

User Guide

# Rat Leptin ELISA Kit

## 96-Well Plate

### EZRL-83K

Intended Use.....	2	Graph of Typical Reference Curve ..	13
Principles of Assay.....	2	Correlation Graph .....	14
Reagents Supplied.....	3	Assay Characteristics.....	15
Storage and Stability .....	4	Sensitivity .....	15
Reagent Precautions .....	4	Specificity .....	15
Sodium Azide.....	4	Precision .....	15
Hydrochloric Acid .....	4	Spike and Recovery of Rat Leptin	
Symbol Definitions.....	5	in Rat Serum .....	16
Materials Required.....	6	Linearity .....	17
Sample Collection and Storage .....	7	Quality Controls .....	18
Rat Leptin ELISA Assay Procedure.....	8	Troubleshooting .....	18
Assay Procedure for Rat Leptin		Product Ordering.....	19
ELISA Kit .....	10	Replacement Reagents.....	19
Microtiter Plate Arrangement.....	11	Notice .....	20
Calculations.....	12	Technical Assistance .....	20
Interpretation .....	12	Terms and Conditions of Sale .....	20
		Contact Information.....	20

---

## Intended Use

This kit is used for the non-radioactive quantification of leptin in rat sera. Plasma samples may also be used but application to samples of other biological fluids may need validation by the user. One kit is sufficient to measure 37 unknown samples in duplicate.

**This kit is for Research Use Only. Not for use in Diagnostic Procedures.**

## Principles of Assay

This assay is a Sandwich ELISA based, sequentially, on:

- Binding of leptin in the sample by a pre-titered antiserum and immobilization of the resulting complexes in the wells of a microtiter plate
- After washing purified biotinylated detection antibody is allowed to bind to the immobilized leptin
- Binding of horseradish peroxidase to the immobilized biotinylated antibodies after free detection antibodies are washed off
- Wash away of free enzyme conjugates
- Quantification of immobilized antibody-enzyme conjugates by monitoring horseradish peroxidase activities in the presence of the substrate 3,3',5,5'-tetramethylbenzidine.

The enzyme activity is measured spectrophotometrically by the increased absorbency at 450 nm, corrected from the absorbency at 590nm, after acidification of formed products. Since the increase in absorbency is directly proportional to the amount of captured leptin in the unknown sample, the latter can be derived by interpolation from a reference curve generated in the same assay with reference standards of known concentrations of rat leptin.

## Reagents Supplied

Each kit is sufficient to run one 96-well plate and contains the following reagents:

**Note:** Store all reagents at 2-8 °C

Reagents Supplied	Volume	Quantity	Cat. No.
Rat/Mouse Leptin ELISA Plate with 2 plate sealers <b>Note:</b> Unused strips should be resealed in the foil pouch with the desiccant provided and stored at 2-8 °C	-	1 plate 2 sealers	EP83
Rat/Mouse Leptin Antiserum	6 mL	1 vial	EAS83
Rat Leptin Standards Rat leptin in buffer: 0.2, 0.5, 1, 2, 5, 10, 20, and 30 ng/mL	0.5 mL	1 vial per standard point	E8083-K
Rat Leptin Quality Controls 1 and 2	0.5 mL	1 vial each	E6083-K
Rat/Mouse Leptin Matrix Solution	0.5 mL	1 vial	EPS0016
Assay Buffer	40 mL	1 bottle	EAB-PTR
10X HRP Wash Buffer Concentrate	50 mL	2 bottles	EWB-HRP
Rat/Mouse Leptin Detection Antibody	12 mL	1 bottle	E1083
Enzyme Solution	12 mL	1 bottle	EHRP-4
Substrate Solution <b>Note:</b> Minimize light exposure.	12 mL	1 bottle	ESS-TMB2
Stop Solution <b>Caution:</b> Corrosive Solution	12 mL	1 bottle	ET-TMB

---

## Storage and Stability

Recommended storage for kit components is 2-8 °C.

All components are shipped and stored at 2-8 °C. Once opened, liquid standards and controls can be stored up to 30 days at 2-8 °C. Refer to expiration dates on all reagents prior to use. Do not mix reagents from different kits unless they have the same lot numbers.

## Reagent Precautions

### Sodium Azide







Sodium azide or Proclin™ has been added to some reagents as a preservative. Although the concentrations are low, Sodium azide and Proclin™ may react with lead and copper plumbing to form highly explosive metal azides. Dispose of unused contents and waste in accordance with international, federal, state, and local regulations.

### Hydrochloric Acid

Hydrochloric acid is corrosive and can cause eye and skin burns. Harmful if swallowed. Causes respiratory and digestive tract burns. Avoid contact with skin and eye. Do not swallow or ingest.

**Note:** See next page for full hazardous labels for the components of this kit.

## Symbol Definitions

Ingredient	Cat. No.	Full Label
Rat Leptin Detection Antibody	E1083	 <b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
Rat Leptin Quality Controls 1 & 2	E6083-K	 <b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
Rat Leptin Standard	E8083-K	 <b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
Antiserum for Rat/Mouse Leptin	EAS83	 <b>Warning:</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
Stop Solution	ET-TMB	 <b>Warning:</b> May be corrosive to metals.
10X HRP Wash Buffer Concentrate	EWB-HRP	 <b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.

---

## Materials Required (Not Provided)

- Multi-channel Pipettes and pipette tips: 0  $\mu$ L-50  $\mu$ L and 50  $\mu$ L-300  $\mu$ L
- Pipettes and pipette tips: 10  $\mu$ L-20  $\mu$ L or 20  $\mu$ L-100  $\mu$ L
- Buffer and Reagent Reservoirs
- Vortex Mixer
- De-ionized water
- Microtiter Plate Reader capable of reading absorbency at 450 nm and 590 nm
- Orbital Microtiter Plate Shaker
- Absorbent Paper or Cloth

---

## Sample Collection and Storage

1. To prepare serum samples, whole blood is directly drawn into a centrifuge tube that contains no anti-coagulant. Let blood clot at room temperature for 30 min.
2. Promptly centrifuge the clotted blood at 2,000 to 3,000  $\times g$  for 15 minutes at  $4 \pm 2^\circ\text{C}$ .
3. Transfer and store serum samples in separate tubes. Date and identify each sample.
4. Use freshly prepared serum or aliquot and store samples at  $\leq -20^\circ\text{C}$  for later use. Avoid multiple ( $> 3$ ) freeze/thaw cycles.
5. To prepare plasma sample, whole blood should be collected into centrifuge tubes containing enough  $\text{K}_3\text{EDTA}$  to achieve a final concentration of 1.735 mg/mL and centrifuge immediately after collection. Observe the same precautions in the preparation of serum samples.
6. If heparin is to be used as anti-coagulant, the effect on the assay outcome at the dose of heparin used should be pre-determined.
7. Avoid using samples with gross hemolysis or lipemia.

## Rat Leptin ELISA Assay Procedure

Warm all reagents to room temperature before setting up the assay.

1. Dilute the 10X concentrated HRP Wash Buffer 10-fold by mixing the entire contents of both buffer bottles with 900 mL de-ionized or distilled water.
2. Remove the required number of strips from the Microtiter Assay Plate. Unused strips should be resealed in the foil pouch and stored at 2-8 °C. Assemble strips in an empty plate holder and wash each well 3 times with 300  $\mu$ L of diluted Wash Buffer per wash. Decant Wash Buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times. **Do not let wells dry before proceeding to the next step. If automated machine is used for assay, follow the manufacturer's instructions for all washing steps described in this protocol.**
3. Add 30  $\mu$ L Assay Buffer to Background wells, Standard wells, and QC1 and QC2 wells. Add 40  $\mu$ L Assay Buffer to sample wells.
4. If samples to be assayed are serum or plasma, add 10  $\mu$ L Matrix Solution to the Background wells, Standard wells, and QC1 and QC2 wells. If samples are free of significant serum matrix components, add 10  $\mu$ L Assay Buffer instead.
5. Add 10  $\mu$ L Assay Buffer to the Background wells and add in duplicates 10  $\mu$ L Rat Leptin Standards in the order of ascending concentration to the appropriate wells.
6. Add 10  $\mu$ L QC1 and 10  $\mu$ L QC2 to the appropriate wells.
7. Add sequentially 10  $\mu$ L of the unknown samples in duplicate to the remaining wells.
8. Transfer Antiserum Solution to a reagent reservoir and add 50  $\mu$ L of this solution to each well with a multi-channel pipette. Cover the plate with plate sealer and incubate at room temperature for 2 hours on an orbital microtiter plate shaker set to rotate at moderate speed, about 400 to 500 rpm.
9. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in well.
10. Wash wells 3 times with diluted Wash Buffer, 300  $\mu$ l per well per wash. Decant and tap after each wash to remove residual buffer.
11. Add 100  $\mu$ L Detection Antibody to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 1 hour on an orbital microtiter plate shaker set to rotate at moderate speed, approximately 400-500 rpm.
12. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in well.
13. Wash wells 3 times with diluted Wash Buffer, 300  $\mu$ l per well per wash. Decant and tap after each wash to remove residual buffer.



- 
14. Add 100  $\mu$ l Enzyme Solution to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 30 minutes on the micro-titer plate shaker.
  15. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in well.
  16. Wash wells 6 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap after each wash to remove residual buffer.
  17. Add 100  $\mu$ l of Substrate solution to each well, cover plate with sealer and shake in the plate shaker for approximately 5 to 20 minutes. Blue color should be formed in wells of Leptin standards with intensity proportional to increasing concentrations of Leptin.

**Note:** Please be aware that the color may develop more quickly or more slowly than the recommended incubation time depending on the localized room temperature. Please visually monitor the color development to optimize the incubation time. One can monitor color development using 370 nm filter, if available on the spectrophotometer. When the absorbance is between 1.2 and 1.8 at 370 nm, the stop solution can be added to terminate the color development.

18. Remove sealer and add 100  $\mu$ L Stop Solution (**Caution:** Corrosive Solution) and shake plate by hand to ensure complete mixing of solution in all wells. The blue color should turn into yellow after acidification. Wipe the bottom of the microtiter plate to remove any residue prior to reading on plate reader. Read absorbance at 450 nm and 590 nm in a plate reader within 5 minutes and ensure that there are no air bubbles in any well. Record the difference of absorbance units.

## Assay Procedure for Rat Leptin ELISA Kit

	Step 1	Step 2	Step 3	Step 4	Step 5-7	Step 8	Step 9-10	Step 11	Step 12-13	Step 14	Step 15-16	Step 17		Step 18	
Well#	Dilute both bottles of 10X Wash Buffer with 900 mL Deionized Water	Wash plate 3X with 300 µL Wash Buffer. Remove residual buffer by tapping smartly on absorbent towels	Assay Buffer	Matrix	StdS/QCs/ Samples	Anti serum	Seal, Agitate, Incubate 2 hours at Room Temperature. Wash 3X with 300 µL Wash Buffer.	Detect. Anti.	Seal, Agitate, Incubate 1 hour at Room Temperature. Remove residual buffer by tapping smartly on absorbent towels. Wash 3X with 300 µL Wash Buffer.	Enzyme Solution	Seal, Agitate, Incubate 30 minutes at Room Temperature. Wash 6X with 300 µL Wash Buffer.	Substrate	Seal, Agitate, Incubate 10-15 minutes at Room Temperature.	Stop Solution	Shake by hand. Read Absorbance at 450 nm and 590 nm within 5 mins
A1 B1			30 µL	10 µL	-	50 µL		100 µL		100 µL		100 µL		100 µL	
C1 D1			30 µL	10 µL	10 µL of 0.2 ng/mL Std.										
E1 F1			30 µL	10 µL	10 µL of 0.5 ng/mL Std.										
G1 H1			30 µL	10 µL	10 µL of 1 ng/mL Std.										
A2 B2			30 µL	10 µL	10 µL of 2 ng/mL Std.										
C2 D2			30 µL	10 µL	10 µL of 5 ng/mL Std.										
E2 F2			30 µL	10 µL	10 µL of 10 ng/mL Std.										
G2 H2			30 µL	10 µL	10 µL of 20 ng/mL Std.										
A3 B3			30 µL	10 µL	10 µL of 30 ng/mL Std.										
C3 D3			30 µL	10 µL	10 µL of QC2										
E3 F3			30 µL	10 µL	10 µL of QC1										
G3 H3			40 µL	-	10 µL of Sample										
A4 B4			40 µL	-	10 µL of Sample										

For research use only. Not for use in diagnostic procedures.

## Microtiter Plate Arrangement

### Rat Leptin ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	2.0 ng/mL	30 ng/mL	Sample 2								
B	Blank	2.0 ng/mL	30 ng/mL	Sample 2								
C	0.2 ng/mL	5.0 ng/mL	QC 1	Etc.								
D	0.2 ng/mL	5.0 ng/mL	QC 1									
E	0.5 ng/mL	10 ng/mL	QC 2									
F	0.5 ng/mL	10 ng/mL	QC 2									
G	1.0 ng/mL	20 ng/mL	Sample 1									
H	1.0 ng/mL	20 ng/mL	Sample 1									

For research use only. Not for use in diagnostic procedures.

---

## Calculations

Graph a reference curve by plotting the absorbance unit of 450nm, less unit at 590nm, on the Y-axis against the concentrations of rat leptin standard on the X-axis. The dose-response curve of this assay fits best to a sigmoidal 4- or 5-parameter logistic equation. The results of unknown samples can be calculated with any computer program having a 4- or 5-parameter logistic function.

**Note:** When sample volumes assayed differ from 10  $\mu\text{L}$ , an appropriate mathematical adjustment must be made to accommodate for the dilution factor (for example, if 5  $\mu\text{L}$  of sample is used, then calculated data must be multiplied by 2). When sample volume assayed is less than 10  $\mu\text{L}$ , compensate the volume deficit with either matrix solution or assay buffer, whichever is appropriate.

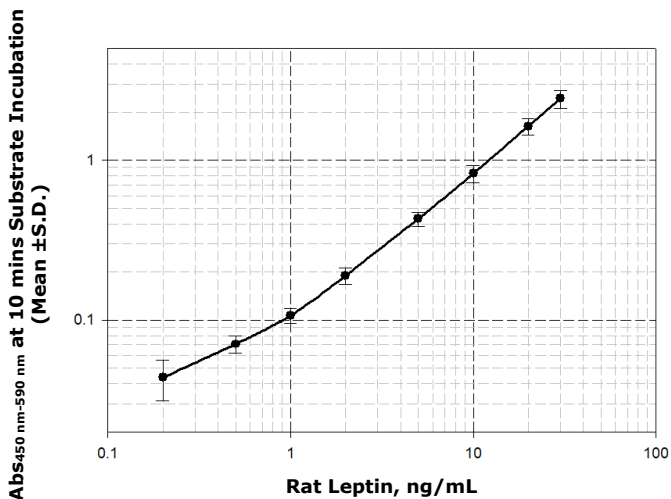
## Interpretation

- The assay will be considered accepted when all Quality Control values fall within the calculated Quality Control Range; if any QC's fall outside the control range, review results with the supervisor.
- If the difference between duplicate results of a sample is  $>15\%$  CV, repeat the sample.
- The limit of sensitivity of this assay is 0.08 ng/mL ( $\sim 5$  pM) leptin (10  $\mu\text{L}$  sample size).
- The appropriate range of this assay is 0.2 ng/mL to 30 ng/mL leptin (10  $\mu\text{L}$  sample size). Any result greater than 30 ng/mL in a 10  $\mu\text{L}$  sample assayed should be repeated on dilution using either matrix solution or assay buffer, whichever is appropriate, as diluent until it falls within range.

## Graph of Typical Reference Curve

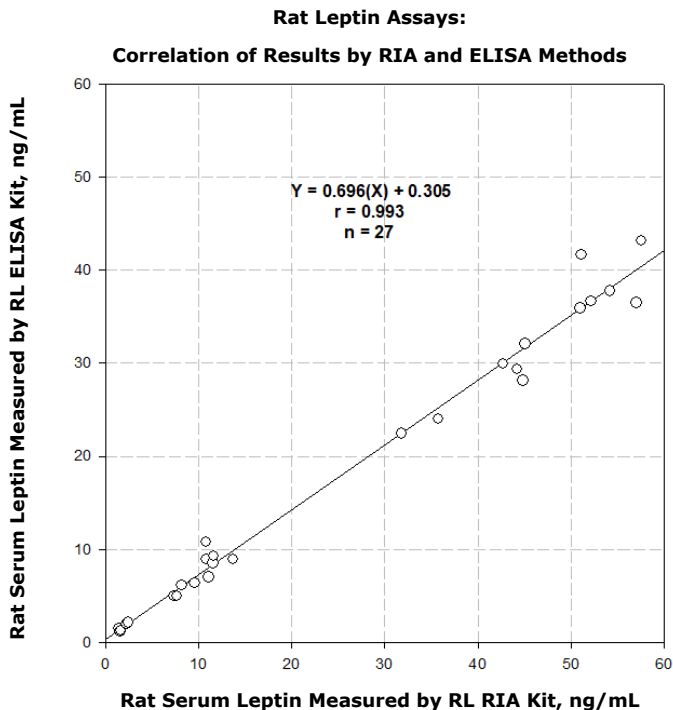
### Rat Leptin ELISA

(n = 9 assays)



For Demonstration Only—Do not use for calculations

## Correlation Graph



Serum samples from 27 rats were assayed for leptin using both Rat Leptin RIA Kit (RL-83K) and Rat Leptin ELISA Kit (EZRL-83K). Correlation of the two kits are derived by linear regression analysis of paired results from each sample.

## Assay Characteristics

### Sensitivity

The lowest level of rat leptin that can be detected by this assay is 0.08 ng/mL using a 10  $\mu$ L sample size.

### Specificity

The specificity (also known as selectivity) of the analytical test is its ability to selectively measure the analytes in the presence of other like components in the sample matrix.

Rat Leptin	100%
Mouse Leptin	143%
Human Leptin	15%
Porcine Leptin	< 0.1%
Ovine Leptin	< 0.1%
Chicken Leptin	< 0.1%
Rat Insulin	0%
Rat C-Peptide	0%
Human Proinsulin	0%
Porcine Proinsulin	0%
Bovine Proinsulin	0%
Glucagon	0%
Human Ghrelin	0%

### Precision

Sample Number	Mean Leptin Level (ng/mL)	Assay Variation (CV)	
		Intra-assay	Inter-assay
1	1.26	2.49%	3.93%
2	5.19	1.88%	3.31%
3	16.56	2.13%	2.95%

The assay variations of Rat Leptin ELISA kit were studied on three rat serum samples with varying concentrations of spiked analyte. The intra-assay variations are calculated from eight duplicate determinations in an assay. The inter-assay variations are calculated from results of 6 separate assays with duplicate samples in each assay.

## Spike and Recovery of Rat Leptin in Rat Serum

Serum Sample #	Rat Leptin		Recovery (%) of Spiked Insulin
	Added (ng/mL)	Observed (ng/mL)	
<b>Rat Serum #1</b>	0	5.00	-
	0.5	5.61	122
	2.0	7.29	115
	10.0	16.17	112
<b>Rat Serum #2</b>	0	5.85	-
	0.5	6.36	102
	2.0	8.20	118
	10.0	16.93	111

Rat leptin at indicated levels was added to two rat serum samples and the resulting leptin content of each sample was assayed by ELISA. The % of recovery = [(observed leptin level after spike - observed leptin level before spike) / spiked level of leptin] x 100%. Mean recovery rate at spiked leptin level of 0.5, 2, and 10 ng/mL is 112%, 117%, and 112%, respectively.



## Linearity

### Effect of Serum Dilution

Serum Sample #	Dilution Factor	Leptin Level		
		Observed (ng/mL)	Expected (ng/mL)	% Of Expected
Rat Serum #1	-	21.18		100
	2x	19.56		92
	5x	19.20	21.18	91
	10x	18.40		87
	20x	19.20		91
Rat Serum #2	-	21.39		100
	2x	20.92		98
	5x	21.00	21.39	98
	10x	21.10		99
	20x	20.80		97

Two rat serum samples are diluted each with matrix solution to various degrees as indicated and assayed for leptin levels along with neat samples of each serum. Measured leptin levels are corrected for dilution factors and reported as observed leptin level.

---

## Quality Controls

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert, or available at our website [SigmaAldrich.com](http://SigmaAldrich.com).

## Troubleshooting

- To obtain reliable and reproducible results the operator should carefully read this manual and fully understand all aspects of each assay step before attempting to run the assay.
- Throughout the assay the operator should adhere strictly to the procedures with good laboratory practice.
- Have all necessary reagents and equipment ready on hand before starting. Once the assay has been started all steps should be completed with precise timing and without interruption.
- Avoid cross contamination of any reagents or samples to be used in the assay.
- Make sure all reagents and samples are added to the bottom of each well.
- Careful and complete mixing of solutions in the well is critical. Poor assay precision will result from incomplete mixing or cross well contamination due to inappropriate mixing.
- Remove any air bubbles formed in the well after acidification of substrate solution because bubbles interfere with spectrophotometric readings.
- Do not let the absorbance reading of the highest standard fall beyond the limit of your microtiter plate reader's capacity. Adjust the length of substrate incubation time accordingly.
- High absorbance in the background or blank wells could be due to:
  - Cross-well contamination by standard solution or sample and
  - Inadequate washing of wells with wash buffer.

## Product Ordering

Products are available for online ordering at [SigmaAldrich.com](https://www.sigmaaldrich.com).

### Replacement Reagents

<b>Reagents</b>	<b>Cat. No.</b>
Rat/Mouse Leptin ELISA Plates	EP83
Rat/Mouse Leptin Antiserum	EAS83
10X HRP Wash Buffer Concentrate	EWB-HRP
Rat Leptin Standards	E8083-K
Rat Leptin Quality Controls 1 and 2	E6083-K
Rat/Mouse Leptin Matrix Solution	EPS0016
Assay Buffer	EAB-PTR
Rat/Mouse Leptin Detection Antibody	E1083
Enzyme Solution	EHRP-4
Substrate	ESS-TMB2
Stop Solution	ET-TMB
10-pack of Rat Leptin ELISA Kits	EZRL-83BK

---

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma, Millipore, Milliplex and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2007-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

For research use only. Not for use in diagnostic procedures.