

GSH/GSSG Ratio Detection Assay Kit II (Fluorometric - Green) ab205811

[32 References](#) [3 图像](#)

概述

产品名称	GSH/GSSG Ratio Detection Assay试剂盒II (Fluorometric - Green)
检测方法	Fluorescent
样品类型	Urine, Plasma, Tissue Extracts, Cell Lysate
检测类型	Quantitative
灵敏度	10 nM
检测时间	0h 30m
产品概述	GSH/GSSG Ratio Detection Assay Kit II (Fluorometric - Green) (ab205811) provides an ultrasensitive assay to quantitate glutathione in the sample.

The GSH/GSSG assay protocol uses a proprietary non-fluorescent water-soluble dye that becomes strongly fluorescent upon reacting with GSH. With a one-step fluorimetric method, the kit can detect as little as 1 picomole of GSH or GSSG in a 100 μ L assay volume.

The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and readily adapted to automation without a separation step. Its signal can be easily read by a fluorescence microplate reader at Ex/Em = 490/520 nm.

GSH/GSSG assay protocol summary:

- add samples (deproteinized) and standards to wells
 - for GSH assay add Thiol Green in assay buffer, or for total glutathione (GSH + GSSG) assay also add GSSG probe
 - incubate for 10 - 60 min at room temp
 - analyze with microplate reader
- GSSG levels can be calculated by subtracting GSH from total glutathione levels.

说明 NOTE: For measuring GSH Standard only, there is enough reagent provided to perform 200 tests.

This product contains water-soluble Thiol green indicator. We recommend this kit as a replacement for **GSH/GSSG Ratio Detection Assay Kit (Fluorometric - Green) (ab138881)**, which uses a DMSO-soluble probe.

Background information on GSH/GSSG

Glutathione (GSH) is a tripeptide that contains L-cysteine, L-glutamic acid and glycine. It is the smallest intracellular protein thiol molecule in the cells, which prevents cell damage caused by reactive oxygen species such as free radicals and peroxides. Glutathione exists in reduced (GSH) and oxidized (GSSG) states.

Reduced glutathione (GSH) is a major tissue antioxidant that provides reducing equivalents for the glutathione peroxidase (GPx) catalyzed reduction of lipid hydroperoxides to their corresponding alcohols and hydrogen peroxide to water. In the GPx catalyzed reaction, the formation of a disulfide bond between two GSH molecules generates oxidized glutathione (GSSG).

Glutathione reductase (GR) recycles GSSG to GSH with the simultaneous oxidation of β -nicotinamide adenine dinucleotide phosphate (β -NADPH₂).

In healthy cells, >90% of the total glutathione pool is in the reduced form (GSH). When cells are exposed to increased levels of oxidative stress, GSSG accumulates and the ratio of GSSG to GSH increases. An increased ratio of GSSG-to-GSH is an indication of oxidative stress.

平台

Microplate reader

性能

存放说明

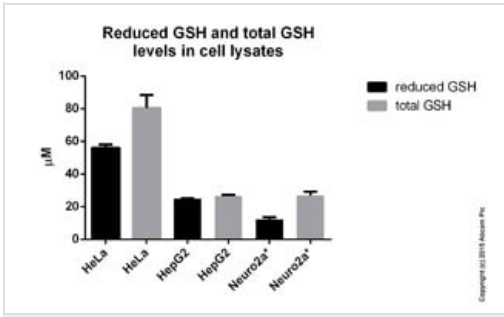
Store at -20°C. Please refer to protocols.

组件	100 tests
Assay Buffer	1 x 25ml
GSH Standard	1 x 62 μ g
GSSG Probe	1 vial
GSSG Standard	1 x 124 μ g
Thiol Green Indicator WS	1 vial

相关性

Glutathione is a small peptide composed of three amino acids: cysteine, glutamic acid, and glycine and is present in tissues in concentrations as high as one millimolar. Glutathione is the principal intracellular low-molecular-weight thiol that plays a critical role in the cellular defense against oxidative and nitrosative stress in mammalian cells. Diminished glutathione levels have been observed in the early stages of apoptosis.

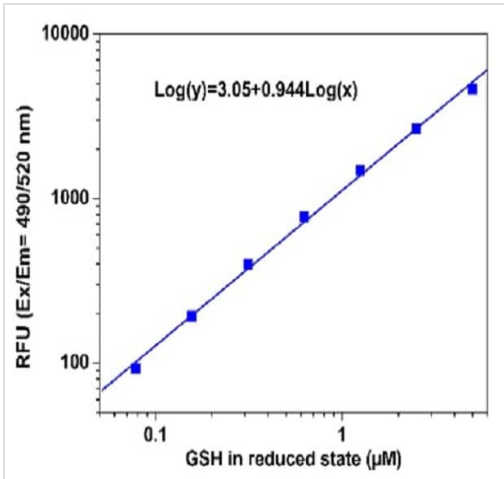
图片



GSH/GSSG Ratio Detection Assay Kit

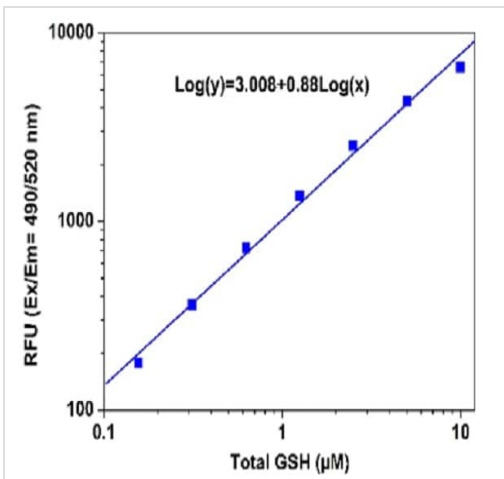
Reduced GSH and total GSH levels in cell lysates. Cells lysed to the concentration of 1×10^7 cells per mL and tested diluted 6-54 fold.

* Cells serum starved for 24 hours.



Sample Calibration Curve - Reduced GSH

Reduced GSH dose responses were measured in a black 96-well plate with ab205811 using a fluorescence microplate reader. 50 μL of GSH standards (0.01 to 5 μM) or blank control was added into each well, and then 50 μL of GSH Assay Mixture was added. The fluorescence intensity was measured at Ex/Em = 490/520 nm after 30 minutes incubation.



Sample Calibration Curve - Total GSH

Total GSH dose responses were measured with ab205811 in a black 96-well plate using a fluorescence microplate reader. 50 μL of GSSG standards (0.01 to 5 μM), GSH-containing samples or blank control were added into each well, and then 50 μL of Total GSH Reaction Mixture was added. Fluorescence intensity was measured at Ex/Em = 490/520 nm after 30 minutes incubation.

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