

## Product datasheet

# GST Assay Kit (Colorimetric) ab65326

[11 References](#) [2 Images](#)

### Overview

---

<b>Product name</b>	GST Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate
<b>Assay type</b>	Enzyme activity
<b>Sensitivity</b>	< 4 mU/well
<b>Assay time</b>	1h 30m
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	GST Assay Kit (Colorimetric) (ab65326) is used to detect GST activity in crude cell lysates and purified protein fractions, and also to quantitate GST-tagged fusion proteins.

The GST assay protocol is based on the GST-catalyzed reaction between GSH and the GST substrate CDNB.

CDNB has the broadest range of isozyme detectability (e.g., alpha-, mu-, pi-, and other GST isoforms). The GST-catalyzed formation of GS-DNB produces a dinitrophenyl thioether which can be detected by spectrophotometer at 340 nm.

One unit of GST activity is defined as the amount of enzyme producing 1  $\mu$ mol of GS-DNB conjugate/min under the conditions of the assay. Detection limit of assay: active GST < 1 mU.

GST assay protocol summary:

- add samples and positive control to wells
- add glutathione to wells
- add reaction mix
- analyze with a microplate reader for at least 10 min

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K263 GST Colorimetric Activity Assay Kit. K263-100 is the same size as the 100 test size of ab65326.

Glutathione S-transferase (GST, EC 2.5.1.13) 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.

**Platform**

Microplate reader

**Properties****Storage instructions**

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

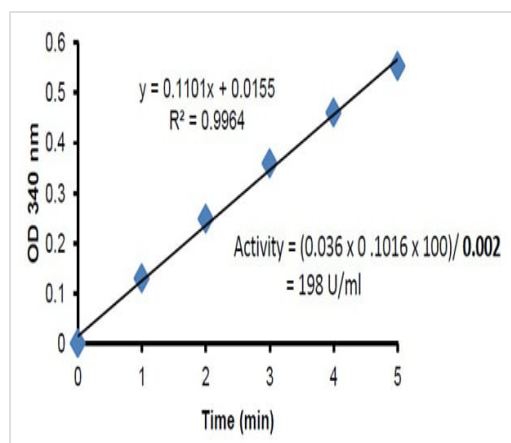
Components	100 tests
Reducing Agent I	2 x 17mg
GST Assay Buffer	1 x 25ml
GST Positive Control	1 x 10µl
GST Substrate	1 x 0.1ml

**Relevance**

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.

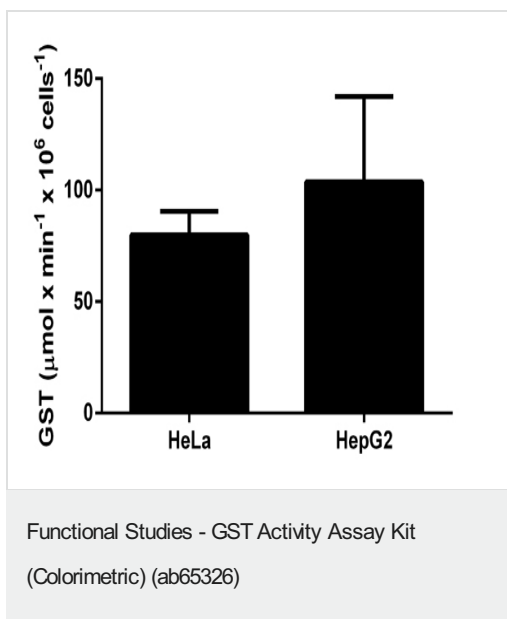
**Cellular localization**

Cytoplasmic

**Images**

GST Kinetic Assay performed according to the protocol.

Functional Studies - GST Activity Assay Kit  
(Colorimetric) (ab65326)



GST activity calculated in cell lysates, with background signal subtracted (duplicates +/- SD), using equation in protocol.

**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.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors