

Zinc Assay Kit ab102507

★★★★★ [1 Abreviews](#) [13 References](#) [画像数 5](#)

製品の概要

製品名	Zinc Assay Kit
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
アッセイタイプ	Quantitative
検出感度	1 µg/ml
全工程の試験時間	0h 15m
製品の概要	Zinc Assay Kit ab102507 is a convenient colorimetric assay in which Zinc binds to a ligand with development of absorbance at 560 nm.

The zinc assay can be used with biological samples such as serum, plasma, csf or urine with detection sensitivity 0.2 µg/ml (~1-3 µM).

Zinc assay protocol summary:
 - add samples and standards to wells
 - add reaction mix and incubate for 10 min
 - analyze with a microplate reader

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K387 Zinc Colorimetric Assay Kit. K387-100 is the same size as the 100 test size of ab102507.</p> <p>Zinc, a metallic chemical element, symbol Zn and atomic number 30 is chemically similar to Magnesium due to its similar size and sole oxidation state of 2+. Zinc is an essential mineral of great biological significance, because many enzymes require it as an essential cofactor.</p>
------	---

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

保存方法	Store at +4°C. Please refer to protocols.
------	---

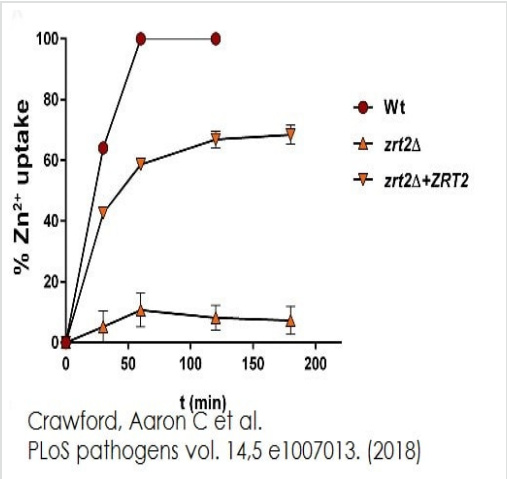
内容	100 tests
7% TCA	1 x 5ml

内容	100 tests
Zinc Reagent 1	1 x 16ml
Zinc Reagent 2	1 x 4ml
Zinc Standard	1 x 0.1ml

関連性

Zinc, a metallic chemical element, symbol Zn and atomic number 30 is chemically similar to Magnesium due to its similar size and sole oxidation state of +2. Zinc is an essential mineral of great biological significance since many enzymes require it as an essential cofactor. Examples of zinc’s biological roles include signal transduction, gene expression, regulation of apoptosis, synaptic plasticity and prostate gland function.

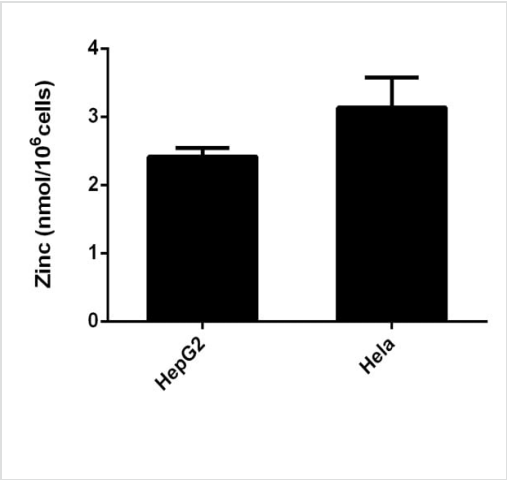
画像



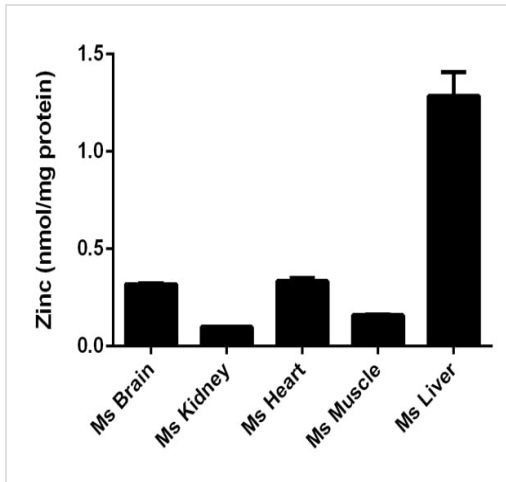
Functional Studies - Zinc Assay Kit (ab102507)

Crawford, Aaron C et al., PLoS pathogens vol. 14,5 e1007013., Fig 2, doi:10.1371/journal.ppat.1007013

Indicated strains were cultured in low zinc medium (SD0, pH ~4.7), exposed to 25 μM ZnSO4 and zinc acquisition determined at indicated time points by measuring how much zinc remained in the cell free supernatant. *C. albicans* wild type acquires all measurable zinc within 60 minute; *zrt2Δ* does not; complementation restored zinc acquisition to 68%. Experiment performed three times

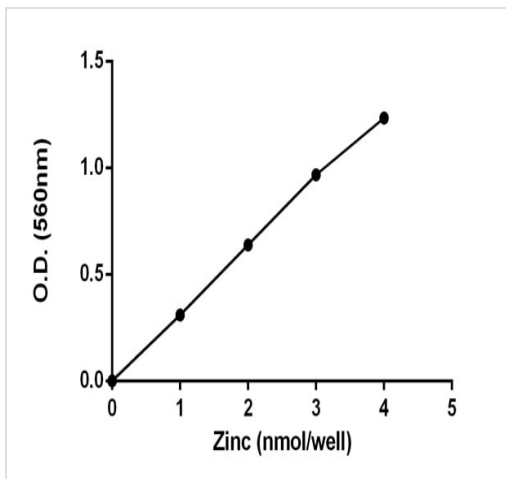


Zinc measured in cell lysates showing quantity (nmol) per 1 mln of tested cells



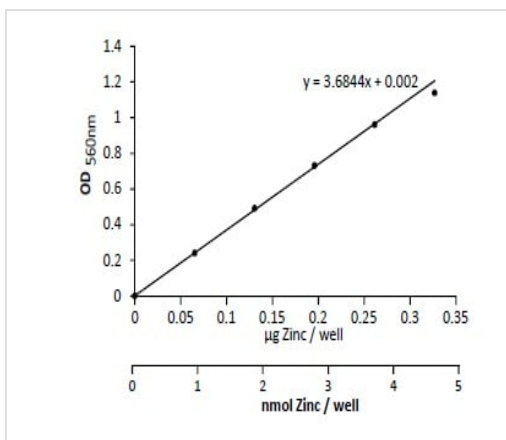
Zinc measured in mouse tissue lysates showing quantity (nmol) per mg protein of tested sample

Functional Studies - Zinc Quantification Kit
(ab102507)



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

Functional Studies - Zinc Quantification Kit
(ab102507)



Representative Standard Curve using ab102507.

Example of Standard Curve

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