



Canine C-Peptide

125 Tubes

Cat. # CCP-24HK

CANINE C-PEPTIDE RIA KIT
125 TUBES (Cat. # CCP-24HK)

| | | |
|-------|-------------------------------------|----|
| I. | Intended Use | 2 |
| II. | Principles Of Procedure | 2 |
| III. | Reagents Supplied | 2 |
| IV. | Storage and Stability | 3 |
| V. | Reagent Precautions | 3 |
| VI. | Materials Required But Not Provided | 5 |
| VII. | Specimen Collection And Storage | 5 |
| VIII. | Assay Procedure | 6 |
| IX. | Calculation | 9 |
| X. | Interpretation | 9 |
| XI. | Assay Characteristics | 10 |
| XII. | Normal Fasting Range | 12 |
| XIII. | Quality Controls | 13 |
| XIV. | Replacement Reagents | 13 |
| XV. | Ordering Information | 14 |
| XVI. | References | 14 |

CANINE C-PEPTIDE RIA KIT
125 TUBES (Cat. # CCP-24HK)

I. INTENDED USE

EMD Millipore's Canine C-Peptide Radioimmunoassay kit is for the quantitative determination of Canine C-Peptide in serum, plasma, and other biological media. It is a completely homologous assay since the antibody was raised against purified Canine C-Peptide and both the tracer and the standard are prepared with Canine C-Peptide.

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 Canine C-Peptide assay utilizes ¹²⁵I-labeled Canine C-Peptide and a Canine C-Peptide antiserum to determine the level of C-Peptide in serum, plasma or tissue culture media by the double antibody/PEG technique.

III. REAGENTS SUPPLIED

Each kit is sufficient to run 125 tubes and contains the following reagents:

A. Assay Buffer

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

Quantity: 40 mL/vial.

Preparation: Ready to use

B. Canine C-Peptide Antibody

Guinea Pig anti-Canine C-Peptide Antibody in Assay Buffer

Quantity: 13 mL/vial

Preparation: Ready to use

C. ¹²⁵I-Canine C-Peptide

¹²⁵I-Canine C-Peptide Label, HPLC purified (specific activity 628 μCi/μg)

Lyophilized for stability. Freshly iodinated label contains 1.5 μCi (56 kBq), calibrated to the 1st Monday of each month.

Quantity: 13.5 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.

III. REAGENTS SUPPLIED (continued)

D. Label Hydrating Buffer

Assay Buffer containing Normal Guinea Pig Serum as a carrier. Used to hydrate ¹²⁵I-Canine C-Peptide.

Quantity: 13.5 mL/vial

Preparation: Ready to use

E. Canine C-Peptide Calibrator

Purified Canine C-Peptide in Assay Buffer

Lyophilized for stability

Quantity: 2 mL/vial upon hydration

Preparation: Hydrate with 2 mL deionized/distilled water.

Note: The actual calibrator concentration of Canine C-Peptide present in the vial will be lot dependent.

Please refer to the Canine C-Peptide analysis sheet for exact Canine C-Peptide calibrator concentration present in a specific lot.

F. Quality Controls 1 & 2

Purified Canine C-Peptide in Assay Buffer

Quantity: Lyophilized, 1 mL/vial upon hydration

Preparation: Hydrate with 1 mL deionized/distilled water.

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: 130 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 $\leq -20^{\circ}\text{C}$. Reconstituted Canine C-Peptide Calibrator should be stored 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.

Note: See Full Labels of Hazardous components on next page.

V. REAGENT PRECAUTIONS (continued)

Full labels of hazardous components in this kit:

| Ingredient, Cat # | | Full Label | |
|--|---------|--|--|
| Quality Controls 1 and 2 (lyophilized) | 6000L-K |  | Warning. Harmful if swallowed. Harmful to aquatic life with long lasting effects. Avoid release to the environment. |
| Precipitating Reagent | PR-UVHK |  | Warning. 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. |
| Canine C-Peptide Antibody | 1024-HK |  | Warning. 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. |
| Canine C-Peptide Standard | 8024-K |   | Danger. 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. |
| ¹²⁵ I-Canine C-Peptide Tracer | 9024-HK |   | Danger. 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. |

V. REAGENT PRECAUTIONS (continued)

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.

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 μ L & 1.0 mL pipet with disposable tips
3. 100 μ L & 1.0 mL repeating dispenser
4. Refrigerated swing bucket centrifuge capable of developing 2,000 - 3,000 xg. (Use of fixed-angle buckets are not recommended.)
5. Absorbent paper
6. Vortex mixer
7. Refrigerator
8. Gamma Counter

VII. SPECIMEN COLLECTION AND STORAGE

1. A maximum of 100 μ L per assay tube of serum or plasma can be used, although, 50 μ 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 values². Use no more than 10 IU heparin per mL of blood collected.
3. Canine C-Peptide must be protected from proteolysis during assay procedures and sample storage. Trasylol (Aprotinin) at a concentration of 500 KIU per mL of serum or plasma should be added to samples to protect from proteolysis.
4. 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.
5. Avoid using samples with gross hemolysis or lipemia.

VIII. ASSAY PROCEDURE

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

A. Canine C-Peptide Calibrator Preparation

Use care in opening the lyophilized calibrator vial. Using a pipette, reconstitute the Canine C-Peptide Calibrator with 2 mL distilled or deionized water to give a concentration described in the analysis sheet. Invert and mix gently, let sit for five minutes then mix well.

Label seven glass tubes 1, 2, 3, 4, 5, 6, and 7. Add 0.5 mL Assay Buffer to each of the seven tubes. Prepare serial dilutions by adding 0.5 mL of the reconstituted calibrator to tube 1, mix well and transfer 0.5 mL from tube 1 to tube 2, mix well and transfer 0.5 mL from tube 2 to tube 3, mix well and transfer 0.5 mL from tube 3 to tube 4, mix well and transfer 0.5 mL from tube 4 to tube 5, mix well and transfer 0.5 mL from tube 5 to tube 6, mix well and transfer 0.5 mL from tube 6 to tube 7 and 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 $\leq -20^{\circ}\text{C}$. Avoid multiple freeze/thaw cycles.

| | Volume of Deionized Water to Add | Volume of Calibrator to Add | Calibrator Concentration ng/mL |
|--|----------------------------------|-----------------------------|---|
| | 2 mL | 0 | X (Refer to calibrator insert sheet for exact concentration) |

| Tube # | Volume of Assay Buffer to Add | Volume of Calibrator to Add | Calibrator Concentration ng/mL |
|--------|-------------------------------|------------------------------------|--------------------------------|
| 1 | 0.5 mL | 0.5 mL of reconstituted calibrator | X/2 |
| 2 | 0.5 mL | 0.5 mL of tube 1 | X/4 |
| 3 | 0.5 mL | 0.5 mL of tube 2 | X/8 |
| 4 | 0.5 mL | 0.5 mL of tube 3 | X/16 |
| 5 | 0.5 mL | 0.5 mL of tube 4 | X/32 |
| 6 | 0.5 mL | 0.5 mL of tube 5 | X/64 |
| 7 | 0.5 mL | 0.5 mL of tube 6 | X/128 |

B. Canine C-Peptide Quality Control 1 and 2 Preparation

Use care in opening the lyophilized Quality Control vials. Using a pipette, reconstitute Quality Control 1 and Quality Control 2 with **1 mL** distilled or deionized water. Invert and mix gently, let sit for five minutes or until completely dissolved then mix well.

III. ASSAY PROCEDURE (continued)

C. Assay Set-Up, Day One

1. Pipet 300 μ L of Assay Buffer to the Non-Specific Binding (NSB) tubes (3-4), 200 μ L to Reference (Bo) tubes (5-6), and 100 μ L in tubes 7 through the end of the assay.
2. Pipet 100 μ L of Calibrators and Quality Controls in duplicate (see flow chart).
3. Pipet 100 μ L of each sample in duplicate. (NOTE: Smaller volumes of sample may be used when Canine C-Peptide 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 μ L, e.g., when using 50 μ L of sample, add 50 μ L of Assay Buffer). Refer to Section IX for calculation modification.
4. Pipet 100 μ L of hydrated 125 I- Canine C-Peptide to all tubes. Important: For preparation, see Section III, Part C.
5. Pipet 100 μ L of Canine C-Peptide 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.

D. 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.
Conversion of rpm to xg:
$$xg = (1.12 \times 10^{-5}) (r) (rpm)^2$$
$$r = \text{radial distance in cm (from axis of rotation to the bottom of the tube)}$$
$$rpm = \text{revolutions per minute}$$
10. Immediately decant the supernatant 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 each. Calculate the ng/mL of Canine C-Peptide in unknown samples using automated data reduction procedures.

III. ASSAY PROCEDURE (continued)

Assay Procedure Flow Chart

| Day One | | | | | | Day Two | |
|---------|------------------|------------------------------------|-----------------------------------|-------------------------------|--|---------------------------|--|
| Set-up | Step 2 | Steps 3-4 | Step 5 | Step 6 | Step 7 | Step 8 | Steps 9-12 |
| Tube # | Add Assay Buffer | Add Calibrator / QC / Sample | Add I-125 Canine C-Peptide Tracer | Add Canine C-Peptide Antibody | Vortex, Cover, and Incubate 20-24 hrs at 4°C | Add Precipitating Reagent | Vortex and incubate 20 min. at 4°C, Centrifuge at 4°C for 20 min., Decant, and Count |
| 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 tube 7 | 100 µL | 100 µL | | 1.0 mL | |
| 9,10 | 100 µL | 100 µL of tube 6 | 100 µL | 100 µL | | 1.0 mL | |
| 11,12 | 100 µL | 100 µL of tube 5 | 100 µL | 100 µL | | 1.0 mL | |
| 13,14 | 100 µL | 100 µL of tube 4 | 100 µL | 100 µL | | 1.0 mL | |
| 15,16 | 100 µL | 100 µL of tube 3 | 100 µL | 100 µL | | 1.0 mL | |
| 17,18 | 100 µL | 100 µL of tube 2 | 100 µL | 100 µL | | 1.0 mL | |
| 19,20 | 100 µL | 100 µL of tube 1 | 100 µL | 100 µL | | 1.0 mL | |
| 21,22 | 100 µL | 100 µL of reconstituted calibrator | 100 µL | 100 µL | | 1.0 mL | |
| 23,24 | 100 µL | 100 µL of QC 1 | 100 µL | 100 µL | | 1.0 mL | |
| 25,26 | 100 µL | 100 µL of QC 2 | 100 µL | 100 µL | | 1.0 mL | |
| 27, n | 100 µL | 100 µL of unknown | 100 µL | 100 µL | | | |

IX. CALCULATIONS

A. Explanation

The calculations for Canine C-Peptide 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
$$\frac{\text{Total Binding Counts}}{\text{Total Counts}} \times 100$$

This should be 30-50%.
4. Calculate the percentage of total binding (%B/Bo) for each standard and sample
$$\%B/Bo = \frac{\text{Sample or Standard}}{\text{Total Binding}} \times 100$$
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 ng/mL of Canine C-Peptide in the unknown samples and controls by interpolation of the reference curve.

[NOTE: When sample volumes assayed differ from 100 μ L, an appropriate mathematical adjustment must be made to accommodate for the dilution factor (e.g., if 50 μ 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.

XI. ASSAY CHARACTERISTICS

A. Sensitivity

The limit of sensitivity for the Canine C-Peptide assay is 0.15 ng/mL (100 µL sample size).

B. Performance

The following parameters of assay performance are expressed as Mean \pm 3 Standard deviations.

ED₈₀=0.56 \pm .15 ng/mL

ED₅₀=2.03 \pm .54 ng/mL

ED₂₀=7.21 \pm 2.76 ng/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.

| | |
|--------------------|------|
| Canine C-Peptide | 100% |
| Human C-Peptide | 15% |
| Rat C-Peptide | ND |
| Human Proinsulin | ND |
| Porcine Proinsulin | ND |
| Bovine Proinsulin | ND |
| Human Insulin | ND |
| Glucagon | ND |

ND- Not Detectable

D. Precision

Within and Between Assay Variation

| Sample number | Mean Canine C-Peptide level (ng/mL) | Assay Variation (%CV) |
|---------------|-------------------------------------|-----------------------|
| | | Intra-assay |
| 1 | 1.07 | 3.06 |
| 2 | 4.68 | 3.33 |
| | | Inter-assay |
| 3 | 1.26 | 6.91 |
| 4 | 4.84 | 5.61 |

XI. ASSAY CHARACTERISTICS (continued)

E. Recovery

Spike and Recovery of Canine C-Peptide in Canine serum

| Serum Sample # | Canine C-Peptide Added (ng/mL) | Observed (ng/mL) | % of Recovery |
|----------------|--------------------------------|------------------|---------------|
| 1 | 0 | 0.16 | 100 |
| | 1.2 | 1.32 | 97.06 |
| | 5 | 5.24 | 101.55 |
| | 10 | 10.12 | 99.61 |
| 2 | 0 | 0.20 | 100 |
| | 1.2 | 1.40 | 100 |
| | 5 | 5.24 | 100.77 |
| | 10 | 10.92 | 107.06 |
| 3 | 0 | 0.08 | 100 |
| | 1.2 | 1.28 | 100 |
| | 5 | 5.2 | 102.36 |
| | 10 | 11.2 | 111.11 |

Varying concentrations of Canine C-Peptide were added to three different canine serum samples and the Canine C-Peptide content was determined by RIA. Mean of the observed levels from duplicate determinations in one assay are shown. Percent recovery was calculated as the observed over expected multiplied by 100.

F. Linearity

Effect of Serum Dilution

| Sample No. | Volume sampled | Observed ng/mL | Expected ng/mL | % Expected |
|------------|----------------|----------------|----------------|------------|
| 1 | 100 μ L | 9.8 | 9.8 | 100 |
| | 50 μ L | 4.96 | | 101 |
| | 25 μ L | 2.32 | | 95 |
| 2 | 100 μ L | 11.04 | 11.04 | 100 |
| | 50 μ L | 5.56 | | 101 |
| | 25 μ L | 2.68 | | 97 |
| 3 | 100 μ L | 11.28 | 11.28 | 100 |
| | 50 μ L | 5.68 | | 101 |
| | 25 μ L | 2.68 | | 95 |

Canine Serum containing varying concentrations of Canine C-Peptide were analyzed in the volumes indicated. Dilution factors of 1, 2, and 4 representing 100 μ L, 50 μ L, and 25 μ L, respectively, were applied in calculating observed concentrations.

XI. ASSAY CHARACTERISTICS (continued)**G. 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 CPM | %B/Bo | ng/mL |
|--------------------------|-------------|-------|---------|-------------|-------|-------|
| 1 | Totals | 16031 | 16131 | | | |
| 2 | " | 16231 | | | | |
| 3 | NSB | 426 | 413 | | | |
| 4 | " | 399 | | | | |
| 5 | Bo | 6854 | 6892 | 6479 | | |
| 6 | " | 6930 | | | | |
| <u>Standards</u> | | | | | | |
| 7 | 0.156 ng/mL | 6394 | 6499 | 6086 | 93.9 | |
| 8 | | 6603 | | | | |
| 9 | 0.313 ng/mL | 6279 | 6264 | 5851 | 90.3 | |
| 10 | | 6248 | | | | |
| 11 | 0.625 ng/mL | 5438 | 5442 | 5029 | 77.6 | |
| 12 | | 5445 | | | | |
| 13 | 1.25 ng/mL | 4740 | 4678 | 4265 | 65.8 | |
| 14 | | 4616 | | | | |
| 15 | 2.5 ng/mL | 3494 | 3464 | 3051 | 47.1 | |
| 16 | | 3433 | | | | |
| 17 | 5.0 ng/mL | 2344 | 2251 | 1838 | 28.4 | |
| 18 | | 2157 | | | | |
| 19 | 10.0 ng/mL | 1505 | 1440 | 1027 | 15.8 | |
| 20 | | 1374 | | | | |
| 21 | 20.0 ng/mL | 899 | 896 | 483 | 7.5 | |
| 22 | | 893 | | | | |
| <u>Controls/Unknowns</u> | | | | | | |
| 23 | | 4543 | 4472 | 4059 | 62.7 | 1.38 |
| 24 | | 4401 | | | | |
| 25 | | 2237 | 2283 | 1870 | 28.9 | 4.95 |
| 26 | | 2329 | | | | |

XII. NORMAL FASTING RANGE

Normal fasting range:

To Be Established

XIII. QUALITY CONTROLS

Good Laboratory Practice (GLP) requires that Quality Control 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 using the catalog number as the keyword.

Recommended batch analysis decision using two controls (Westgard Rules³):

1. When both controls are within ± 2 SD.
Decision: Approve batch and release analytical results.
2. When one control is outside ± 2 SD and the second control is within ± 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

| Reagents | Cat. # |
|---|---------|
| ¹²⁵ I-Canine C-Peptide (<1.5 μ Ci, 56 kBq) | 9024-HK |
| Label Hydrating Buffer (13.5mL) | LHB-PHK |
| Canine C-Peptide Calibrator (Lyophilized) | 8024-K |
| Canine C-Peptide Antibody (13 mL) | 1024-HK |
| Precipitating Reagent (130 mL) | PR-UVHK |
| QC 1&2 (Lyophilized) | 6000L-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

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.

XVI. REFERENCES

1. 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.
2. Thorell, J.I. *Scand J. Clin. Lab. Invest.* 31:187, 1973
3. Westgard, J.O., et.al. Amulti Shewhart chart for quality for quality control in clinical chemistry. *Clin. Chem.* 27:493-501, 1981.