

# AMP'D<sup>®</sup> GLP-1 ELISA

## kit

Ultra-sensitive AMP'D<sup>®</sup> 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 »](#)

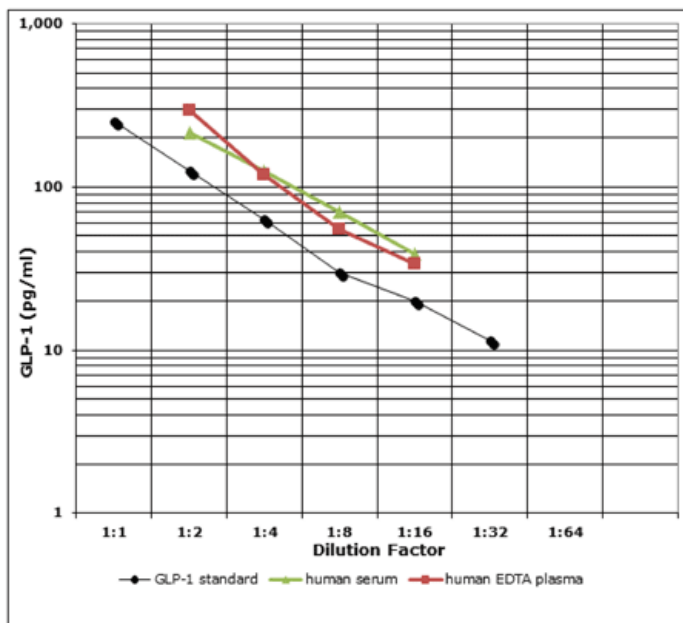
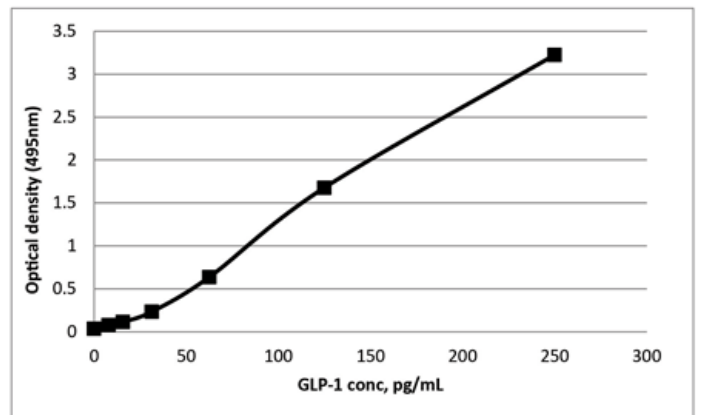
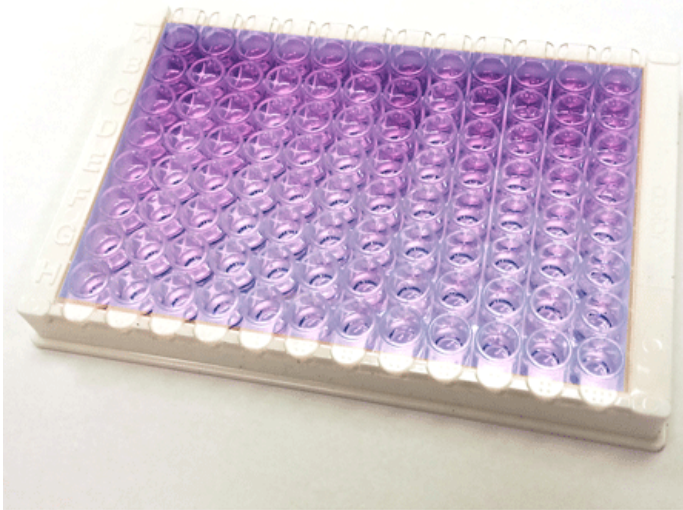
ENZ-KIT104-0001

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 <a href="#">Absorbance 96 Plate Reader</a> .
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) (≤2.1%) and ≤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



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)