



## **Canine C-Peptide ELISA**

### **96-Well Plate Assay**

**Cat. # EZCCP-47K,  
EZCCP47BK**

## **Canine C-Peptide ELISA KIT**

### **96-Well Plate**

**Cat. # EZCCP-47K and EZCCP47BK**

<b><u>TABLE OF CONTENTS</u></b>	<b><u>PAGE</u></b>
I. Intended Use	2
II. Principles Of Procedure	2
III. Reagents Supplied	3
IV. Storage and Stability	3
V. Reagent Precautions	3
VI. Materials Required But Not Provided	5
VII. Sample Collection And Storage	5
VIII. Reagent Preparation	6
IX. Assay Procedure	8
X. Microtiter Plate Arrangement	11
XI. Graph of Typical Reference Curve	12
XII. Assay Characteristics	12
XIII. Quality Controls	16
XIV. Troubleshooting Guide	16
XV. Replacement Reagents	16
XVI. Ordering Information	17

## I. INTENDED USE

This Canine C-Peptide ELISA kit is used for the non-radioactive quantification of C-peptide in canine serum and plasma. One kit is sufficient to measure 39 unknown samples in duplicate. ***This kit is for Research Use Only. Not for Use in Diagnostic Procedures.***

## II. PRINCIPLES OF ASSAY

This assay is a competitive ELISA based, sequentially, on: 1) capture of canine C-peptide molecules from samples/standard by a pre-titered limited amount of anti-canine C-peptide serum whose IgG fraction can be immobilized on the wells of a microtiter plate, 2) subsequent competition of binding to the immobilized anti-canine C-peptide IgG by a pre-titered biotinylated canine C-peptide, 3) washing of unbound materials from samples, 4) binding of streptavidin-horseradish peroxidase conjugate to the immobilized biotinylated canine C-peptide, 5) washing of excess free enzyme conjugates, and 7) quantification of immobilized 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 absorbance at 450 nm – 590 nm after acidification of formed products. Since the decrease in absorbance is directly proportional to the amount of captured canine C-peptide in the unknown sample because of the fixed limited amount of anti-canine C-peptide IgG, the concentration can be derived by interpolation from a reference curve generated in the same assay with reference standards of known concentrations of canine C-peptide.

### III. 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	Catalog Number	Volume	Quantity
Microtiter Plate with 2 plate sealers	EP83	-----	1 plate 2 sealers
Canine C-Peptide Standard	E8047-K	Lyophilized	1 vial
Canine C-Peptide Quality Control 1 and 2	E6047-K	Lyophilized	2 vials
Matrix Solution	EMTX-1	Lyophilized	1 vial
Assay Buffer	EAB-GLU	25 mL	1 bottle
10X Wash Buffer	EWB-HRP	50 mL	2 bottles
Biotinylated Canine C-Peptide	E1047-D	3 mL	1 bottle
Canine C-Peptide Antibody	E1047-P	3 mL	1 bottle
Enzyme Solution	EHRP-5	12 mL	1 bottle
Substrate Solution	ESS-TMB2	12 mL	1 bottle
Stop Solution	ET-TMB	12 mL	1 bottle

### IV. STORAGE AND STABILITY

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







All components are shipped and stored at 2-8°C. Reconstituted standards and controls can be frozen for future use but repeated freeze/thaw cycles should be avoided. Refer to expiration dates on all reagents prior to use. Do not mix reagents from different kits unless they have the same lot numbers.

### V. REAGENT PRECAUTIONS

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.

**See next page for Full Hazardous Labels for components of this kit.**

## Full Labels of Hazardous components:

Ingredient, Cat #		Full Label	
Canine C-Peptide ELISA Standard	E8047-K	 	<b>Warning.</b> Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Canine C-Peptide Quality Control 1 and 2	E6047-K	 	<b>Warning.</b> Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.
Matrix Solution	EMTX-1	No Symbol Required	Harmful to aquatic life with long lasting effects. Avoid release to the environment.
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.
Stop Solution	ET-TMB		<b>Warning.</b> May be corrosive to metals.

## VI. MATERIALS REQUIRED BUT NOT PROVIDED

1. Multi-channel Pipettes and pipette tips: 5-50  $\mu$ L and 50-300  $\mu$ L
2. Pipettes and pipette tips: 10  $\mu$ L-20  $\mu$ L or 20  $\mu$ L-100  $\mu$ L
3. Reagent Reservoirs
4. Polypropylene Microfuge Tubes
5. Vortex Mixer
6. De-ionized water
7. Microtiter Plate Reader capable of reading absorbency at 450 nm
8. Orbital Microtiter Plate Shaker
9. Absorbent Paper or Cloth

## VII. SAMPLE COLLECTION AND STORAGE

### A. Preparation of Serum Samples:

- Whole blood is directly drawn into a Vacutainer® serum tube that contains no anti-coagulant. Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000xg.
- Use freshly prepared serum or aliquot and store samples at  $\leq -20^{\circ}\text{C}$ .
- Avoid multiple [ $>2$ ] freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.

### B. Preparation of Plasma Samples:

- Whole blood is directly drawn into a Vacutainer® EDTA-plasma tube. Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000xg.
- Use freshly prepared plasma or aliquot and store samples at  $\leq -20^{\circ}\text{C}$ .
- Avoid multiple [ $>2$ ] freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.

### NOTE:

- A maximum of 20  $\mu$ L per well of neat serum or plasma can be used.
- All samples must be stored in polypropylene tubes. **DO NOT STORE SAMPLES IN GLASS.**
- Avoid debris, lipids and cells when using samples with gross hemolysis or lipemia.
- Care must be taken when using heparin as an anti-coagulant since an excess of heparin will provide falsely high values. Use no more than 10 IU heparin per mL of blood collected.

## VIII. REAGENT PREPARATION

### A. Canine C-Peptide Standard Preparation

1. Use care in opening the lyophilized Standard vial. Using a pipette, reconstitute the canine C-peptide Standard with 0.5 mL distilled or de-ionized water. Invert and mix gently, let sit for 5 minutes then mix well.
2. Label 5 polypropylene microfuge tubes as 1, 2, 3, 4, and 5. Add 100  $\mu$ L Assay Buffer to each of the tubes. Prepare serial dilutions by adding 100  $\mu$ L of the reconstituted standard to Tube 5, mix well and transfer 100  $\mu$ L of Tube 5 to Tube 4, mix well and transfer 100  $\mu$ L of Tube 4 to Tube 3, mix well and transfer 100  $\mu$ L of Tube 3 to Tube 2, mix well and transfer 100  $\mu$ L of the Tube 2 to Tube 1, mix well. The 0 ng/mL Canine C-Peptide standard (Background) will be Assay Buffer.

Note: Change tip for every dilution. Wet tip with standard before dispensing. Unused portions of reconstituted standard should be stored in small aliquots at  $\leq -20^{\circ}\text{C}$ . Avoid multiple freeze/thaw cycles.

Tube #	Volume of Deionized Water to Add	Volume of Standard to Add	Standard Stock Concentration
Reconstituted standard/Tube 6	0.5 mL	0	X (refer to analysis sheet for exact concentration)

Tube #	Volume of Assay Buffer to Add	Volume of Standard to Add	Standard Concentration (ng/mL)
5	100 $\mu$ L	100 $\mu$ L of reconstituted standard	X/2
4	100 $\mu$ L	100 $\mu$ L of Tube 5	X/4
3	100 $\mu$ L	100 $\mu$ L of Tube 4	X/8
2	100 $\mu$ L	100 $\mu$ L of Tube 3	X/16
1	100 $\mu$ L	100 $\mu$ L of Tube 2	X/32

## **VIII. REAGENT PREPARATION (continued)**

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

Use care in opening the lyophilized Quality Control vials. Reconstitute each Canine C-Peptide Quality Control 1 and Quality Control 2 with 0.5 mL distilled or de-ionized water and gently invert to ensure complete hydration. Unused portions of the reconstituted Quality Controls should be stored in small aliquots at  $\leq -20^{\circ}\text{C}$ . Avoid further freeze/thaw cycles.

### **C. Preparation of Wash Buffer**

Bring the 10X Wash Buffer to room temperature and mix to bring all salts into solution. Dilute 50 mL of 10X Wash Buffer with 450 mL deionized water. Store unused portion at 2-8°C for up to one month.

### **D. Preparation of Matrix Solution**

Add 1 mL distilled or de-ionized water to the bottle containing lyophilized Matrix Solution. Mix well. Allow at least 5 minutes for complete reconstitution. Any unused portion of the reconstituted Matrix Solution should be stored at  $\leq -20^{\circ}\text{C}$  for up to one month.



## IX. CANINE C-PEPTIDE ELISA ASSAY PROCEDURE

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

1. 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 the strips in an empty plate holder. Add 300 µL diluted Wash Buffer to each well of the plate. Decant Wash Buffer and remove the residual volume by inverting the plate and tapping it smartly onto absorbent towels several times. Repeat wash procedure 2 additional times. **Do not let wells dry before proceeding to the next step.** If an automated machine is used for the assay, follow the manufacturer's instructions for all washing steps described in this protocol.
2. Add 30 µL Assay Buffer to B<sub>0</sub> wells, 10 µL to Standard/Quality Control wells and 30 µL to sample wells.
3. Add 20 µL of appropriate Matrix Solution to B<sub>0</sub>, Standards and Quality Control wells (refer to Microtiter Plate Arrangement section for suggested sample order placement).
4. Add 20 µL Standards and Quality Controls to the appropriate wells.
5. Add 20 µL of sample to the appropriate wells.
6. Add 25 µL Canine C-Peptide Antibody to all wells.
7. Cover the plate with plate sealer and incubate at room temperature for 3 hours on an orbital microtiter plate shaker set to rotate at moderate speed, about 400 to 500 rpm.
8. Remove plate sealer and add 25 µL Biotinylated Canine C-Peptide to all wells. Re-cover plate with sealer and incubate at room temperature for 1 hour on an orbital microtiter plate shaker set to rotate at a moderate speed, about 400 to 500 rpm.
9. Remove plate sealer and decant reagents from the plate. Tap as before to remove residual volume in well. Wash wells 3 times with diluted Wash Buffer, 300 µL per well per wash. Decant and tap after each wash to remove residual buffer.
10. Add 100 µL Enzyme Solution to each well. Cover plate with sealer and incubate at room temperature for 30 minutes on an orbital microtiter plate shaker set to rotate at a moderate speed, about 400 to 500 rpm.

## IX. CANINE C-PEPTIDE ELISA ASSAY PROCEDURE (continued)






11. Remove sealer, decant reagents from the plate and tap plate to remove the residual volume. 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.
12. Add 100  $\mu$ L of Substrate Solution to each well, cover plate with sealer and shake on the plate shaker for approximately 5-20 minutes. Blue color should be formed in wells of the canine C-peptide standards and in samples with intensity inversely proportional to increasing concentrations of canine C-peptide.

**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.**

13. Remove sealer and add 100  $\mu$ L Stop Solution [**CAUTION: CORROSIVE SOLUTION**] and gently shake plate by hand to ensure complete mixing of solution in all wells. The blue color should turn to 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 590nm in a plate reader within 5 minutes and ensure that there are no air bubbles in any well. Record the difference of absorbance units. The absorbance of the lowest Canine C-Peptide standard should be approximately 2.0 - 3.0, or not to exceed the capability of the plate reader used.

Note: When sample volumes assayed differ from 20  $\mu$ L, an appropriate mathematical adjustment must be made to accommodate for the dilution factor (e.g., if 10  $\mu$ L of sample is used, then calculated data must be multiplied by 2). When sample volume assayed is less than 20  $\mu$ L, compensate for the volume deficit with the Matrix Solution provided.

### Assay Procedure for Canine C-Peptide ELISA Kit (Cat. # EZCCP-47K and EZCCP47BK)

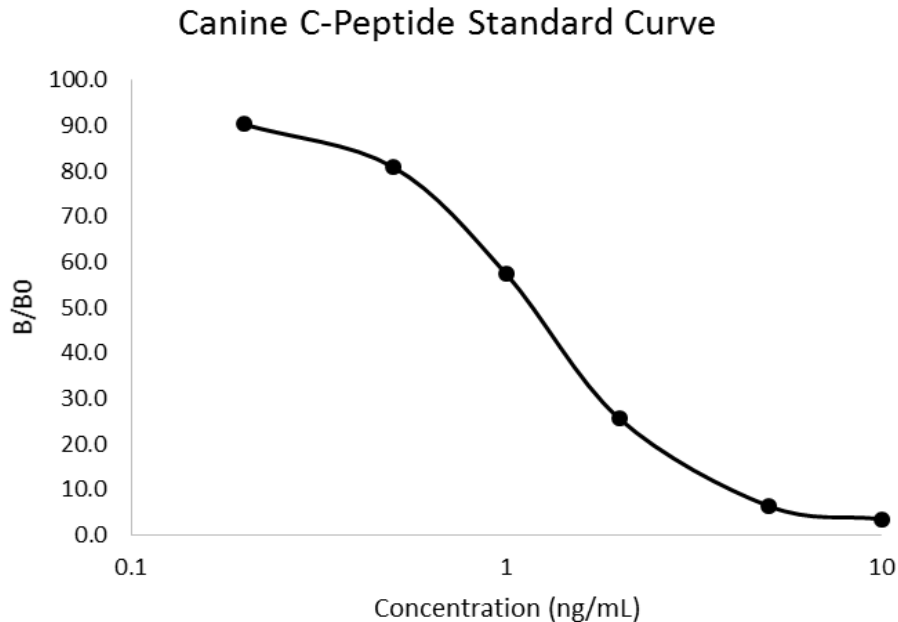
	Step 1	Step 2	Step 3	Step 4-5	Step 6	Step 7	Step 8	Step 9	Step 10	Step 11	Step 12-13			
Well #	Wash plate 3X with 300 µL 1X Wash Buffer. Remove residual buffer by tapping smartly on absorbent towels.	Assay Buffer	Matrix Solution	Standards/ QCs/ Samples	Canine C-Peptide Antibody	Seal, Agitate, Incubate 3 hrs at Room Temperature.	Biotinylated Canine C-Peptide	Seal, Agitate, Incubate 1 hour at Room Temperature. Wash 3X with 300 µL Wash Buffer.	Enzyme Solution	Seal, Agitate, Incubate 30 minutes at Room Temperature. Wash 3X with 300 µL Wash Buffer.	Substrate Solution	Seal, Agitate, Incubate 5 – 20 minutes at Room Temperature.	Stop Solution	Read Absorbance at 450 nm and 590 nm.
A1, B1		30 µL	20 µL	--	25 µL		25 µL		100 µL		100 µL			
C1, D1		10 µL	20 µL	20 µL of Tube 1										
E1, F1		10 µL	20 µL	20 µL of Tube 2										
G1, H1		10 µL	20 µL	20 µL of Tube 3										
A2, B2		10 µL	20 µL	20 µL of Tube 4										
C2, D2		10 µL	20 µL	20 µL of Tube 5										
E2, F2		10 µL	20 µL	20 µL of Tube 6										
G2, H2		10 µL	20 µL	20 µL of QC1										
A3, B3		10 µL	20 µL	20 µL of QC2										
C3, D3		30 µL	--	20 µL of Sample										
E3, F3		30 µL	--	20 µL of Sample										
G3, H3 Etc.		30 µL	--	20 µL of Sample										

## X. MICROTITER PLATE ARRANGEMENT

### Canine C- Peptide ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	B <sub>0</sub>	Tube 4	QC2									
B	B <sub>0</sub>	Tube 4	QC2									
C	Tube 1	Tube 5	Sample 1									
D	Tube 1	Tube 5	Sample 1									
E	Tube 2	Reconstituted Standard	Sample 2									
F	Tube 2	Reconstituted Standard	Sample 2									
G	Tube 3	QC1	Sample 3									
H	Tube 3	QC1	Sample 3									

## XI. GRAPH OF TYPICAL REFERENCE CURVE



Typical Standard Curve, not to be used to calculate data.

## XII. ASSAY CHARACTERISTICS

### A. Sensitivity

The Minimum Detectable Concentration (MinDC) of Canine C-Peptide is 0.24 ng/mL. It is calculated by using MILLIPLEX® Analyst 5.1. It measures the true limits of detection for an assay by mathematically determining what the empirical MinDC would be if an infinite number of standard concentrations were run for the assay under the same conditions. This reported value is the mean plus 2 standard deviations of the MinDC of multiple assays (n= 8).

### B. Specificity

Canine C-Peptide	100%
Human C-Peptide	0.8%
Porcine C-Peptide	ND
Rat C-peptide 1	ND
Rat C-peptide 2	ND
Human Proinsulin	ND
Porcine Proinsulin	ND
Bovine Proinsulin	ND
Human Insulin	ND
Glucagon	ND

## XII. ASSAY CHARACTERISTICS (continued)

### C. Precision

#### Intra-Assay Variation

	Mean Canine C-Peptide Levels (ng/mL)	Intra-Assay %CV
1	0.39	<15
2	1.95	<10

#### Inter-Assay Variation

	Mean Canine C-Peptide Levels (ng/mL)	Inter-Assay %CV
1	0.39	<10
2	1.96	<10

The assay variations of EMD Millipore's Canine C-Peptide ELISA kit was studied on two samples at two levels on the Canine C-Peptide standard curve. The mean intra-assay variation was calculated from results of eight determinations of the indicated samples. The mean inter-assay variations of each sample were calculated from results of 6 separate assays with duplicate samples in each assay.

## XII. ASSAY CHARACTERISTICS (continued)

### D. Spike Recovery of Canine C-Peptide in Assay Samples

Sample	Canine C-Peptide Added (ng/mL)	Expected (ng/mL)	Observed (ng/mL)	Recovery
1	0.5	1.17	1.12	96
	1	1.67	1.54	92
	2	2.67	2.38	89
2	0.5	0.84	0.87	104
	1	1.34	1.26	94
	2	2.34	2.11	90
3	0.5	1.07	1.06	99
	1	1.57	1.47	94
	2	2.57	2.4	93
4	0.5	0.83	0.86	104
	1	1.33	1.31	98
	2	2.33	2.28	98
5	0.5	1.01	0.99	98
	1	1.51	1.41	93
	2	2.51	2.2	88
Average				95

Varying amounts of C-Peptide were added to individual canine serum and plasma samples and the resulting C-peptide content of each sample was assayed by the Canine C-Peptide ELISA. The recovery = [(observed C-peptide/ (spiked C-peptide concentration + basal C-peptide level)) x 100%.

## XII. ASSAY CHARACTERISTICS (continued)

### E. Linearity of Sample Dilution

Sample	Volume (µL)	Expected (ng/mL)	Observed (ng/mL)	Expected
1	20	5.53	5.53	
	10	2.77	2.84	103
	5	1.39	1.55	112
	2.5	0.69	0.79	114
2	20	4.84	4.84	
	10	2.42	2.7	112
	5	1.21	1.43	118
	2.5	0.61	0.66	108
3	20	4.8	4.8	
	10	2.4	2.63	110
	5	1.2	1.42	118
	2.5	0.6	0.62	103
4	20	4.95	4.95	
	10	2.48	2.67	108
	5	1.24	1.42	115
	2.5	0.62	0.65	105
5	20	5.49	5.49	
	10	2.75	2.78	101
	5	1.37	1.45	106
	2.5	0.68	0.65	96
Average				109

Five canine serum and plasma samples with the indicated sample volumes were assayed. Required amounts of serum matrix were added to compensate for lost volumes below 20 µL. The resulting dilution factors of neat, 2, 4 and 8 representing 20 µL, 10 µL, 5 µL and 2.5 µL sample volumes assayed, respectively, were applied in the calculation of observed Canine C-Peptide concentrations. % expected = (observed/expected) x 100%.



### XIII. QUALITY CONTROLS

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).

### XIV. TROUBLESHOOTING GUIDE

1. 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.
2. Throughout the assay the operator should adhere strictly to the procedures with good laboratory practice.
3. 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.
4. Avoid cross contamination of any reagents or samples to be used in the assay.
5. Make sure all reagents and samples are added to the bottom of each well.
6. 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.
7. Remove any air bubbles formed in the well after acidification of substrate solution because bubbles interfere with spectrophotometric readings.
8. High signal in background or blank wells could be due to 1.) cross well contamination by standard solution or sample or 2.) inadequate washing of wells with Wash Buffer or 3.) overexposure to light after substrate has been added.

### XV. REPLACEMENT REAGENTS

Reagents	Cat. #
ELISA Plate	EP83
10X HRP Wash Buffer Concentrate	EWB-HRP
Canine C-Peptide ELISA Standard	E8047-K
Canine C-Peptide Quality Control 1 and 2	E6047-K
Matrix Solution	EMTX-1
Assay Buffer	EAB-GLU
Canine C-Peptide Antibody	E1047-P
Biotinylated Canine C-Peptide	E1047-D
Enzyme Solution	EHRP-5
Substrate Solution	ESS-TMB2
Stop Solution	ET-TMB
10 pack of Canine C-Peptide ELISA Kits	EZCCP47BK

## **XVI. 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).