

### GST His-tag ELISA Kit ab302758

Recombinant SimpleStep ELISA<sup>®</sup>

[4 Images](#)

#### Overview

**Product name** GST His-tag ELISA Kit

**Detection method** Colorimetric

#### Precision

Intra-assay

Sample	n	Mean	SD	CV%
Supernatant	8			7.8%

Inter-assay

Sample	n	Mean	SD	CV%
Supernatant	3			9.7%

**Sample type** Cell culture supernatant, Cell culture media

**Assay type** Sandwich (quantitative)

**Sensitivity** 3.476 pg/ml

**Range** 35.156 pg/ml - 2250 pg/ml

#### Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	108	92% - 120%
Cell culture media	91	81% - 110%

**Assay time** 1h 30m

**Assay duration** One step assay

#### Product overview

GST His-tag SimpleStep ELISA<sup>®</sup> kit is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of GST His-tag protein. Quantitate GST His-tag with 3.476 pg/ml sensitivity.

SimpleStep ELISA<sup>®</sup> technology employs capture antibodies conjugated to an affinity tag that is

recognized by the monoclonal antibody used to coat our SimpleStep ELISA<sup>®</sup> plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA<sup>®</sup> protocol summary in the image section for further details. Our SimpleStep ELISA<sup>®</sup> technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA<sup>®</sup> microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA<sup>®</sup> kits.

**Notes**

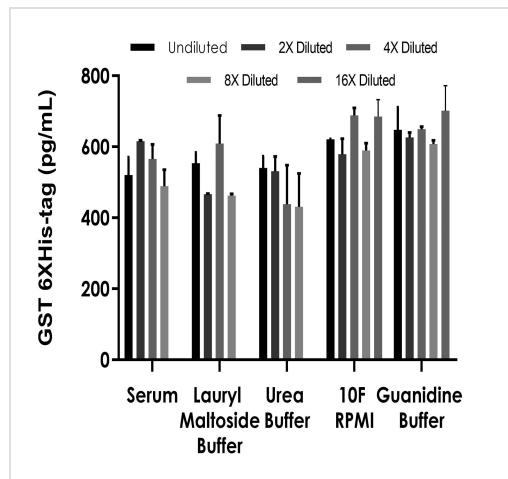
Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

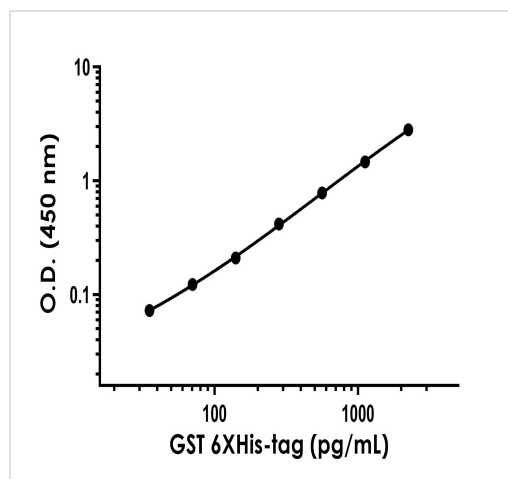
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	96 tests
10X GST His-tag Capture Antibody	1 x 600µl
10X GST His-tag Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BR	1 x 6ml
GST His-tag Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml



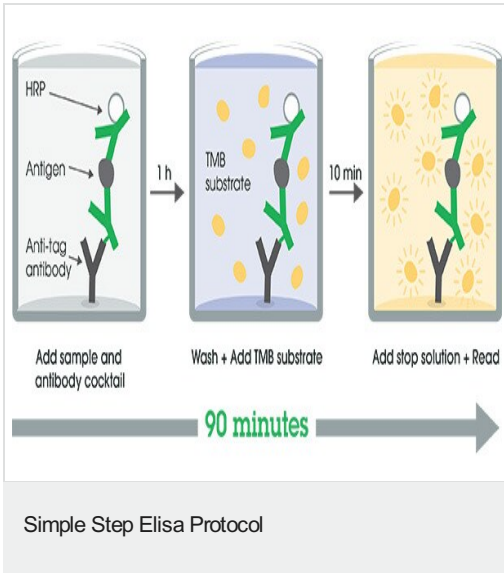
Interpolated concentrations of spiked species GST His-tag in serum, lauryl maltoside buffer, urea buffer, 10F RPMI, and Guanidine buffer.

Recombinant GST His-tag was spiked into the following samples and diluted in a 2-fold dilution series in 1X Cell Extraction Buffer PTR. Undiluted samples are as follows: serum 10%, lauryl maltoside buffer 50%, urea buffer 10%, 10F RPMI 50%, and guanidine buffer 0.1%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).



Example of species GST His-tag standard curve in 1X Cell Extraction Buffer PTR.

Example of species GST His-tag standard curve. Background-subtracted data values (mean +/- SD) are graphed.



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

### Why choose a recombinant antibody?

<p><b>Research with confidence</b> Consistent and reproducible results</p>	<p><b>Long-term and scalable supply</b> Recombinant technology</p>
<p><b>Success from the first experiment</b> Confirmed specificity</p>	<p><b>Ethical standards compliant</b> Animal-free production</p>

**Benefits of Recombinant Antibodies**

**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