted protein), background signal subtracted (duplicates +/- SD).



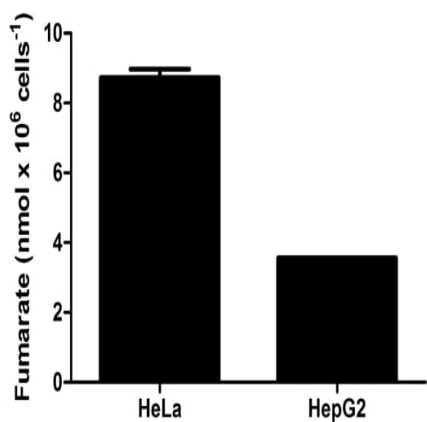
Fumarate measured in biological fluids, background signal subtracted (duplicates +/- SD).

Functional Studies - Fumarate Detection Kit
(ab102516)



Fumarate measured in cell culture medium and control medium, background signal subtracted (duplicates +/- SD).

Functional Studies - Fumarate Detection Kit
(ab102516)



Fumarate measured in cell lysates, background signal subtracted (duplicates +/- SD).

Functional Studies - Fumarate Detection Kit
(ab102516)

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