



**Porcine Insulin**

**250 Tubes**

**Cat. # PI-12K**

**PORCINE INSULIN RIA KIT  
250 TUBES (Cat. # PI-12K)**

I.	Intended Use	2
II.	Principles Of Procedure	2
III.	Reagents Supplied	3
IV.	Storage and Stability	3
V.	Reagent Precautions	4
VI.	Materials Required But Not Provided	5
VII.	Specimen Collection And Storage	5
VIII.	Assay Procedure	5
IX.	Calculations and Transformations	8
X.	Interpretation	8
XI.	Normal Fasting Range	8
XII.	Assay Characteristics	9
XIII.	Quality Controls	11
XIV.	Replacement Reagents	11
XV.	Ordering Information	12
XVI.	References	12

**PORCINE INSULIN RIA KIT  
250 TUBES (Cat. # PI-12K)**

**I. INTENDED USE**

EMD Millipore's Porcine Insulin Radioimmunoassay (RIA) Kit has been developed for the measurement of Porcine Insulin in plasma, serum, or tissue culture media. The tracer is prepared with human insulin. ***This kit is for Research Use Only. Not for Use in Diagnostic Procedures.***

**II. PRINCIPLES OF PROCEDURE**

In radioimmunoassay, a fixed concentration of labeled tracer antigen is incubated with a constant dilution of antiserum such that the concentration of antigen binding sites on the antibody is limited, for example, only 50% of the total tracer concentration may be bound by antibody. If unlabeled antigen is added to this system, there is competition between labeled tracer and unlabeled antigen for the limited and constant number of binding sites on the antibody. Thus, the amount of tracer bound to antibody will decrease as the concentration of unlabeled antigen increases. This can be measured after separating antibody-bound from free tracer and counting one or the other, or both fractions. A calibration or standard curve is set up with increasing concentrations of standard unlabeled antigen and from this curve the amount of antigen in unknown samples can be calculated. Thus, the four basic necessities for a radioimmunoassay system are: a specific antiserum to the antigen to be measured, the availability of a radioactive labeled form of the antigen, a method whereby antibody-bound tracer can be separated from the unbound tracer, and finally, an instrument to count radioactivity.

The EMD Millipore Porcine Insulin assay utilizes <sup>125</sup>I-labeled Insulin and an Insulin antiserum to determine the level of Insulin in serum, plasma or tissue culture media by the double antibody/PEG technique.

### III. REAGENTS SUPPLIED

Each kit is sufficient to run 250 tubes and contains the following reagents.

#### A. Assay Buffer

0.05M Phosphosaline pH 7.4 containing 0.025M EDTA, 0.08% Sodium Azide, and 1% RIA Grade BSA

Quantity: 40 mL/vial

Preparation: Ready to use

#### B. Insulin Antibody

Insulin Serum in Assay Buffer

Quantity: 26 mL/vial

Preparation: Ready to use

#### C. <sup>125</sup>I-Insulin

<sup>125</sup>I-Insulin Label, HPLC purified (specific activity 367  $\mu$ Ci/ $\mu$ g)

Lyophilized for stability. Freshly iodinated label contains <5  $\mu$ Ci, <185 kBq calibrated to the 1st Monday of each month.

Quantity: 27 mL/vial upon hydration

Preparation: Contents Lyophilized. Hydrate with entire contents of Label Hydrating Buffer. Allow to sit at room temperature for 30 minutes, with occasional gentle mixing.

#### D. Label Hydrating Buffer

Assay Buffer containing Normal Guinea Pig IgG as a carrier. Used to hydrate <sup>125</sup>I-Insulin

Quantity: 27 mL/vial

Preparation: Ready to use

#### E. Human Insulin Standards

Purified Recombinant Human Insulin in Assay Buffer at the following concentration: 200  $\mu$ U/mL

Quantity: 2 mL/vial

Preparation: Ready to use

#### F. Quality Controls 1 & 2

Purified Recombinant Porcine Insulin in Assay Buffer

Quantity: 1 mL/vial

Preparation: Ready to use

#### G. Precipitating Reagent

Goat anti Guinea Pig IgG Serum, 3% PEG and 0.05% Triton X-100 in 0.05M Phosphosaline, 0.025M EDTA, 0.08% Sodium Azide

Quantity: 260 mL/vial

Preparation: Ready to use; chill to 4°C.

### IV. STORAGE AND STABILITY

Refrigerate all reagents between 2 and 8°C for short term storage. For prolonged storage (>2 weeks), freeze at  $\leq -20^\circ\text{C}$ . Avoid multiple (>5) freeze/thaw cycles. Refer to date on bottle for expiration when stored at  $\leq -20^\circ\text{C}$ . Do not mix reagents from different kits unless they have the same lot number.

## V. REAGENT PRECAUTIONS

### A. Radioactive Materials

This radioactive material may be received, acquired, possessed and used only by research personnel or clinical laboratories for in vitro research tests not involving internal or external administration of the material, or the radiation there from, to human beings or animals. Its receipt, acquisition, possession, use and transfer are subject to the regulations of the U. S. Nuclear Regulatory Commission (NRC) or of a State with which the Commission has entered into an agreement for the exercise of regulatory authority.




The following are suggested general rules for the safe use of radioactive material. The customer's Radiation Safety Officer is ultimately responsible for the safe handling and use of radioactive material.

1. Wear appropriate personal devices at all times while in areas where radioactive materials are used or stored.
2. Wear laboratory coats, disposable gloves, and other protective clothing at all times.
3. Monitor hands, shoes, and clothing and immediate area surrounding the work station for contamination after each procedure and before leaving the area.
4. Do not eat, drink, or smoke in any area where radioactive materials are stored or used.
5. Never pipette radioactive material by mouth.
6. Dispose of radioactive waste in accordance with NRC rules and regulations.
7. Avoid contaminating objects such as telephones, light switches, doorknobs, etc.
8. Use absorbent pads for containing and easily disposing of small amounts of contamination.
9. Wipe up all spills immediately and thoroughly and dispose of the contaminated materials as radioactive waste. Inform Radiation Safety Officer.

### B. Sodium Azide

Sodium Azide has been added to all reagents as a preservative at a concentration of 0.08%. Although it is at a minimum concentration, sodium azide may react with lead and copper plumbing to form highly explosive metal azides. On disposal, flush with a large volume of water to prevent azide build up.

#### Full labels of hazardous components in this kit:

Ingredient, Cat #		Full Label	
<sup>125</sup> I-Insulin Tracer	9011		<p><b>Danger.</b> Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/ attention.</p>
Insulin Antibody	1013-K		<p><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.</p>
Precipitating Reagent	PR-UV		<p><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.</p>

## VI. MATERIALS REQUIRED BUT NOT PROVIDED

1. Borosilicate glass tubes, 12 x 75 mm. (NOTE: Polypropylene or polystyrene tubes may be used if the investigator finds that the pellet formation is acceptably stable in their system.)
2. 100  $\mu$ L pipet with disposable tips
3. 100  $\mu$ L & 1.0 mL repeating dispenser
4. Refrigerated swing bucket centrifuge capable of developing 2,000 - 3,000xg. (Use of fixed-angle buckets is not recommended.)
5. Absorbent paper
6. Vortex mixer
7. Refrigerator
8. Gamma Counter

## VII. SPECIMEN COLLECTION AND STORAGE

1. A maximum of 100  $\mu$ L per assay tube of serum or plasma can be used, although, 50  $\mu$ L per assay tube is adequate for most applications. Tissue culture and other media may also be used.
2. Care must be taken when using heparin as an anticoagulant, since an excess will provide falsely high values.<sup>2</sup> Use no more than 10 IU heparin per mL of blood collected.
3. Specimens can be stored at 4°C if they will be tested within 24 hours of collection. For longer storage, specimens should be stored at  $\leq$  -20°C. Avoid multiple (>5) freeze/thaw cycles.
4. Avoid using samples with gross hemolysis or lipemia.

## VIII. ASSAY PROCEDURE

### Standard Preparation

Use care in opening the Standard vial.

Label six glass tubes 1, 2, 3, 4, 5, and 6. Add 1.0 mL Assay Buffer to each of the six tubes. Prepare serial dilutions by adding 1.0 mL of the 200uU/mL standard to tube 1, mix well and transfer 1.0 mL of tube 1 to tube 2, mix well and transfer 1.0 mL of tube 2 to tube 3, mix well and transfer 1.0 mL of tube 3 to tube 4, mix well and transfer 1.0 mL of tube 4 to tube 5, mix well and transfer 1.0 mL of tube 5 to tube 6, mix well.

Note: Do not use a Repeater pipette. Change tip for every dilution. Wet tip with Standard before dispensing. Unused portions of standard should be stored at  $\leq$  -20°C. Avoid multiple freeze/thaw cycles.

Tube #	Standard Concentration uU/mL	Volume of Assay Buffer to Add	Volume of Standard to Add
1	100 uU/mL	1.0 mL	1.0 mL of 200 uU/mL
2	50 uU/mL	1.0 mL	1.0 mL of 100 uU/mL
3	25 uU/mL	1.0 mL	1.0 mL of 50 uU/mL
4	12.5 uU/mL	1.0 mL	1.0 mL of 25 uU/mL
5	6.25 uU/mL	1.0 mL	1.0 mL of 12.5 uU/mL
6	3.125 uU/mL	1.0 mL	1.0 mL of 6.25 uU/mL

For optimal results, accurate pipetting and adherence to the protocol are recommended.

## VIII. ASSAY PROCEDURE (continued)

### A. Assay Set-Up, Day One

1. Pipet 300  $\mu$ L of Assay Buffer to the Non-Specific Binding (NSB) tubes (3-4), 200  $\mu$ L to Reference (Bo) tubes (5-6), and 100  $\mu$ L to tubes 7 through the end of the assay.
2. Pipet 100  $\mu$ L of Standards and Quality Controls in duplicate (see flow chart).
3. Pipet 100  $\mu$ L of each sample in duplicate. (NOTE: Smaller volumes of sample may be used when Insulin concentrations are anticipated to be elevated or when sample size is limited. Additional Assay Buffer should be added to compensate for the difference so that the volume is equivalent to 100  $\mu$ L (e.g., when using 50  $\mu$ L of sample, add 50  $\mu$ L of Assay Buffer). Refer to Section IX for calculation modification.
4. Pipet 100  $\mu$ L of  $^{125}$ I-Insulin to all tubes. Important: For preparation, see Section III, Part C.
5. Pipet 100  $\mu$ L of Insulin antibody to all tubes except Total Count tubes (1-2) and NSB tubes (3-4).
6. Vortex, cover, and incubate overnight (20-24 hours) at 4°C.

### B. Day Two

7. Add 1.0 mL of cold (4°C) Precipitating Reagent to all tubes (except Total Count tubes).
8. Vortex and incubate 20 minutes at 4°C.
9. Centrifuge, 4°C, all tubes [except Total Count tubes (1-2)] for 20 minutes at 2,000-3,000 xg. NOTE: If less than 2,000 xg is used or if slipped pellets have been observed in previous runs, the time of centrifugation must be increased to obtain a firm pellet (e.g., 40 minutes). Multiple centrifuge runs within an assay must be consistent.  
$$xg = (1.12 \times 10^{-5}) (r) (rpm)^2$$

r = radial distance in cm (from axis of rotation to the bottom of the tube)  
rpm = rotational velocity of the rotor
10. Immediately decant the supernate of all tubes except Total Count tubes (1-2), drain tubes for at least 15-60 seconds (be consistent between racks), and blot excess liquid from lip of tubes. NOTE: Invert tubes only one time. Pellets are fragile and slipping may occur.
11. Count all tubes in a gamma counter for 1 minute. Calculate the  $\mu$ U/mL of Porcine Insulin in unknown samples using automated data reduction procedures.

VIII. ASSAY PROCEDURE (continued)

Assay Flow Chart

Day One						Day Two		
	Step 1	Step 2-3	Step 4	Step 5	Step 6	Step 7	Step 8	Step 9-11
Tube #	Add Assay Buffer	Add Standard / QC Sample	Add <sup>125</sup> I-Insulin Tracer	Add Insulin Antibody	<b>Vortex, Cover, and Incubate 20-24 hrs at 4°C</b>	Add Precipitating Reagent	<b>Vortex, and Incubate 20 min. at 4°C</b>	<b>Centrifuge at 4°C for 20 min., Decant, and Count pellets</b>
1,2	----	----	100 µL	----		----		
3,4	300 µL	----	100 µL	----		1.0 mL		
5,6	200 µL	----	100 µL	100 µL		1.0 mL		
7,8	100 µL	100 µL of 3.125 µU/mL	100 µL	100 µL		1.0 mL		
9,10	100 µL	100 µL of 6.25 µU/mL	100 µL	100 µL		1.0 mL		
11,12	100 µL	100 µL of 12.5 µU/mL	100 µL	100 µL		1.0 mL		
13,14	100 µL	100 µL of 25 µU/mL	100 µL	100 µL		1.0 mL		
15,16	100 µL	100 µL of 50 µU/mL	100 µL	100 µL		1.0 mL		
17,18	100 µL	100 µL of 100 µU/mL	100 µL	100 µL		1.0 mL		
19,20	100 µL	100 µL of 200 µU/mL	100 µL	100 µL		1.0 mL		
21,22	100 µL	100 µL of QC 1	100 µL	100 µL		1.0 mL		
23,24	100 µL	100 µL of QC 2	100 µL	100 µL		1.0 mL		
25-n	100 µL	100 µL of unknown	100 µL	100 µL	1.0 mL			



## IX. CALCULATIONS AND TRANSFORMATIONS

### A. Explanation

The calculations for Porcine Insulin can be automatically performed by most gamma counters possessing data reduction capabilities or by independent treatment of the raw data using a commercially available software package. Choose weighted 4-parameter or weighted log/logit for the mathematical treatment of the data. [NOTE: Be certain the procedure used subtracts the NSB counts from each average count, except Total Counts, prior to final data reduction.]

### B. Manual Calculation

1. Average duplicate counts for Total Count tubes (1-2), NSB tubes (3-4), Total Binding tubes (reference, Bo) (5-6), and all duplicate tubes for standards and samples to the end of the assay.
2. Subtract the average NSB counts from each average count (except for Total Counts). These counts are used in the following calculations.
3. Calculate the percentage of tracer bound  
(Total Binding Counts/Total Counts) X 100  
This should be 35-50%.
4. Calculate the percentage of total binding (%B/Bo) for each standard and sample.  
 $\%B/Bo = (\text{Sample or Standard}/\text{Total Binding}) \times 100$
5. Plot the % B/Bo for each standard on the y-axis and the known concentration of the standard on the x-axis using log-log graph paper.
6. Construct the reference curve by joining the points with a smooth curve.
7. Determine the  $\mu\text{U}/\text{mL}$  of Porcine Insulin in the unknown samples and controls by interpolation of the reference curve.

[NOTE: When sample volumes assayed differ from 100  $\mu\text{L}$ , an appropriate mathematical adjustment must be made to accommodate for the dilution factor (e.g., if 50  $\mu\text{L}$  of sample is used, then calculated data must be multiplied by 2).]

## X. INTERPRETATION

### A. Acceptance Criteria

1. The run 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.
2. If the difference between duplicate results of a sample is >10% CV, repeat the sample.
3. The limit of sensitivity for the Porcine Insulin assay is 1.611  $\mu\text{U}/\text{mL}$  (100  $\mu\text{L}$  sample size).
4. The limit of linearity for the Porcine Insulin assay is 200  $\mu\text{U}/\text{mL}$  (100  $\mu\text{L}$  sample size). Any result greater than 200  $\mu\text{U}/\text{mL}$  should be repeated on dilution using Assay Buffer as a diluent.

## XI. NORMAL FASTING RANGE

5-15  $\mu\text{U}/\text{mL}$

## XII. ASSAY CHARACTERISTICS

### A. Sensitivity

The lowest level of insulin that can be detected by this assay is 1.611  $\mu\text{U/mL}$  when using a 100  $\mu\text{l}$  sample size.

### B. Performance

The following parameters of assay performance are expressed as Mean  $\pm$  Standard Deviation.

$$\text{ED}_{80} = 5 \pm 2 \mu\text{U/mL}$$

$$\text{ED}_{50} = 22 \pm 6 \mu\text{U/mL}$$

$$\text{ED}_{20} = 98 \pm 24 \mu\text{U/mL}$$

### C. Specificity

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

Porcine Insulin	100%
Human Insulin	100%
Bovine Insulin	90%
Intact Human Proinsulin	38%
Des 31,32 HPI	47%
Des 64,65 HPI	72%
Glucagon	*
Somatostatin	*
Pancreatic Polypeptide	*

\* Not Detectable

## XII. ASSAY CHARACTERISTICS (continued)

### D. Example of Assay Results

This data is presented as an example only and should not be used in lieu of a standard curve prepared with each assay.

Tube #	ID	CPM	Ave CPM	Ave Net	%	μU/mL
				CPM	B/Bo	
1	Totals	16371				
2	"	16333	16352			
3	NSB	569				
4	"	570	570			
5	Bo	7537				
6	"	7638	7588	7018		
<u>Standards</u>						
7	3.125 μU/mL	6741				
8		6646	6694	6124	87.3	
9	6.25 μU/mL	6037				
10		6117	6077	5508	78.5	
11	12.5 μU/mL	5156				
12		4911	5034	4464	63.6	
13	25 μU/mL	3993				
14		3728	3861	3291	46.9	
15	50 μU/mL	2727				
16		2721	2724	2155	30.7	
17	100 μU/mL	1866				
18		1860	1863	1294	18.4	
19	200 μU/mL	1345				
20		1229	1287	718	10.2	
<u>Controls/Unknown</u>						
21	QC 1	4903				
22		4947	4925	4356	62.1	13.48
23	QC 2	2751				
24		2754	2753	2183	31.1	49.39
25-n	Unknown					

### XIII. QUALITY CONTROLS

Good Laboratory Practice (GLP) requires that Quality Control (QC) specimens be run with each standard curve to check the assay performance. Two levels of controls are provided for this purpose. These and any other control materials should be assayed repeatedly to establish mean values and acceptable ranges. Each individual laboratory is responsible for defining their system for quality control decisions and is also responsible for making this system a written part of their laboratory manual. The ranges for Quality Control 1 and 2 are provided on the card insert or can be located at the EMD Millipore website [emdmillipore.com](http://emdmillipore.com) using the catalog number as the keyword.

Recommended batch analysis decision using two controls (Westgard Rules):<sup>4</sup>

1. When both controls are within  $\pm 2$  SD. Decision: Approve batch and release analyte results.
2. When one control is outside  $\pm 2$  SD and the second control is within  $\pm 2$  SD. Decision: Hold results, check with supervisor. If no obvious source of error is identified by the below mentioned check of systems, the supervisor may decide to release the results.

Technician check of systems:

1. Check for calculation errors
2. Repeat standards and controls
3. Check reagent solutions
4. Check instrument

### XIV. REPLACEMENT REAGENTS

Reagent	Cat #
<sup>125</sup> I-Insulin (<5 $\mu$ Ci, <185 kBq)	9011
Label Hydrating Buffer (27 mL)	LHB-P
Human Insulin Standards (2 mL each)	8014-K
Insulin Antibody (26 mL)	1013-K
Precipitating Reagent (260 mL)	PR-UV
Quality Control 1 & 2 (1 mL each)	6000-K
Assay Buffer (40 mL)	AB-P

## XV. ORDERING INFORMATION

To place an order or to obtain additional information about our immunoassay products, please contact your Customer Service or Technical Support Specialist.

Contact information for each region can be found on our website:

[emdmillipore.com/contact](http://emdmillipore.com/contact)

### Conditions of Sale

For Research Use Only. Not for Use in Diagnostic Procedures.

### Safety Data Sheets (SDS)

Safety Data Sheets for EMD Millipore products may be ordered by fax or phone or through our website at [emdmillipore.com/msds](http://emdmillipore.com/msds).

## XVI. REFERENCES

1. Morgan, C.R., and Lazarow, A. Immunoassay of insulin: Two antibody system. Plasma insulin levels in normal, subdiabetic and diabetic rats. *Diabetes*. 12: 115-126, 1963.
2. Thorell, J.I. *Scand. J. Clin. Lab. Invest.* 31:187, 1973.
3. Feldman, H. and Rodbard, D. "Mathematical Theory of Radioimmunoassay," in: W.D Odell and Doughaday, W.H. (Ed.), Principles of Competitive Protein-Binding Assays. Philadelphia: J.B. Leppincott Company; pp 158-203, 1971.
4. Westgard, J.O., et. al. A multi-rule Shewhart chart for quality control in clinical chemistry. *Clin. Chem.* 27:493-501, 1981.