



# Mercodia

# Glucagon ELISA - 10 $\mu$ L

Directions for Use

**10-1281-01**

**Reagents for 96 determinations**

For Research Use Only  
Not for Use in Diagnostic Procedures

Manufactured by

**Mercodia AB**  
Sylveniusgatan 8A  
SE-754 50 Uppsala  
Sweden

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063





Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) info@eaglebio.com



**EAGLE**  
BIOSCIENCES

## Explanation of symbols used on labels

 $\Sigma = 96$	Reagents for 96 determinations
	Expiry date
	Store between 2-8°C
	Lot No.

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
BIOSCIENCES

## Intended Use

Mercodia Glucagon ELISA - 10  $\mu$ L provides a method for the quantitative determination of glucagon in rat, mouse and non-human primate (NHP) serum, EDTA plasma and cell culture media samples.

## Summary and explanation of the test

Glucagon is a 29 amino acid polypeptide processed from proglucagon in pancreatic alpha cells. In intestinal L-cells proglucagon is cleaved into glicentin, corresponding to proglucagon residues no 1-69. Glicentin can further be processed into oxyntomodulin, corresponding to proglucagon residues no 33-69. These peptides are released simultaneously upon stimulation. Moreover, a fragment of glucagon corresponding to its C-terminal part (residues no 19-29), also designated mini-glucagon, is reported to be present in the pancreas in low amounts compared to the total glucagon content.

In general, glucagon has an effect opposite that of insulin, i.e. it raises blood glucose levels. It causes the liver to convert glycogen into glucose, which is then released into the blood stream. With longer stimulation, glucagon action in the liver results in a glucose-sparing activation of free fatty acid oxidation and production of ketones. During hypoglycaemia, glucagon secretion offers a protective feedback mechanism, defending the organism against damaging effects of glucose deficiency in the brain and nerves.

## Principle of the procedure

Mercodia Glucagon ELISA - 10  $\mu$ L is a solid phase two-site enzyme immunoassay. It is based on the direct sandwich technique in which two monoclonal antibodies are directed against separate antigenic determinants on the glucagon molecule. During incubation glucagon in the sample reacts with peroxidase-conjugated anti-glucagon antibodies (clone E6A11K) and anti-glucagon antibodies (clone M5F9S) bound to microplate wells. A simple washing step removes unbound enzyme labelled antibody. The bound conjugate is detected by reaction with 3,3',5,5'-tetramethylbenzidine (TMB). The reaction is stopped by adding acid to give a colorimetric endpoint that is read spectrophotometrically.

## Warnings and precautions

- For research use only. Not for use in diagnostic procedures.
- Not for internal or external use in humans or animals.
- The Stop Solution in this kit contains 0.5M H<sub>2</sub>SO<sub>4</sub>. Follow routine precautions for handling hazardous chemicals.
- All samples should be handled as capable of transmitting infections.
- Each well can only be used once.
- The Stop Solution contains <5% Sulphuric acid.  
The Stop Solution is labeled:



### Danger

H318 – Causes serious eye damage.

H315 – Causes skin irritation.

P280 – Wear protective gloves. Wear eye or face protection.

P264 – Wash hands thoroughly after handling.

P302 + P352 + P362 + P364 – IF ON SKIN: Wash with plenty of soap and water.

Take off contaminated clothing and wash it before reuse.

P332 + P313 – If skin irritation occurs: Get medical attention.

P305 + P351 + P338 + P310 – IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Immediately call a POISON CENTER or physician.

- The Enzyme Conjugate Buffer, Cal 0, 1, 2, 3, 4, 5 and Wash Buffer contain <0.06% 5-chloro-2-methyl-4-isothiazolin-3-one and 2-methyl-2H-isothiazol-3-one (3:1).

The Enzyme Conjugate Buffer, the Calibrators and Wash Buffer are labeled:



### Warning

H317 – May cause an allergic skin reaction.

P280 – Wear protective gloves.

P261 – Avoid breathing vapour.

P272 – Contaminated work clothing should not be allowed out of the workplace.

P302 + P352 – IF ON SKIN: Wash with plenty of soap and water.

P333 + P313 – If skin irritation or rash occurs: Get medical attention.

P501 – Dispose of contents and container in accordance with all local, regional, national and international regulations.

## Material required but not provided

- Pipettes with appropriate volumes (repeating pipettes preferred for addition of enzyme conjugate 1X solution, Substrate TMB and Stop Solution)
- Tubes, beakers and cylinders for reagent preparation
- Redistilled water
- Magnetic stirrer
- Vortex mixer
- Microplate reader with 450 nm filter
- Microplate shaker (700–900 cycles per minute, orbital movement)
- Refrigerator (2–8°C) with room for microplate shaker
- Microplate washing device with overflow function (recommended but not required)

## Reagents 1 X 96

Each Mercodia Glucagon ELISA – 10 µL kit (10-1281-01) contains reagents for 96 wells, sufficient for 42 samples and one Calibrator curve in duplicate. For larger series of assays, use pooled reagents from packages bearing identical lot numbers. The expiry date for the complete kit is stated on the outer label. The recommended storage temperature is 2–8°C.

<b>Coated Plate</b> Mouse monoclonal anti-glucagon For unused microplate strips, reseal the bag using adhesive tape, store at 2–8°C and use within 2 months.	1 plate	96 wells 8-well strips	Ready for Use
<b>Calibrators 1, 2, 3, 4, 5</b> Synthetic glucagon Color coded yellow Concentration stated on vial label. Storage after reconstitution: 2–8°C for 1 month. For storage of reconstituted Calibrators for more than 1 month, store at -20°C.	5 vials	1000 µL	Lyophilized Add 1000 µL redistilled water per vial.
<b>Calibrator 0</b> Color coded yellow	1 vial	5 mL	Ready for Use
<b>Enzyme Conjugate 11X</b> Mouse monoclonal anti-glucagon	1 vial	0.6 mL	Preparation, see below
<b>Enzyme Conjugate Buffer</b> Color coded blue	1 vial	6 mL	Ready for Use
<b>Wash Buffer 21X</b> Storage after dilution: 2–8°C for 2 months.	1 bottle	50 mL	Dilute with 1000 mL redistilled water to make wash buffer 1X solution.
<b>Substrate TMB</b> Colorless solution <i>Note! Light sensitive!</i>	1 bottle	22 mL	Ready for Use
<b>Stop Solution</b> 0.5 M H <sub>2</sub> SO <sub>4</sub>	1 vial	7 mL	Ready for Use

## Preparation of enzyme conjugate 1X solution

Prepare the needed volume of enzyme conjugate 1X solution by dilution of Enzyme Conjugate 11X (1+10) in Enzyme Conjugate Buffer according to the table below. When preparing enzyme conjugate 1X solution for the whole plate, pour all of the Enzyme Conjugate Buffer into the Enzyme Conjugate 11X vial. Mix gently. Use within 1 week.

Number of strips	Enzyme Conjugate 11X	Enzyme Conjugate Buffer
12 strips	1 vial	1 vial
8 strips	0.36 mL	3.6 mL
4 strips	0.18 mL	1.8 mL

## Specimen collection and handling

Serum or EDTA plasma can be used. However, glucagon in serum or EDTA plasma samples will be sensitive to storage conditions and freeze-thaw cycles. Addition of aprotinin to samples will not improve sample stability. It is recommended to keep samples on ice when handling them at room temperature.

### Serum

Collect blood by venipuncture, allow to clot, and separate the serum by centrifugation. Avoid storing samples at room temperature or 2-8°C for longer than 2 hours. Store samples at -80°C and avoid freeze-thaw cycles.

### Plasma

#### EDTA plasma

Collect blood by venipuncture into tubes containing EDTA as anticoagulant, and separate the plasma fraction by centrifugation. Avoid storing samples at room temperature or 2-8°C for longer than 2 hours. Store samples at -80°C and avoid freeze-thaw cycles.

#### EDTA plasma in BD P800 tubes

For studies in which very low levels of glucagon need to be detected, it may be beneficial to use BD (Becton Dickinson) P800 tubes for sample collection, since this will prevent the degradation of glucagon. Avoid storing samples at room temperature or 2-8°C for longer than 6 hours. Store samples at -80°C and avoid freeze-thaw cycles.

### Cell culture medium

Note that different chemicals used in cell culture media can interfere with the assay (such as sodium azide ( $\text{NaN}_3$ ) and beta-mercaptoethanol).

### Preparation of samples

No dilution is normally required, however, samples above the obtained value of Calibrator 5 should be diluted with Calibrator 0. Samples that appear cloudy benefit from centrifugation before pipetting. *Note!* Buffers containing sodium azide ( $\text{NaN}_3$ ) can not be used for sample dilution.

## Test procedure

Prepare a calibrator curve for each assay run.

1. Prepare enzyme conjugate 1X solution and wash buffer 1X solution.
2. Prepare sufficient microplate wells to accommodate Calibrators, controls and samples in duplicate.
3. Pipette 10  $\mu$ L each of Calibrators, controls and samples into appropriate wells.
4. Add 50  $\mu$ L enzyme conjugate 1X solution to each well and attach the plate sealer.
5. Incubate on a plate shaker (700-900 rpm) over night (18-22h) at 2-8°C.
6. Wash 6 times with 700  $\mu$ L wash buffer 1X solution per well using an automatic plate washer with overflow-wash function. After final wash, invert and tap the plate firmly against absorbent paper. Do not include soak step in washing procedure.  
Or manually:  
Discard the reaction volume by inverting the microplate over a sink. Add 350  $\mu$ L wash solution to each well. Discard the wash buffer 1X solution, tap firmly several times against absorbent paper to remove excess liquid. Repeat 5 times. Avoid prolonged soaking during washing procedure.
7. Add 200  $\mu$ L Substrate TMB.
8. Incubate on the bench for 30 minutes at room temperature (18-25°C).
9. Add 50  $\mu$ L Stop Solution to each well.  
Place plate on a shaker for approximately 5 seconds to ensure mixing.
10. Read optical density at 450 nm and calculate results.  
Read within 30 minutes.

*Note!* Be extra careful not to contaminate the Substrate TMB with enzyme conjugate solution.

## Internal quality control

Commercial controls and/or internal serum pools with low, intermediate and high glucagon concentrations should routinely be assayed as samples, and results charted from day to day. It is good laboratory practice to record the following data for each assay: kit lot number, dilution and/or reconstitution dates of kit components, OD values for the blank, Calibrators and controls.

## Calculation of results

The concentration of glucagon is obtained by plotting the absorbance of the calibrators, except for Calibrator 0, versus their concentration. It is important to use an appropriate curve fitting model that represent the true dose-response relationship to get accurate results. It is every laboratory's responsibility to try out the functionality of the chosen curve fitting model and used software. Note that weighting of the curve fit is important to get a proper fit.

The Mercodia Glucagon ELISA - 10  $\mu$ L is validated with Four parameter logistic with weighting  $1/s^2$ , using Magellan (Tecan) Software.

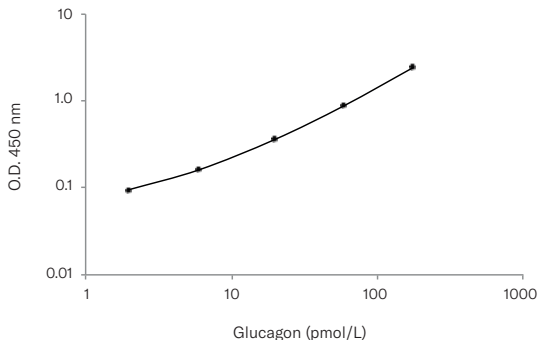
## Example of results

Wells	Identity	A <sub>450</sub> nm	Mean conc. pmol/L
1A-B	Calibrator 0	0.069/0.076	
1C-D	Calibrator 1*	0.095/0.095	
1E-F	Calibrator 2*	0.154/0.169	
1G-H	Calibrator 3*	0.360/0.363	
2A-B	Calibrator 4*	0.841/0.925	
2C-D	Calibrator 5*	2.468/2.408	
2E-F	Sample 1	0.178/0.176	7.218
2G-H	Sample 2	0.355/0.373	20.51
3A-B	Sample 3	0.916/1.007	64.36

\* Concentration stated on vial label.

## Example of calibrator curve

A typical calibrator curve is shown here. Do not use this curve to determine actual assay results.





## Conversion factor

1 pmol/L = 3.5 pg/mL

## Limitations of the procedure

Grossly lipemic or icteric samples do not interfere in the assay. High levels of hemoglobin (>500 mg/dL) can interfere in the assay.

## Expected values

Good practice dictates that each laboratory establishes its own expected range of values.

## Performance characteristics

### Detection limit

Detection limit is defined as the Capability of Detection according to ISO11843-Part 1. Capability of Detection should be seen as part of a method validation, rather than the lowest concentration that can be measured.

The detection limit is 1.5 pmol/L as determined with the methodology described in ISO11843-Part 4.

Concentrations of samples with absorbance below Calibrator 1 should not be calculated, but expressed as less than or equal to ( $\leq$ ) the concentration indicated on the vial for Calibrator 1.

### Recovery

Species	Addition			Dilution		
	Min	Max	Mean	Min	Max	Mean
NHP*	96	112	103	85	106	95
Mouse	109	126	118	78	101	85
Rat	114	121	118	71	125	92

Figures in %

\*Non-human primate samples kindly provided by Professor Barbara C. Hansen at University of South Florida, USA.

## Hook effect

Samples with a concentration up to at least 8 µmol/L can be measured without giving falsely low results.

## Precision

Each sample was analyzed in 4 replicates on at least 7 different occasions.

Species	Sample	Mean value pmol/L	Coefficient of variation	
			Repeatability %*	Within laboratory %**
NHP	1	12.6	7.2	12.3
	2	36.9	7.4	12.3
	3	44.7	7.7	10.6
Rat	1	7.3	5.1	10.1
	2	8.8	6.9	16.1
	3	9.6	4.4	17.9

\*Within assay variation

\*\*Total assay variation

## Specificity

	Crossreaction	Highest concentrations tested (pmol/L)
Mini-glucagon	n.d.	120
Glicentin, human	1.0%	300
Glicentin, mouse	7.0%	135
Glicentin, rat	4.0%	160
Oxyntomodulin, human/rat/mouse	2.0%	400
Oxyntomodulin, bovine/canine/porcine	n.d.	400
GLP-1 (7-36)	n.d.	100
GLP-1 (9-36)	n.d.	600
GLP-2	n.d.	100

n.d. = not detectable

## Calibration

Mercodia Glucagon ELISA – 10 µL is calibrated against WHO 1<sup>st</sup> International reference preparation 69/194.

## Warranty

The performance data presented here was obtained using the procedure indicated. Any change or modification in the procedure not recommended by Mercodia AB may affect the results, in which event Mercodia AB disclaims all warranties expressed, implied or statutory, including the implied warranty of merchantability and fitness for use.

Mercodia AB and its authorized distributors, in such event, shall not be liable for damages indirect or consequential.

## References

Bagger JI *et al* (2011) Glucagon antagonism as a potential therapeutic target in type 2 diabetes. *Diabetes Obes Metab* 13:965-971.

Holst J *et al* (2004) Role of incretin hormones in the regulation of insulin secretion in diabetic and nondiabetic humans. *Am J Physiol Endocrinol Metab* 287:E199-206.

Holst J (2010) Glucagon and Glucagon-Like Peptides 1 and 2. *Results Probl Cell Differ* 50:121-135.

Young A (2005) Inhibition of Glucagon Secretion. *Adv Pharmacol* 52:151-171.

Further references can be found on our website: [www.mercodia.com](http://www.mercodia.com)

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
BIOSCIENCES









**Summary of protocol sheet**  
**Mercodia Glucagon ELISA - 10 µL**

Add Calibrators, controls* and samples	10 µL
Add enzyme conjugate 1X solution and attach plate sealer	50 µL
Incubate	Over night (18-22 h) at 2-8°C on a plate shaker, 700-900 rpm
Wash plate with wash buffer 1X solution	700 µL, 6 times
Add Substrate TMB	200 µL
Incubate	30 minutes at 18-25°C
Add Stop Solution	50 µL Shake for 5 seconds to ensure mixing
Measure $A_{450}$ nm	Evaluate results

\*not provided

For full details see page 7

For technical support please contact: support@merckodia.com

distributed in the US/Canada by:  
**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
 BIOSCIENCES

31-3189  
 Version 6.0  
 2020-01-20