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 SecAMP'D®, GaPn1ajd\$Arkitica reporterio immunoassay kit with results in 3 hours. 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 »

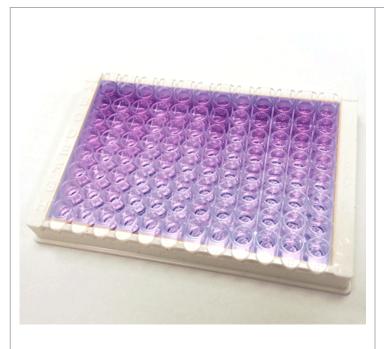
ENZ-KIT104-0001

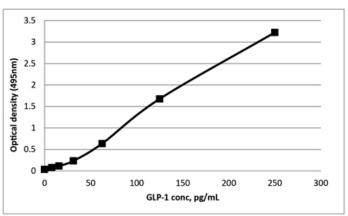
96 wells

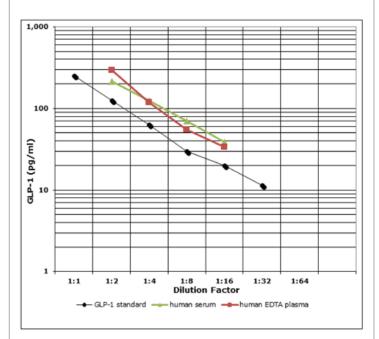
Manuals, SDS & CofA

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) (≤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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com