

Heme Assay Kit ab272534

画像数 1

製品の概要

製品名 Heme Assay Kit

検出方法 Colorimetric

サンプルの種類 Urine, Serum, Plasma, Whole Blood

アッセイタイプ Quantitative

検出範囲 0.6 μ M - 125 μ M

製品の概要

Heme Assay Kit (ab272534) is a simple, direct and automation-ready procedure for measuring heme concentration. This Assay Kit is based on an improved aqueous alkaline solution method, in which the heme is converted into a uniform colored form. The intensity of color, measured at 400 nm, is directly proportional to the heme concentration in the sample. The optimized formulation substantially reduces interference by substances in the raw samples and exhibits high sensitivity.

Sensitive and accurate: Linear detection range 0.6 – 125 μ M heme in 96-well plate assay.

Simple and high-throughput: The “mix-and-read” procedure involves addition of a single working reagent and reading the optical density. Can be readily automated as a high-throughput assay in 96-well plates for thousands of samples per day.

Safety: Reagents are non-toxic.

Versatility: Assays can be executed in 96-well plate or cuvette.

特記事項

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

アプリケーション

適用あり: Functional Studies

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

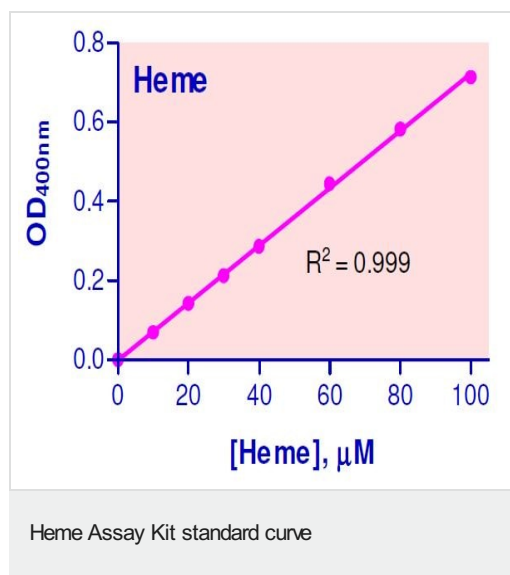
内容	250 tests
Calibrator	1 x 10ml
Reagent	1 x 50ml

アプリケーション

The Abpromise guarantee **Abpromise保証は、次のテスト済みアプリケーションにおけるab272534の使用に適用されず**
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

画像



Standard curve with freshly prepared Heme - Dilution series for freshly prepared Heme for visualization of the assay linearity - please note standard curve generation is not required for this assay.

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