

GST Activity Assay Kit (Fluorometric) ab65325

[2 References](#) [画像数 1](#)

製品の概要

製品名	GST Activity Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Saliva, Urine, Serum, Plasma, Other biological fluids, Cell Lysate, Purified protein, Tissue Lysate
アッセイタイプ	Enzyme activity
検出感度	< 0.5 mU/well
全工程の試験時間	2h 00m

製品の概要

Abcam's GST Activity Assay Kit (Fluorometric) provides a simple, fluorescence-based *in vitro* assay for detecting the GST activity using fluorescence plate reader. The assay utilizes monochlorobimane (MCB), a dye that reacts with glutathione. The free form of MCB is almost nonfluorescent, whereas the dye fluoresces in blue (Ex/Em = 380/461 nm) after reaction with glutathione catalysed by GST. The change in fluorescence overtime allows for an easy measurement of the sample GST level.

The kit can detect GST activity in crude cell lysate or purified protein fraction, and also quantitate GST-tagged fusion protein.

Visit our [FAQs page](#) for tips and troubleshooting.

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K260 GST Fluorometric Activity Assay Kit. K260-100 is the same size as the 100 test size of ab65325.

Glutathione S-transferase (GST) is a family of enzymes that play an important role in detoxification of xenobiotics. GST catalyzes the formation of the thiol group of glutathione to electrophilic xenobiotics. It utilizes glutathione to scavenge potentially toxic compounds including those produced as a result of oxidative stress and is part of the defense mechanism against the mutagenic, carcinogenic and toxic effects of such compounds.

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.

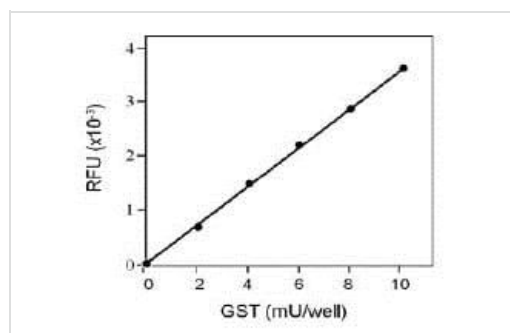
内容	100 tests
GSH Substrate	2 vials

内容	100 tests
GST Assay Buffer	1 x 25ml
GST Enzyme Standard	1 x 10 μ l
MCB Probe	1 x 200 μ l

関連性 Glutathione S-transferase (GST) is a family of enzymes that play an important role in detoxification of xenobiotics. GST catalyzes the formation of the thiol group of glutathione to electrophilic xenobiotics. It utilizes glutathione to scavenge potentially toxic compounds including those produced as a result of oxidative stress and is part of the defense mechanism against the mutagenic, carcinogenic and toxic effects of such compounds. Based on their biochemical, immunological, and structural properties, the soluble human GSTs are categorized into 4 main classes: alpha, mu, pi, and theta.

細胞内局在 Cytoplasmic

画像



Standard Calibration Curve of GST Measured by Fluorometry. Various amounts of standard GST were incubated with GSH and MCB according to the kit instructions. Fluorescence was measured at ex/em 380/460 nm.

Functional Studies - GST Activity Assay Kit
(Fluorometric) (ab65325)

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