

Hemin Assay Kit ab65332

[17 References](#) [3 图像](#)

概述

产品名称	Hemin Assay试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Quantitative
检测时间	0h 30m
产品概述	<p>Hemin Assay Kit (ab65332) utilizes peroxidase activity in the presence of hemin to provide a simple, exquisitely sensitive assay which causes the conversion of a colorless probe to a strongly colored ($\lambda = 570$) compound. Trace amounts of Hemin can be quantitated in the 5-160 pg (10-250 fmol) range.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>

说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K672 Heme Colorimetric Assay Kit. K672-100 is the same size as the 100 test size of ab65332.</p> <p>Free hemin results from the breakdown of hemin-containing proteins such as hemoglobin and myoglobin. It can be detected in various body fluids such as saliva, urine and csf under various pathological states. Free hemin exists in cells at a very minute concentrations ($< 1\mu\text{M} \sim 650 \text{ ng/ml}$) exerting regulatory functions such as repression of nonspecific d-aminolevulinate synthase expression and induction of microsomal hemin oxygenase-1. Hemin can stimulate growth of oral bacteria involved in gingivitis and is an indicator of possible pathological conditions when found in the urine or feces.</p>
----	---

平台	Microplate reader
----	-------------------

性能

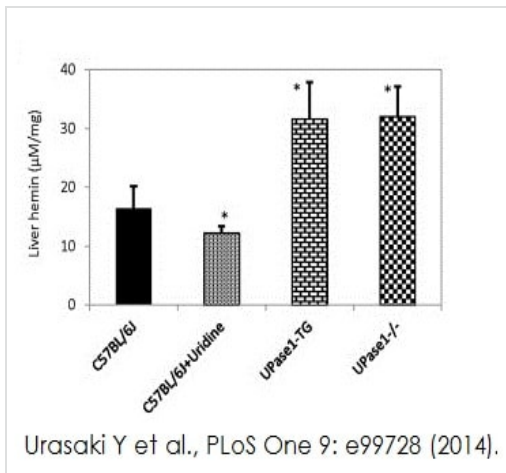
存放说明 Store at +4°C. Please refer to protocols.

组件	100 tests	2000 tests
Apoperoxidase Enzyme Mix	1 vial	20 vials
Assay Buffer V	1 x 25ml	20 x 25ml
OxiRed Probe II	1 x 0.2ml	20 x 0.2ml

组件	100 tests	2000 tests
Hemin Standard	1 vial	20 vials
Substrate V	1 x 1ml	20 x 1ml

相关性 Free hemin results from the breakdown of hemin-containing proteins such as hemoglobin and myoglobin. Hemin level has been used as an indicator of possible pathological conditions when found in the urine or faeces. It plays regulatory functions for enzyme activities.

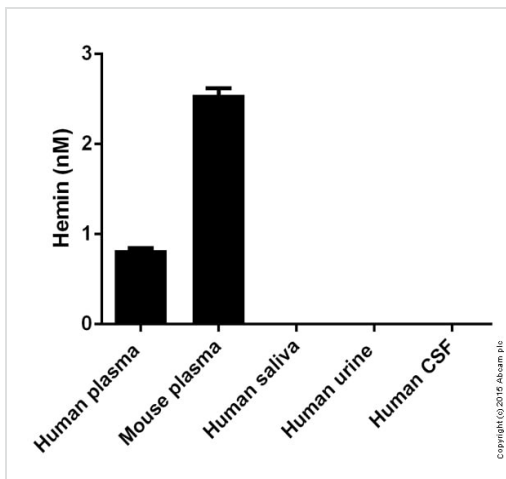
图片



Urasakai et al investigates the effect of perturbations to uridine homeostasis on heme-deficiency response pathway. Mouse liver hemin concentration was measured using ab65332.

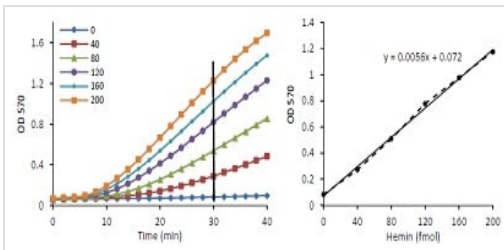
Functional studies - ab65332

Urasaki Y., PLoS One 9(6), Fig 4D. doi: 10.1371/journal.pone.0099728 Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>



Hemin measured in various biofluids showing concentration (nM)

Functional studies - ab65332 Hemin Assay Kit



A: Time course of development of color and B: Standard Curve at 30 minutes for Hemin standard.

Functional Studies - Hemin Assay Kit (ab65332)

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors