

Glucagon-Like Peptide 2 ELISA Kit ab222863

★★★★★ [1 Abreviews](#) [1 Image](#)

Overview

Product name Glucagon-Like Peptide 2 ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
Overall	60			4.4%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	60			10.5%

Sample type Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.14 ng/ml

Range 0.563 ng/ml - 4.5 ng/ml

Recovery 97 %

Assay time 4h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse, Rat, Horse, Human, Pig, Monkey  
**Does not react with:** Rabbit, Cow

Product overview The Glucagon-Like Peptide 2 ELISA (Enzyme-Linked Immunosorbent Assay) Kit (ab222863) kit is designed for detection of GLP-2 in plasma and serum samples.

The kit has been validated using rat, mouse, equine, human, swine, and monkey samples. This assay employs a quantitative sandwich enzyme immunoassay technique, which measures GLP-2 in approximately 4 hours. A polyclonal antibody specific for GLP-2 has been pre-coated onto a 96-well microplate with removable strips. GLP-2 in standards and samples is sandwiched by the immobilized antibody and biotinylated polyclonal antibody specific for GLP-2, which is recognized by a streptavidin-peroxidase conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

**The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.**

**Notes** Glucagon-like peptide 2 (GLP-2) is a thirty-three amino acid peptide that is secreted by intestinal endocrine cells alongside GLP-1. GLP-2 and GLP-1 are co-secreted from enteroendocrine L-cells located in distal intestine in response to enteral nutrient ingestion, particularly fats and carbohydrates. GLP-2 secretion is mediated by direct nutrient stimulation of the L-cells and indirect action from enteroendocrine and neural inputs. GLP-2 acts to enhance nutrient absorption by inhibiting gastric motility and secretion and by stimulating nutrient transport.

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

## Properties

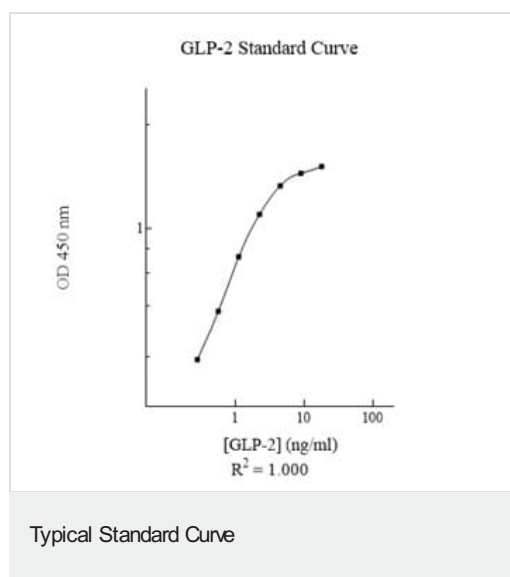
**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 20ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated GLP-2 Antibody	1 x 120µl
Anti-GLP-2 coated Microplate (12 x 8 wells)	1 unit
Chromogen Substrate	1 x 7ml
GLP-2 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

**Relevance** Human GLP2 (glucagon like peptide 2) is a peptide hormone of the glucagon family produced by the L cells of the intestinal mucosa and originates from the same prohormone as glucagon. GLP2 has been shown to be an important growth factor for the intestinal mucosa.

**Cellular localization** Secreted

## Images



Representative Standard Curve using ab222863

**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