

AMP'D[®] GLP-1 ELISA

kit

Ultra-sensitive AMP'D[®] GLP-1 ELISA kit enabling the ability to use less sample to detect levels of human Glucagon-like peptide 1 (7-36) amide, a potent promoter of insulin

secretion, a major incretin hormone, and therapeutic for type 2 diabetes.

GLP-1 functions in the regulation of blood glucose through insulin, and plays a role in obesity and diabetes. It is the most potent promoter of insulin secretion and an important gastrointestinal peptide hormone, or incretin. In response to eating, increased levels of GLP-1 stimulate insulin secretion and inhibit glucagon release leading to lower blood glucose levels. These positive effects of GLP-1 on glucose regulation make it an important target for the management of type 2 diabetes.

Ordering Information

[Order Online »](#)

ENZ-KIT104-0001

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Ultra-sensitive measurement of GLP-1, detecting as little as 5.5 pg/ml
- Negligible reactivity from similar GLP-1 forms
- High throughput format with results in 3 hours for up to 40 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Handling & Storage

Use/Stability Store all components at +4°C, except standard at -20°C.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Glucagon-like peptide-1

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of Glucagon-like peptide 1 (GLP-1) (7-36) amide in serum and plasma of human origin.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter Plate, Assay Buffer 13, Standard, Wash Buffer Concentrate, Antibody, Conjugate, and AMP'D[®] (Signal Amplification) Substrate, Amplifier, Substrate Diluent, Amplifier Diluent, and Stop Solution

Crossreactivity GLP-1 (7-36) amide (100%), GLP-1 (7-37) ($\leq 2.1\%$) and $\leq 0.02\%$: GLP-1 (9-36) amide, GLP-1 (1-36) amide, GLP-2, Glucagon

GenBank ID X03991.1

Gene/Protein Identifier NP_002045.1 (RefSeq), NM_002054.4 (RefSeq), 2641 (Entrez GeneID), 138030 (OMIM)

Sensitivity 5.5 pg/ml (range 7.8-250 pg/ml)

Species Reactivity Human

UniProt ID P01275[98-127]

Wavelength 495 nm

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)