

# Glutathione Peroxidase Assay Kit (Colorimetric) ab102530

★★★★★ [1 Abreviews](#) [111 References](#) [画像数 6](#)

### 製品の概要

製品名	Glutathione Peroxidase Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Platelets, Other biological fluids, Tissue Extracts
アッセイタイプ	Enzyme activity
検出感度	0.5 mU/ml
全工程の試験時間	0h 40m
種交差性	交差種: Mammals, Other species
製品の概要	Glutathione Peroxidase Assay Kit (Colorimetric) ab102530 can be used to quantitate the activity of all of the glutathione dependent peroxidases in plasma, erythrocyte lysates, tissue homogenates, and cell lysates.

In the glutathione peroxidase assay protocol, glutathione peroxidase (GPx) oxidizes GSH to produce GSSG as part of the reaction in which it reduces cumene hydroperoxide. Glutathione reductase (GR) then reduces the GSSG to produce GSH, and in the same reaction consumes NADPH. The decrease of NADPH (measured at OD=340 nm) is proportional to GPx activity.

The assay has a detection sensitivity of ~ 0.5 mU/ml of GPx in samples.

Glutathione peroxidase assay protocol summary:

- add samples and standards to wells
- add reaction mix and incubate for 15 min at room temp to deplete all GSSG in sample
- add cumene hydroperoxide
- analyze with microplate reader immediately, and after at least 5 min

**特記事項** This product is manufactured by BioVision, an Abcam company and was previously called K762 Glutathione Peroxidase Activity Colorimetric Assay Kit. K762-100 is the same size as the 100 test size of ab102530.

The Glutathione Peroxidase (GPx, EC 1.11.1.9) family of enzymes plays an important role in the protection of organisms from oxidative damage. GPx converts reduced glutathione (GSH) to oxidized glutathione (GSSG) while reducing lipid hydroperoxides to their corresponding alcohols or free hydrogen peroxide to water.

**試験プラットフォーム** Microplate reader

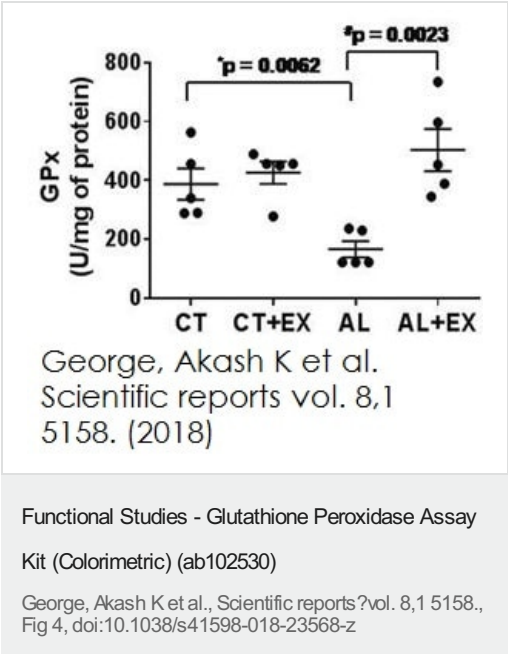
製品の特性

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

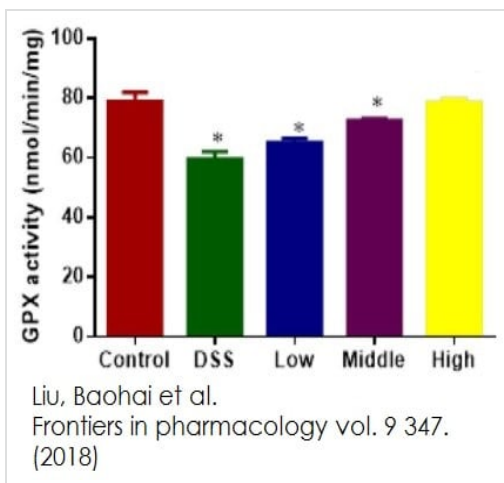
内容	100 tests	2000 tests
Cumene Hydroperoxide	1 vial	20 vials
Glutathione	1 vial	20 vials
Active Glutathione Reductase	1 vial	20 vials
Glutathione Peroxidase Assay Buffer	1 x 50ml	20 x 50ml
Active Glutathione Peroxidase	1 vial	20 vials
NADPH II	1 vial	20 vials

関連性Glutathione Peroxidase (GPx, EC 1.11.1.9) is an enzyme family with peroxidase activity, and plays important role in protecting of organisms from oxidative damage. It converts reduced glutathione (GSH) to oxidized glutathione (GSSG), to reduce lipid hydroperoxides to their corresponding alcohols, or reduce free hydrogen peroxide to water. Several isozymes have been found in different cellular locations and with different substrate specificity. Low levels of GPx have been correlated with free radical related disorders.

画像



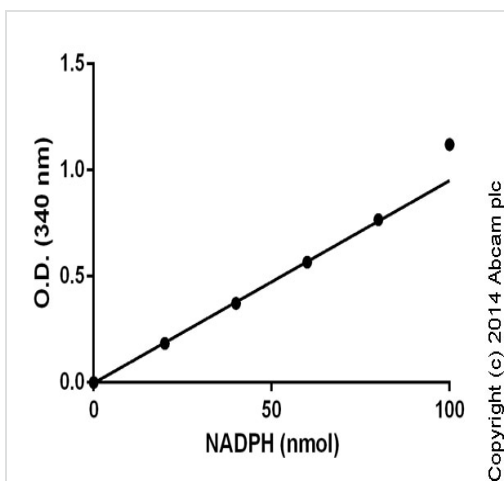
Scatter dot plots representing the levels of glutathione peroxidase (GPx) in brain tissue in different mice groups.



The effects of PCSE on the activity of GPX in a mouse UC model.

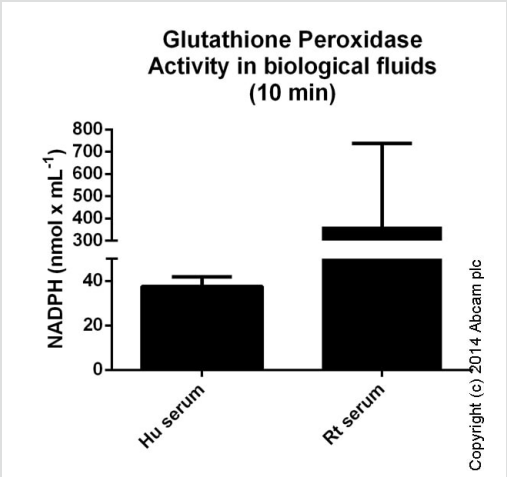
Functional Studies - Glutathione Peroxidase Assay  
Kit (Colorimetric) (ab102530)

Liu, Baohai et al., Frontiers in pharmacology?vol. 9  
347., Fig 8, doi:10.3389/fphar.2018.00347



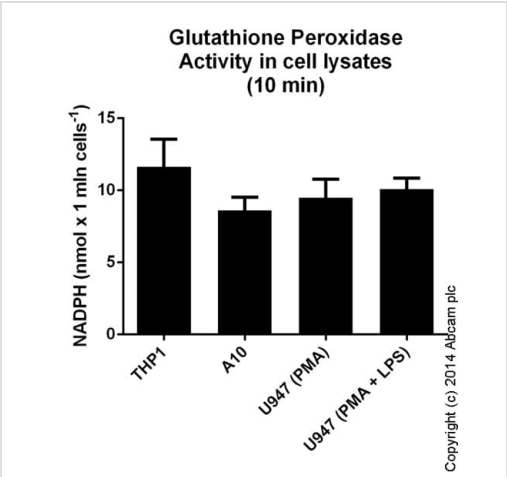
Standard curve: mean of duplicates (+/- SD) with background reads  
subtracted

Functional Studies - Glutathione Peroxidase Assay  
Kit (ab102530)



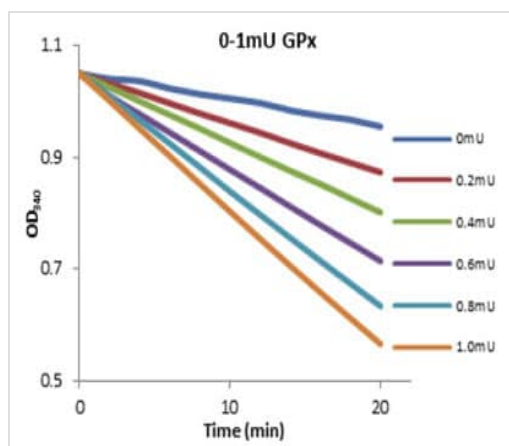
Functional Studies - Glutathione Peroxidase Assay  
Kit (ab102530)

Glutathione Peroxidase Activity measured in biological fluids showing quantity (nmol) per 1 mL after 10 min of incubation. Samples diluted 1-27 fold.



Functional Studies - Glutathione Peroxidase Assay  
Kit (ab102530)

Glutathione Peroxidase Activity measured in cell lysates showing quantity (nmol) per 1 mln cells after 10 min of incubation. Undiluted sample contained 2.5e6 cells per well. Samples were diluted 1-9 fold.



Tests example obtained using ab102530

Functional Studies - Glutathione Peroxidase Assay  
Kit (ab102530)

**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