

NQO1 Activity Assay Kit ab184867

13 References 6 图像

概述

| | | | | |
|-------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|---|------|--------|
| 产品名称 | NQO1 Activity Assay试剂盒 | | | |
| 检测方法 | Colorimetric | | | |
| 精确度 | 批次内 | | | |
| | 样品 | n | Mean | SD |
| | Overall | 3 | | 2.183% |
| | 批次间 | | | |
| | 样品 | n | Mean | SD |
| | Overall | 9 | | 6.035% |
| 样品类型 | Cell culture extracts, Adherent cells, Suspension cells, Tissue Extracts, Cell Lysate, Tissue Homogenate, Tissue Lysate | | | |
| 检测类型 | Enzyme activity | | | |
| 检测时间 | 0h 30m | | | |
| 种属反应性 | 与反应: Mouse, Rat, Cow, Human | | | |
| 产品概述 | <p>NQO1 Activity Assay kit (ab184867) is designed for the sensitive and accurate measurement of NQO1 (NAD(P)H dehydrogenase [quinone] 1) activity.</p> <p>The enzyme activity is determined by following the reduction of Menadione with cofactor NADH and the simultaneous reduction of WST1 which leads to increased absorbance at 440 nm. Dicoumarol is used as an inhibitor of NQO1 as part of the assay.</p> <p>NQO1 assay protocol summary:</p> <ul style="list-style-type: none">- add samples to wells- add reaction buffer + inhibitor and reaction buffer to separate sample wells- analyze with a microplate reader for 5 min | | | |
| 说明 | <p>NQO1 serves as a quinone reductase in connection with conjugation reactions of hydroquinones involved in detoxification pathways, NQO1 protects against quinone-induced damage by competing with potentially toxic one-electron pathways. It also functions in biosynthetic processes such as the vitamin K-dependent gamma-carboxylation of glutamate residues in prothrombin</p> | | | |

| | |
|----|-------------------|
| | synthesis. |
| 平台 | Microplate reader |
| 性能 | |

存放说明

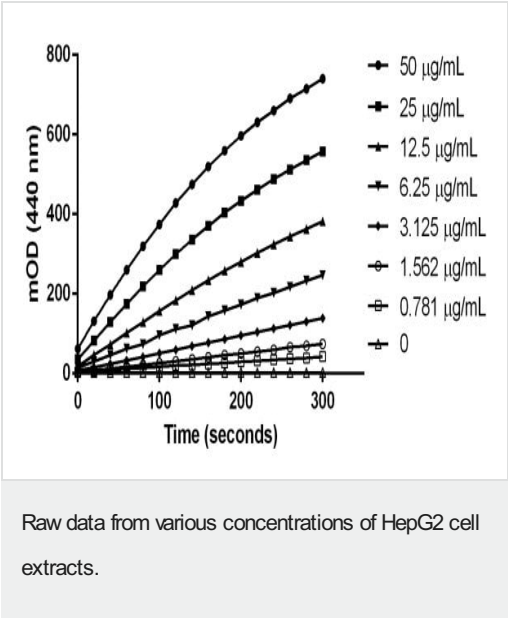
Store at -20°C. Please refer to protocols.

| 组件 | 1 x 96 tests |
|-------------------------------|--------------|
| 1000X Cofactor (lyophilized) | 1 vial |
| 1000X Inhibitor (lyophilized) | 1 vial |
| 100X Dye (lyophilized) | 1 vial |
| 20X Basic Buffer | 1 x 3ml |
| 2X Extraction Buffer | 1 x 15ml |
| 5000X Menadione | 1 x 100µl |
| 500X NADH (lyophilized) | 1 vial |
| 96-well Microplate | 2 units |

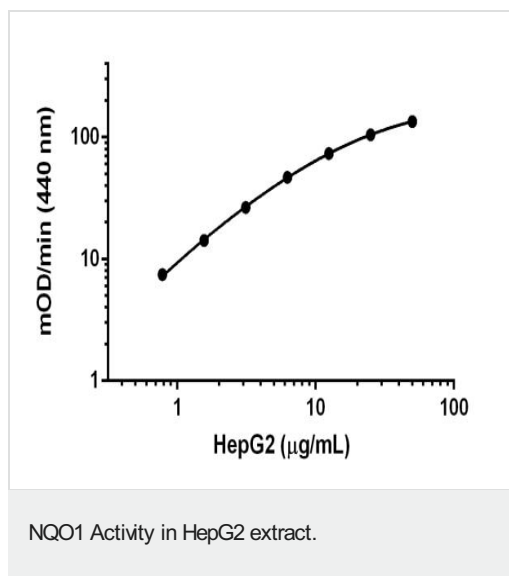
细胞定位

Cytoplasmic

图片



An example is shown below where the rate/slope is calculated between these time points.



This raw data is expressed as rate (mOD/min) per microgram of cell extract added per well as shown above. The molar extinction coefficient factor for the WST1 dye is 25.9/mM/well.

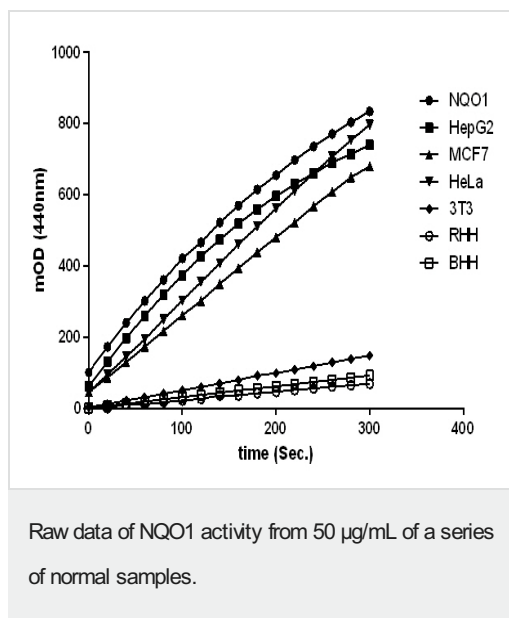


Figure 3a

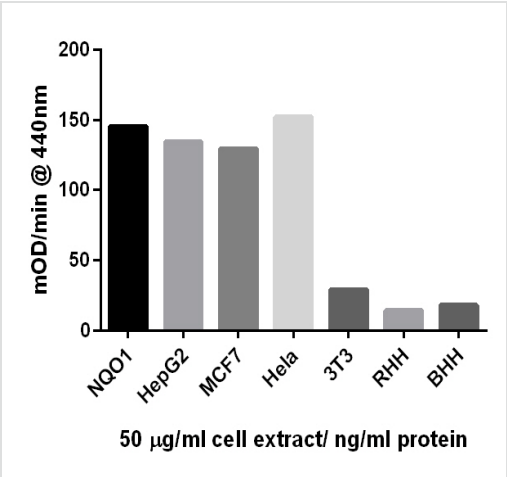


Figure 3b

Comparison of NQO1 activity in different samples at 50 ug/mL

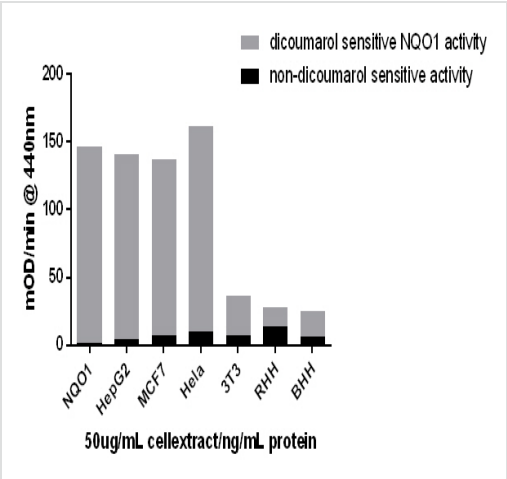


Figure 3c

Dicoumarol sensitive NQO1 activity/total activity in a series of normal cell lysates at 50 ug/mL.

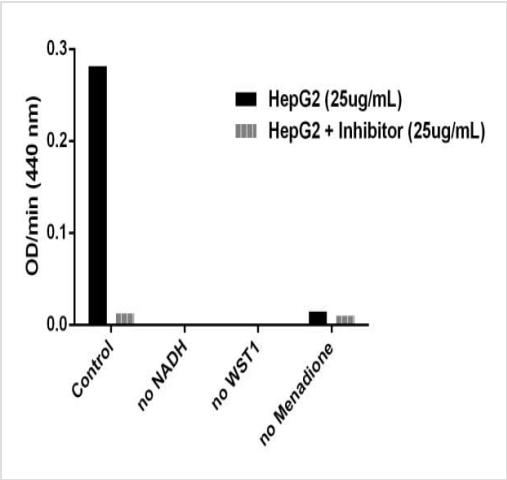


Figure 4

NQO1 activity assay components requirement test.

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