



Northern Lights Mercodia Lispro NL-ELISA

Directions for Use

10-1291-01

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 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.

Intended use

Mercodia Lispro NL-ELISA provides a method for the quantitative determination of insulin lispro in plasma or serum samples.

Summary and explanation of the test

The Mercodia Lispro NL-ELISA is a ligand binding assay able to measure insulin lispro specifically, without cross-reaction to native insulin, native proinsulin, or any of the tested insulin analogues. Also, no interference from insulin autoantibodies (IgG antibodies), which can be present in samples from diabetic patients, has been observed¹.

Mercodia Lispro NL-ELISA is designed to fit the needs of the pharmaceutical industry, by meeting the requirements specified in the EMA²/FDA³ guidelines. Seven serum calibrators (incl. Calibrator 0), two anchor points and three controls are included in the kit.

Principle of the procedure

Mercodia Lispro NL-ELISA is a solid phase two-site enzyme immunoassay based on the sandwich technique, in which two monoclonal antibodies are directed against separate antigenic determinants on the lispro molecule. Insulin lispro in the sample reacts with anti-lispro antibodies bound to microtitration wells and peroxidase-conjugated anti-lispro antibodies in the solution. A simple washing step removes unbound enzyme-labelled antibody. The bound conjugate is detected by reaction with the chemiluminescent substrate. A chemiluminescence plate reader is used to read the intensity of light generated.

Warnings and precautions

- For research use only. Not for use in diagnostic procedures.
- Instrument settings should be optimized according to the manufacturer's instructions.
- Not for internal or external use in humans or animals.
- All samples should be handled as capable of transmitting infections.
- Each well can only be used once.
- The Enzyme Conjugate Buffer, Calibrator 0, 1, 2, 3, 4, 5, and 6, Anchor Points (Low and High), Controls (Low, Medium, High), Assay Buffer 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, Anchor Points (Low and High), Controls (Low, Medium, High), Assay Buffer 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.

Warning! This kit contains reagents that may be infectious!

The serum Calibrators, Controls and Anchor Points are manufactured from human blood components. The source materials have been tested by immunoassay for Hepatitis B surface antigen, antibodies for Hepatitis C virus and for antibodies for HIV virus and found to be negative.

Nevertheless, all recommended precautions for the handling of blood derivatives should be observed as no test method can provide total assurance that these viruses or other contagion are absent. Please refer to HHS Publication No. (CDC) 88-8395 or corresponding local/national guidelines on laboratory safety procedures.

Material required but not provided

- Pipettes with appropriate volumes (multichannel or repeating pipettes preferred for addition of Assay Buffer, enzyme conjugate 1X solution and substrate reagent solution)
- Tubes, beakers and cylinders for reagent preparation
- Redistilled water for resuspension of calibrators and controls
- Magnetic stirrer
- Vortex mixer
- Microplate reader for chemiluminescence (glow)
- Microplate shaker (700–900 cycles per minute. orbital movement)
- Microplate washing device with overflow function (recommended but not required)

Reagents 1 X 96

Each Mercodia Lispro NL-ELISA kit (10-1291-01) contains reagents for 96 wells, sufficient for 36 samples, 3 controls 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.

Coated Plate Mouse monoclonal anti-lispro For unused microplate wells completely reseal the bag using adhesive tape, store at 2–8°C and use within 4 weeks.	1 plate	96 wells 8-well strips	Ready for Use
Calibrators 1, 2, 3, 4, 5, 6 Lispro Serum-based Concentration stated on vial label Reconstituted Calibrators are stable for 1 month at 2–8°C.	6 vials	500 µL	Lyophilized Resuspend with 500 µL redistilled water per vial
Calibrator 0 Serum-based Reconstituted Calibrators are stable for 1 month at 2–8°C.	1 vial	500 µL	Lyophilized Resuspend with 500 µL redistilled water per vial
Anchor Points Low, High Lispro Serum-based Concentration stated on on vial label Reconstituted Anchor Points are stable for 1 month at 2–8°C.	2 vials	500 µL	Lyophilized Resuspend with 500 µL redistilled water per vial
Controls Low, Medium, High Serum-based Concentration stated on on vial label Reconstituted Controls are stable for 1 month at 2–8°C.	3 vials	500 µL	Lyophilized Resuspend with 500 µL redistilled water per vial
Assay Buffer Color coded red	1 vial	14 mL	Ready for Use

Enzyme Conjugate 11X Peroxidase conjugated mouse monoclonal anti-insulin	1 vial	1.3 mL	Preparation see below
Enzyme Conjugate Buffer Color coded blue	1 vial	13 mL	Ready for use
Wash Buffer 21X 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
Substrate NL Ultra A Colorless solution	1 vial	2 mL	Preparation see below
Substrate NL Ultra B Colorless solution	1 vial	4 mL	Preparation see below

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 the blue Enzyme Conjugate Buffer into the Enzyme Conjugate 11X vial. Mix gently. Store at 2-8°C and use within 4 weeks.

Number of strips	Enzyme Conjugate 11X	Enzyme Conjugate Buffer
12 strips	1 vial	1 vial
8 strips	700 µL	7000 µL
4 strips	350 µL	3500 µL

Preparation of substrate working solution

Prepare the needed volume of substrate working solution by mixing Substrate NL Ultra A with Substrate NL Ultra B (1:2) according to the table below. Mix gently. Store at 2-8°C and use within 1 day. *Note!* Protect substrate from light.

Number of strips	Substrate NL Ultra A	Substrate NL Ultra B
12 strips	1 vial	1 vial
8 strips	1.2 mL	2.4 mL
4 strips	0.6 mL	1.2 mL

Specimen collection and handling

Serum, EDTA plasma and P-800 plasma can be used. Store samples at -80°C and avoid freeze-thaw cycles. Up to three freeze-thaw cycles for serum and plasma samples had minimal effect on lispro levels in samples. Avoid long time storage of samples at room temperature or 2-8°C. Lispro in serum and plasma samples was found to be stable at room temperature for 4 hours and at 2-8°C for 24 hours.

Serum

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

EDTA plasma and P-800 plasma

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

Preparation of samples

No dilution is normally required, however, samples above the obtained value of Calibrator 6 should be diluted with Calibrator 0 or a serum pool. *Note!* Buffers containing sodium azide (NaN_3) cannot be used for sample dilution.

Test procedure

All reagents and samples must be brought to room temperature before use. Assay a calibrator curve in each run. Keep substrate and conjugate solution separated from each other during all steps to avoid contamination. The product has been optimized and validated without plate sealer.

1. Prepare enzyme conjugate 1X solution, substrate working solution and wash buffer 1X solution.
2. Resuspend Calibrators, Anchor Points and Controls with 500 μ L redistilled water.
3. Prepare sufficient microplate wells to accommodate Calibrators, Anchor Points, Controls and samples in duplicate.
4. Pipette 10 μ L each of Calibrators, Anchor Points, Controls and samples into appropriate wells.
5. Add 100 μ L of Assay Buffer into each well.
6. Incubate on plate shaker for 1 hour (700-900 rpm) at room temperature (18-25°C).
7. On an automatic plate washer, use plate mode combined with overflow-wash function to wash the plate with 700 μ L wash buffer 1X solution per well for 6 cycles. Invert and tap the plate firmly against absorbent paper after the final wash.
Or manually,
discard the reaction volume by inverting the microplate over a sink. Add 350 μ L wash buffer 1X solution to each well. Discard the wash buffer 1X solution, tap firmly several times against absorbent paper to remove excess liquid. Repeat 5 times for a total of 6 cycles. See TN34-0106 for details.
8. Add 100 μ L enzyme conjugate 1X solution to each well.
9. Incubate on plate shaker for 1 hour (700-900 rpm) at room temperature (18-25 °).
10. Wash as described in step 7.
11. Add 50 μ L substrate working solution into each well.
12. Incubate for 10 minutes (18-25°C).
13. Read the luminosity of the wells using a 1 second integration time, without using a filter. If the reader has manual gain, gain should be set so that the signal of Anchor Point High is within the readers dynamic range. Read within 15 minutes.
14. Concentrations of Controls and samples should be read/calculated using Anchor Points and Calibrator 1-6 with a 5-pl weighted fit (1/Y²). *Note!* Calibrator 0 should not be included in the curve fit.

Internal quality controls

Serum controls included in the kit and/or internal plasma/serum pools with low, intermediate and high insulin lispro concentrations should routinely be assayed as samples and results should be charted from day to day. It is good laboratory practice to record the following data for each assay: kit lot number, preparation dates of kit components, calibrators and controls.

Laboratories should follow government regulations or accreditation requirements for quality control frequency.

The concentration ranges stated on the control vials are based on the nominal value of insulin lispro with acceptance criteria $100\% \pm 20\%$ as recommended by EMA²/FDA³.

Calculation of results

The concentration of lispro is obtained by plotting the relative light units (RLU) of the Calibrators, except for Calibrator 0, versus their concentration. It is important to use an appropriate curve fitting model that represents 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 at the low range of the standard curve, especially when the measuring range is wide.

The Mercodia Lispro NL-ELISA is validated using MARS (BMG Labtech) with Five Parameter Logistic (5PL) and automatic weighting using $1/Y^2$.

Note!

- Include the Anchor Points in the curve fit.
- Do not include Calibrator 0 in the curve fit.

Example of results

These values were obtained using BMG Labtech Clariostar with 1 s integration time, 2500 gain and 11.5 mm focal height.

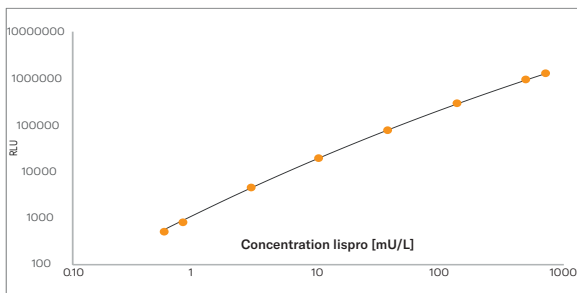
Wells	Identity	Mean RLU**	Mean conc. mU/L
1 A-D	Calibrator 0	66	<Anchor Point Low
	Anchor Point Low*	649	0.71
1 E-H	Calibrator 1*	1109	1.06
2 A-D	Calibrator 2*	4887	3.38
2 E-H	Calibrator 3*	21279	11.6
3 A-D	Calibrator 4*	81047	38.8
3 E-H	Calibrator 5*	306242	143
4 A-D	Calibrator 6*	973581	495
4 E-H	Anchor Point High*	1327796	704
8 A-D	Control Low*	2524	2.0
8 E-H	Control Medium*	46396	23.2
9 A-D	Control High*	806775	401

*Concentration stated on vial label.

**The RLU is affected by reader model and gain settings. Different readers can give different RLU values. Absolute RLU values do not affect calculated results.

Calibrator curve

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



Conversion factor

1 mU/L = 6 pmol/L = 0,035 µg/L = 35 pg/mL

Molecular weight of insulin lispro is: 5808 g/mol

Limitations of the procedure

The assay is specific for insulin lispro with no cross-reactivity to any of the tested insulin analogs, insulin autoantibodies, native insulin and native proinsulin as shown in performance characteristics.

Expected values

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

Performance characteristics

The assay has been validated according to guidelines. Selected studies are presented here. Additional data can be obtained from Mercodia.

Validation of curve fit

The curve fitting was validated with Five Parameter Logistics (weighted fit $1/Y^2$).

Sensitivity and range of quantification

Capability of Detection is 0.36 mU/L as determined by the methodology described in ISO11843-Part 4⁴.

Lower Limit of Quantification, LLOQ, is 1.0 mU/L as determined according to EMA²/FDA³ guidelines.

The Upper Limit of Quantification, ULOQ, is 500 mU/L as determined according to EMA²/FDA³ guidelines.

Precision and accuracy

Human serum

Human serum controls spiked with Humalog® (lispro) were analyzed in 4 replicates on at least 12 different occasions by 2 different technicians on 4 days.

Sample	Mean value mU/L	Between run- Accuracy %	Coefficient of variation	
			Repeatability %*	Within laboratory %**
QC _{LOQ}	0.94	94	6.7	7.7
QC _{Low}	1.89	95	3.5	5.8
QC _{Medium}	21.4	86	2.2	3.7
QC _{High}	364	91	1.9	3.0
QC _{ULOQ}	459	92	2.5	3.4

*Within assay variation

**Total assay variation

Human EDTA Plasma

Human EDTA plasma, human P-800 plasma and porcine EDTA plasma were spiked with EDQM Lispro⁵ and analyzed in 4 replicates on 6 occasions by 2 different technicians on 2 days.

Sample	Mean value mU/L	Between run- Accuracy %	Coefficient of variation	
			Repeatability %*	Within laboratory %**
QC _{LOQ}	0.98	98	2.4	3.6
QC _{Low}	1.95	98	3.7	5.2
QC _{Medium}	22.5	90	2.1	4.0
QC _{High}	333	83	2.1	3.2
QC _{ULOQ}	415	83	3.5	4.7

*Within assay variation

**Total assay variation

Human P-800 Plasma

Sample	Mean value mU/L	Between run- Accuracy %	Coefficient of variation	
			Repeatability %*	Within laboratory %**
QC _{LOQ}	1.07	107	3.3	3.4
QC _{Low}	2.08	104	1.3	4.1
QC _{Medium}	22.2	89	1.4	2.7
QC _{High}	336	84	1.1	1.7
QC _{ULOQ}	417	83	1.4	3.5

*Within assay variation

**Total assay variation

Porcine EDTA Plasma

Sample	Mean value mU/L	Between run- Accuracy %	Coefficient of variation	
			Repeatability %*	Within laboratory %**
QC _{LLOQ}	1.09	109	2.1	4.3
QC _{Low}	2.00	100	1.8	3.85
QC _{Medium}	22.3	89	2.1	2.96
QC _{High}	332	83	0.8	1.9
QC _{ULOQ}	424	85	1.7	3.41

*Within assay variation

**Total assay variation

Selectivity

Selectivity has been validated according to FDA³/EMA² guidelines. At least 10 blank samples from different individuals of each tested matrix (human serum, human EDTA plasma, human P-800 plasma and porcine EDTA-plasma) showed no positive signal in the assay in the absence of lispro. Auto-insulin antibodies, HAMA, rheumatoid factor (RF), grossly lipemic, icteric or hemolyzed samples do not interfere in the assay.

Parallelism

Samples with various concentration of insulin lispro were serially diluted 1/2, 1/4 and 1/8 with Calibrator 0 as dilution matrix. Patient samples (human, porcine) dosed with lispro were parallel to the serum calibrator curve, with acceptance criteria CV < 30 % between back-calculated samples according to EMA² guidelines.

High dose hook effect

Samples with a concentration up to 10 000 mU/L (20X Cal 6) can be measured without giving falsely low results.

Specificity

Cross-reactivity

The following cross-reactions and interferences were studied:

Substance	Concentrations tested	Cross-reaction	Interference (Acceptance criteria 100 ± 25 %)	
			LLOQ (Recovery %)	ULOQ (Recovery %)
Native Human Insulin	50-400 mU/L	N.D.	105-112	97-114
Native Human Proinsulin	50-300 pmol/L	N.D.	104-113	103-110
Glargine	50-600 mU/L	N.D.	96-117	89-107
Glargine M1	50-600 mU/L	N.D.	98-105	100-107
Glargine M2	50-600 mU/L	N.D.	104-123	109-115
Degludec	50-600 mU/L	N.D.	101-113	103-109
Detemir	50-600 mU/L	N.D.	107-116	108-110
Insulin NPH	50-600 mU/L	N.D.	106-116	102-108
Aspart	50-600 mU/L	N.D.	95-111	101-106
Glulisine	50-600 mU/L	N.D.	101-112	99-103
Native Porcine Insulin	50-600 mU/L	N.D.	-	-
Native Porcine Proinsulin	50-300 pmol/L	N.D.	-	-

N.D. = Not Detected

Calibration

Mercodia Lispro NL-ELISA is calibrated against a highly purified, fully validated, commercial European Pharmacopoeia Reference Standards: Insulin Lispro CRS4. Accuracy of reference standard was evaluated by measuring known concentrations of insulin lispro from Humalog® (Eli Lilly and Company).

Warranty

The performance data presented here was obtained using the procedure indicated. Any change or modification in the procedure not recommended by Merckodia AB may affect the results, in which event Merckodia AB disclaims all warranties expressed, implied or statutory, including the implied warranty of merchantability and fitness for use. Merckodia AB and its authorized distributors, in such event, shall not be liable for damages indirect or consequential.

References

1. Bingley, P. J. et al. Combined analysis of autoantibodies improves prediction of IDDM in islet cell antibody-positive relatives. *Diabetes* 43, 1304-10 (1994)
2. Guideline on bioanalytical method validation, European Medicines Agency (EMA), Science Medicines Health, 21 July 2011.
3. Guidance for Industry, Bioanalytical Method Validation, U.S. Department of Health and Human Services, Food and Drug Administration, Center for Drug Evaluation and Research (CDER), Center for Veterinary Medicine (CVM), May 2018, Biopharmaceutics.
4. International Organization for Standardization, www.iso.org/standard, ISO11843-Part 4
5. European Directorate for the Quality of Medicines & HealthCare European Pharmacopoeia (Ph. Eur.). Reference Standard, INSULIN LISPRO CRS batch 2. Catalogue code: Y0000348 (<https://crs.edqm.eu>)

Further references can be found on our website: **www.mercodia.com**.

Summary of protocol sheet
Northern Lights Mercodia Lispro NL-ELISA

Add calibrators, anchor points, controls and samples	10 μ L
Add Assay Buffer to all wells	100 μ L
Incubate	1 h at 18-25°C on a plate shaker (700-900 rpm)
Wash plate with wash buffer 1X solution	700 μ L, 6 times
Add enzyme conjugate solution to all wells	100 μ L
Incubate	1 h at 18-25°C on a plate shaker (700-900 rpm)
Wash plate with wash buffer 1X solution	700 μ L, 6 times
Add substrate solution	50 μ L
Incubate	10 minutes at 18-25°C on the bench
Read chemiluminescence	1 s integration time (glow)

For full details see page 8

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 • info@eaglebio.com



EAGLE
BIOSCIENCES

31-3196
Version 3.0
2021-01-27