

Zinc Quantification Kit (Fluorometric) ab176725

7 References 画像数 1

製品の概要

製品名	Zinc Quantification Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Urine, Serum, Plasma, Cell culture extracts, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	>= 13 ng/ml
種交差性	交差種: Mammals, Other species
製品の概要	Abcam's Zinc Quantification Kit (Fluorometric) (ab176725) provides a robust method for detecting zinc concentration in biological samples using our proprietary Zn Detector, in which Zinc binds to the probe with enhanced fluorescence at Ex/Em= 485/525 nm. The Zinc probe exhibits a large increase in fluorescence in response to Zn ²⁺ (greater than 200~300 folds). It has high Zn ²⁺ - specificity with little responses to other metals, including Ca ²⁺ , Mg ²⁺ , Mn ²⁺ , and Cu ²⁺ . The assay can be used with biological samples such as serum, plasma, and urine with detection sensitivity at 0.2 μM (13 ng/mL).
特記事項	Zinc is an essential trace mineral element that plays an important role in a number of biological processes. It is an essential factor required for many enzymes, protein structures, and control of genetic expression. Zinc status also affects basic processes of cell division, growth, differentiation, development, and aging. Clinical signs of zinc deficiency include acrodermatitis, low immunity, diarrhea, poor healing, stunting, hypogonadism, fetal growth failure, teratology and abortion. Simple, direct and automation-ready procedures for measuring zinc ion are highly desirable in research and drug discovery.
試験プラットフォーム	Microplate reader

製品の特性

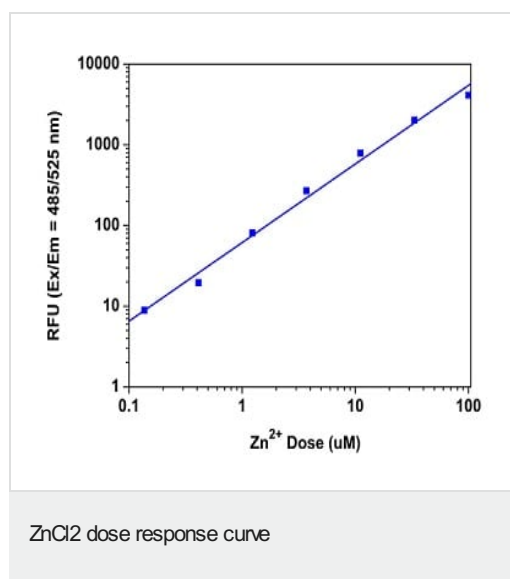
保存方法 Store at -20°C. Please refer to protocols.

内容	200 tests
Assay Buffer	1 x 15ml
Zn Detector	1 vial

内容	200 tests
ZnCl ₂ Standard	1 vial

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



ZnCl₂ dose response was measured on a 96-well black plate with Zinc Quantification Kit (Fluorometric) (ab176725). As low as ~ 0.2 μM Zn²⁺ can be detected with 5 minutes incubation time (n=3).

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