

Cat. # MK149

For Research Use

---

**TAKARA**

**Pig Glu-Osteocalcin  
EIA Kit**

---

Product Manual

v201607Da

---

## Table of Contents

I. Description .....	3
II. Principle .....	3
III. Components .....	4
IV. Materials Required but not Provided .....	4
V. Storage .....	4
VI. Intended Use.....	4
VII. Protocol.....	5
VIII. Performance.....	7
IX. Experimental Examples .....	10
X. Related Products .....	12
XI. Precautions.....	12

## I. Description

The Pig Glu-Osteocalcin EIA Kit is a quantitative kit that enables specific and highly sensitive assay of decarboxylated osteocalcin released from porcine bone tissues by enzymes in osteoclasts or uncarboxylated Glu-osteocalcin (inactive osteocalcin) produced by osteoblasts. The capture antibody (plate-bound antibody) is a plate-bound solid-phased monoclonal antibody that specifically recognizes the Glu residues at positions 21 and 24 on osteocalcin. It is paired with a labeled antibody – a monoclonal antibody for detecting porcine osteocalcin.

The concurrent measurement of porcine active osteocalcin (Gla-osteocalcin) may be achieved with the Pig Gla-Osteocalcin EIA Kit (Cat. #MK139) to monitor both bone formation and bone resorption based on a relative evaluation of Gla/Glu-osteocalcins.

Osteocalcin (OC) comprises 49 amino acids, including 2 to 3  $\gamma$ -carboxyglutamate residues (Gla), and has a molecular weight of approximately 5,900. It is known as a vitamin K-dependent calcium-binding non-collagen protein. Osteocalcin is an osteoblast-specific marker as it is produced only by osteoblasts. The Gla-osteocalcin, in particular, is a marker of osteogenesis.

[Primary amino acid structures of osteocalcins of various species]

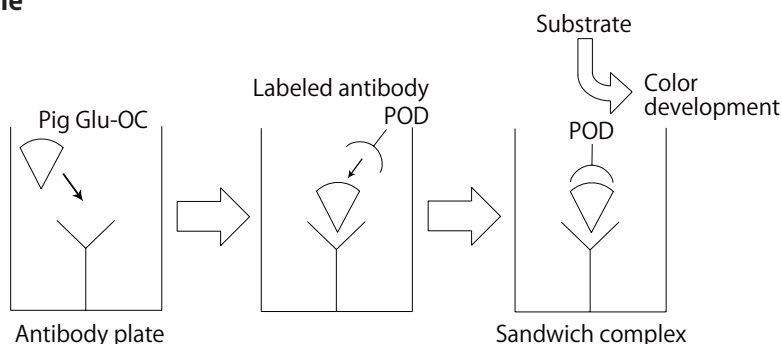
		10	20	30	40	50
Human	1	YLYQWLGAPV	PYPDPLEPRR	EVCELNPDCD	ELADHIGFQE	AYRRFYGP-V
Bovine	1	YLDHWLGAPA	PYPDPLEPKR	EVCELNPDCD	ELADHIGFQE	AYRRFYGP-V
Rat	1	YLNNGLGAPA	PYPDPLEPHR	EVCELNPDCD	ELADHIGFQD	AYKRIYGTTV
Mouse	1	YL----GASV	PSPDPLEPTR	EQCELNPACD	ELSDQYGLKT	AYKRIYGITI
Chicken	1	YAQDSGVAGA	P-PNPLEAQR	EVCELSPDCD	ELADQIGFQE	AYRRFYGP-V
Monkey	1	YLYQWLGAPA	PYPDPLEPKR	EVCELNPDCD	ELADHIGFQE	AYRRFYGP-V
Pig	1	YLDHGLGAPA	PYPDPLEPRR	EVCELNPDCD	ELADHIGFQE	AYRRFYGI-A

In bone formation, the dynamic osteogenesis in growing young animals is called “modeling” (new construction). In matured animals beyond the growing phase, bone morphology undergoes no apparent changes and remains stable, but a certain percentage of the bone are constantly being replaced. This process is called “remodeling” (reconstruction).

Efficacy assessments of osteoporosis drugs require the use of 2 types of animal models to allow evaluation for bone modeling and remodeling. Rodents, such as mouse and rat, are commonly used animal models to evaluate modeling, while monkey and miniature pig are used frequently to evaluate remodeling. These animal models are essential in the developments and efficacy assessments of potential therapeutic agents.

Osteocalcin is also a hormone that plays a role in insulin regulation and in male fertility. The undercarboxylated form (ucOC), in particular, is responsible for regulation of glucose metabolism. This kit enables researchers to study the role that ucOC plays in metabolic disease.

## II. Principle



**III. Components**

(1) Antibody Coated Microtiterplate Anti-Glu-OC monoclonal antibody-coated plate (96 wells: 8 wells x 12 Strips)	1 plate
(2) Antibody-POD Conjugate (lyophilized) Peroxidase-Labeled Anti-Pig-OC Monoclonal Antibody	for 11 ml
(3) Standard (lyophilized) Pig Glu Osteocalcin Full-length Peptide 32 ng	for 1 ml
(4) Sample Diluent BlockAce-containing PBS (with preservative)	11 ml x 2
(5) Substrate Solution (TMBZ) 3, 3', 5, 5' Tetramethylbenzidine Solution	12 ml

**IV. Materials Required but not Provided**

- Wash and Stop solution for ELISA without Sulfuric Acid (Cat. #MK021)  
Contains wash solution (10X PBS, 50 ml x 5 tubes; Tween 20, 3 ml) and reaction stop solution (60 ml).
  - \* This product is a stop solution for peroxidase reactions without 1N sulfuric acid.
  - \* 1N sulfuric acid can be used as a stop solution. Handle 1N sulfuric acid with caution.
- Pipette, micropipette, and tips
- Microplate reader (capable of measuring absorbance of up to 3.5 when set to 450 nm)

**V. Storage**            4°C**VI. Intended Use**

Quantitative determination of Glu-type osteocalcin (Pig Glu-OC) in porcine biological samples.

**VII. Protocol****1. Sample**

- Suitable samples include porcine serum, plasma, ascite fluid, urine, and cell culture supernatant, etc.
- Samples may be stored up to 12 hours at 4°C. If the assay will be performed longer than 12 hours after sample preparation, then store samples frozen at -20°C. Use (4) Sample Diluent for dilution if necessary.
- The recommended dilution for porcine serum samples is 2 - 4-fold. (Investigate the optimum dilution ratio before assaying a sample for the first time.)
- As this product cross-reacts slightly with bovine antigens, any bovine-serum-supplemented medium in a cultured sample may interfere with the assay. Switch to a serum-free medium.
- This product cross-reacts slightly with rat antigens, resulting in low sensitivity. It is therefore unsuitable for assaying samples of rat origin. We recommend Rat Glu-Osteocalcin High Sensitive EIA Kit (Cat. #MK146) for assaying rat samples.
- This product does not react to mouse, rabbit, or human antigens.

**2. Preparation of Solutions**

- Antibody Coated Microtiter Plate  
Allow the (1) Anti-Glu OC Monoclonal Antibody Coated Plate to reach room temperature unopened in its package before use.
- POD-labeled Antibody Solution  
Reconstitute (2) Antibody - POD Conjugate with 11 ml of distilled water.  
Once reconstituted, it is stable for up to 1 week at 4°C. For longer storage, freeze at -20°C, at which it is stable for up to 1 month. Once thawed, it may not be returned to frozen storage.
- Pig Glu-OC standard solution  
Add 1 ml of distilled water to the lyophilized (3) Standard to reconstitute it (32.0 ng/ml). Dilute the Standard with (4) Sample Diluent before use to prepare fresh serial dilutions of Standard Solution at concentrations of 16.0, 8.0, 4.0, 2.0, 1.0, and 0.5 ng/ml. Use Sample Diluent as the 0-concentration standard.  
The dissolved Pig Glu-OC standard solution (32.0 ng/ml) is stable for up to 1 week after preparation when stored at 4°C, or for up to 1 month at -20°C. Once thawed, however, it may not be returned to frozen storage.
- Substrate Solution  
Return (5) Substrate Solution (TMBZ) to room temperature before use. It is supplied ready to use. Check before use that the Substrate Solution has not developed a dark blue color. A reaction with metal ions will result in coloration; make sure it is not contaminated with any tap water.  
If the Substrate Solution will be used for several assays, divide it into aliquots of the required volume in advance.

- Stop solution  
Use the Stop solution included in Wash and Stop Solution for ELISA without Sulfuric Acid (Cat. #MK021) directly.  
\* Because this is highly viscous, mix well using a plate mixer after its introduction.
- PBS with 0.1% Tween 20 for washing  
Dilute the 10X PBS included in Wash and Stop solution for ELISA without Sulfuric Acid (Cat. #MK021) 10 fold with distilled water, and then add Tween 20 to a final concentration of 0.1%.  
For 96 reactions performed with this kit, 300 ml of washing solution is required.

### 3. Procedure

Assay samples in duplicate.

Return each reagent in the kit and samples to room temperature and make sure solutions are mixed uniformly without creating bubbles before use.

1. Prepare reagents and samples (100  $\mu$ l each) in a separate 96-well plate in advance so that they can be added to the (1) Antibody Coated Microtiterplate quickly (within 5 minutes) using an 8-channel pipette or similar apparatus. In order to provide highly reliable results, it is recommended to place serial dilutions of the Standard Solution in the 1st and 12th rows. Perform this reaction at room temperature (20 - 30°C) for 1 hour; incubation at 37°C may compromise antigenicity.  
[First reaction]
2. Discard reaction mixtures, followed by 3 washes with Washing Buffer. Then add 100  $\mu$ l of the POD-labeled Antibody Solution per well using an 8-channel pipette and allow to react for 1 hour at room temperature (20 - 30°C).  
[Second reaction]
3. Discard reaction mixtures, followed by 4 washes with Washing Buffer. Then add 100  $\mu$ l of (5) Substrate Solution (TMBZ) per well using an 8-channel pipette and allow to react at room temperature (20 - 30°C) for 10 - 15 minutes. [Third reaction]
4. Add 100  $\mu$ l of Stop Solution to each well to stop the reaction in the same order as for (5) Substrate Solution (TMBZ). Then mix well.
5. Use distilled water as a control to make zero adjustment and measure absorbance at 450 nm.  
The color is stable for up to 1 hour after reaction termination.
6. Plot a standard curve based on the results obtained from the Standard Solutions (with concentration as x-axis and absorbance as y-axis) and use it to determine the corresponding concentrations of Pig Glu-OC based on the sample's absorbance.

#### Note:

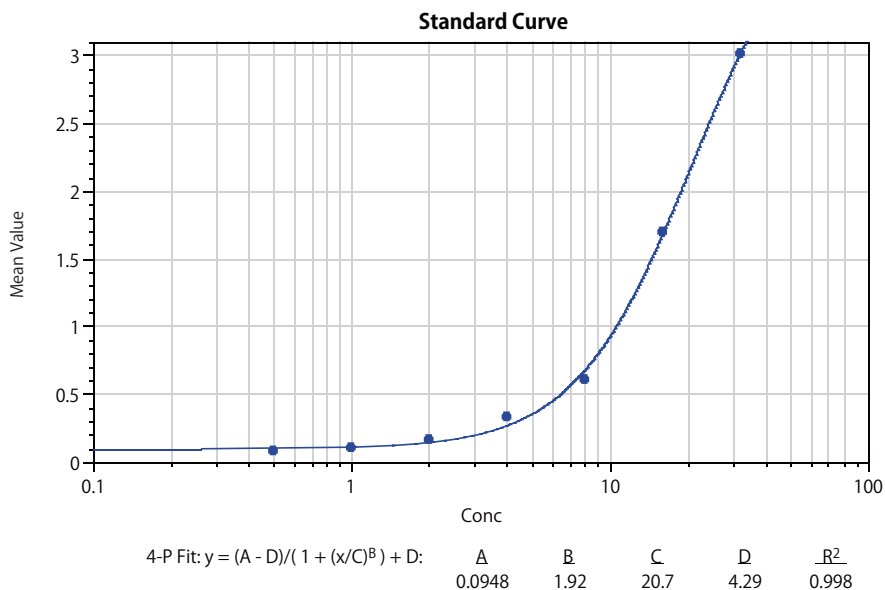
- Cover the plate with film or the like to prevent evaporation of solutions during reactions at room temperature or in an incubator.
- It is recommended that the Washing Buffer be completely discarded to get rid of the residual fluid.

**VIII. Performance**

**1. Standard curve (Pig Glu Osteocalcin EIA Kit)**

The following shows a typical standard curve of this kit as an example.  
The standard curve for calculation needs to be established in each assay.

Limit of Detection: 0.5 ng/ml



Pig Glu-OC concentration (ng/ml)	32.0	16.0	8.0	4.0	2.0	1.0	0.5	0.0
A <sub>450</sub>	3.012	1.701	0.611	0.339	0.172	0.103	0.081	0.057

(Color development time: 15 min)

**2. Reproducibility**

<Intra-assay precision test (n=16)>

A reproducibility test was performed with 16 replicates, using 3 different concentrations of porcine serum.

Sample	Mean (ng/ml)	SD	CV (%)
control A	14.9	0.99	6.8
control B	7.6	0.50	6.5
control C	4.2	0.51	12.1

<Inter-assay precision test (n=3)>

The reproducibility test was performed with triplicates, by assaying 3 different concentrations of sample over 3 days.

Sample	Mean (ng/ml)	SD	CV (%)
control D	17.3	1.50	8.7
control E	8.5	0.40	4.7
control F	4.4	0.21	4.7

**3. Recovery test**

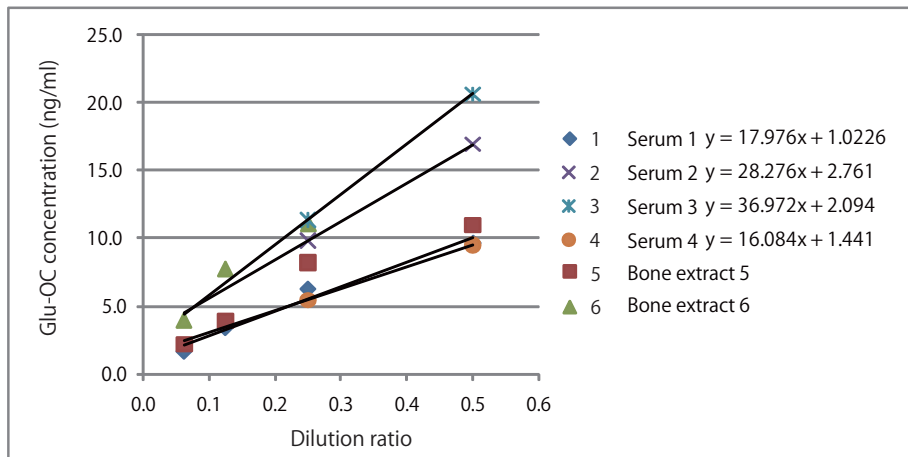
Equal volumes of samples in different concentrations were combined and assayed. The result of each mixture was compared with the theoretical value to determine the recovery rate. The mean recovery rate was 109.1%.

Sample A	Sample B	Theoretical Value (A+B)/2	Assay Result	Recovery Rate (%)
26.8	22.0	24.4	25.1	102.8
13.0	22.0	17.5	19.0	108.7
5.2	22.0	13.6	14.8	108.7
2.3	22.0	12.2	12.1	99.5
26.8	8.2	17.5	16.9	96.5
13.0	8.2	10.6	9.7	91.7
5.2	8.2	6.7	7.3	108.6
2.3	8.2	5.3	6.4	121.4
26.8	1.4	14.1	16.0	113.5
13.0	1.4	7.2	8.2	114.4
5.2	1.4	3.3	4.0	120.9
2.3	1.4	1.9	2.4	129.0
26.8	4.9	15.9	15.6	98.2
13.0	4.9	9.0	8.4	93.8
5.2	4.9	5.1	5.7	111.8
2.3	4.9	3.6	4.6	126.1

Unit: ng/ml

**4. Linearity of Porcine Serum**

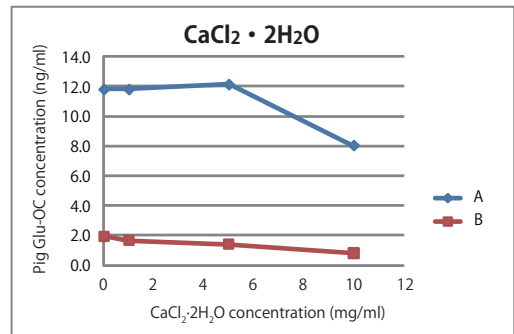
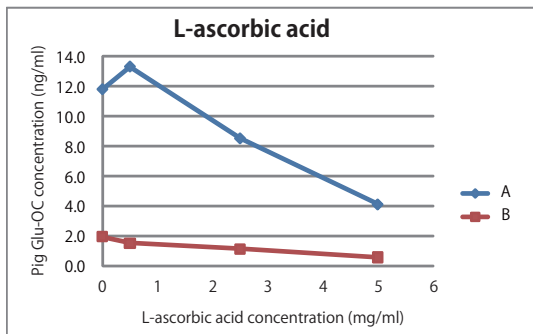
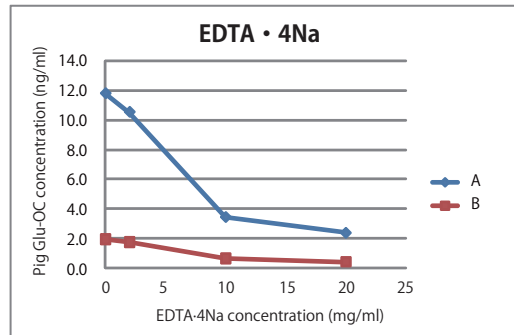
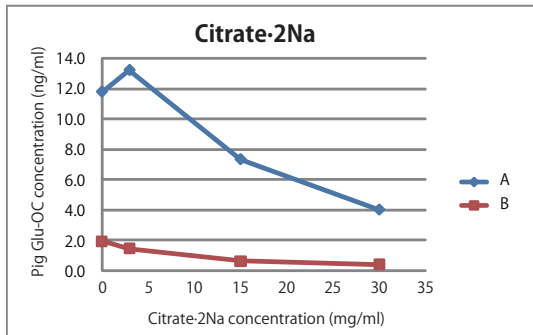
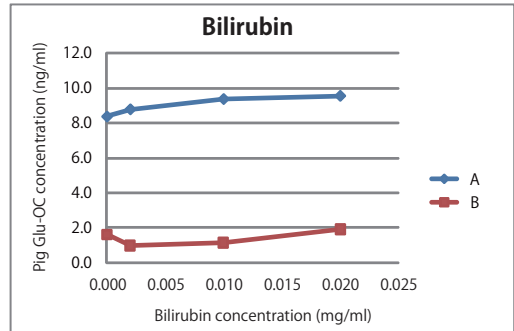
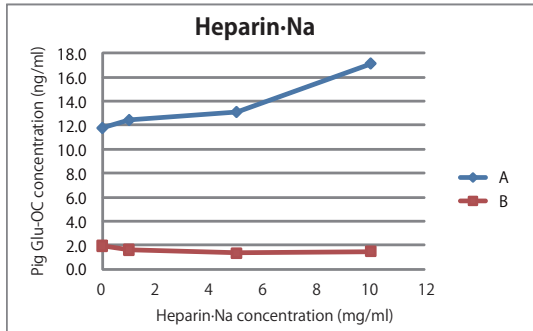
Serum samples from 8-month-old miniature pigs (1 - 4) and porcine bone extract (5, 6) were assayed. For serum samples, it is preferable to assay 2- or 4-fold dilutions based on the range of the standard curve. For porcine bone extracts, linearity was seen only at dilutions of 4-fold or higher because of the high pH in the extractant, which contained sodium carbonate (pH 9.5).





5. Effects of Coexisting Substances

1 part by volume of the substance being tested was added to 9 parts by volume of osteocalcin Standard Solution (2 concentrations: A and B), and the effects on the reaction were examined. The final concentration of the test substance is shown on the horizontal axis of the graph. The concentration of Glu-osteocalcin measured is shown on the vertical axis. (Unit: ng/ml)



**IX. Experimental Examples**

**1. Assay of miniature pig serum samples**

We monitored the individual serum Glu-osteocalcin levels in 4 miniature pigs (2 males, 2 females) aged 8 to 10 months old. Assayed samples were diluted 2- and 4-fold.

Miniature Pig		Dilution Ratio	
		2-Fold	4-Fold
Animal A (female)	8 months	16.0	10.2
	9 months	19.2	11.4
	10 months	9.9	6.5
Animal B (female)	8 months	9.1	5.4
	9 months	4.2	2.9
	10 months	5.0	3.5
Animal C (male)	8 months	22.1	14.4
	9 months	10.9	6.4
	10 months	9.1	5.8
Animal D (male)	8 months	19.2	12.7
	9 months	12.2	7.8
	10 months	10.9	6.7

[Result]

The levels of Glu-osteocalcin detected at 8, 9, and 10 months after birth declined gradually, suggesting the bone metabolic turnover had peaked over the growth phase and begun to settle down.

**2. Monitoring processed food for components of pig bone-origin**

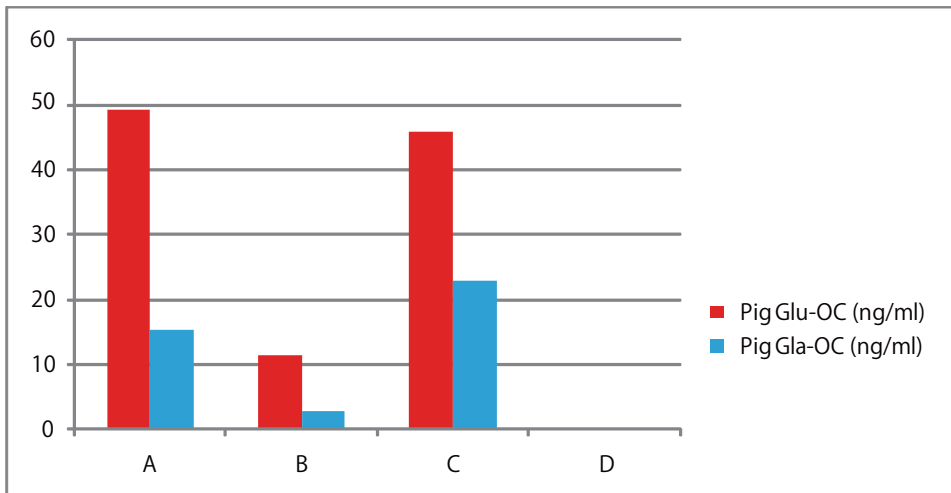
We measured the Gla- and Glu-osteocalcin levels in pig bone extracts contained in processed food.

The powdered soups that came with 4 different types of instant food (A - D) were dissolved according to the recipes to prepare stock solutions, which were assayed for Gla- and Glu-osteocalcin levels.

With heating a part of their manufacturing process, the Gla-osteocalcin in these powdered soups may have become decarboxylated depending on the heating condition. However, the Gla- and Glu-osteocalcin level in powdered soups A and C was detected with this kit.

Gla-OC: Pig Gla-Osteocalcin EIA Kit (Cat. #MK139) was used for assay

Glu-OC: Pig Glu-Osteocalcin EIA Kit (Cat. #MK149) was used for assay



[Result]

The results demonstrated that this kit may be used to periodically monitor the content of pig-bone components during the manufacturing of processed food.

**X. Related Products**

Pig Gla-Osteocalcin EIA Kit (Cat. #MK139)  
Gla-Type Osteocalcin (Gla-OC) EIA Kit (Cat. #MK111)  
Rat Gla-Osteocalcin High Sensitive EIA Kit (Cat. #MK126)  
Rat Glu-Osteocalcin High Sensitive EIA Kit (Cat. #MK146)  
Human Gla-Osteocalcin High Sensitive EIA Kit (Cat. #MK128)  
Anti-Osteocalcin (Bovine) (Cat. #M041 - M044)  
Anti-Osteocalcin (Human) (Cat. #M171, M184)  
Anti-Osteocalcin ((Mouse) (Cat. #M188)  
Anti-Osteocalcin (Mouse), polyclonal (Cat. #M173)  
Anti-Osteocalcin (Rat) (Cat. #M185 - M187)  
Osteoblast Inducer Reagent (Cat. #MK430)  
TRACP & ALP Assay Kit (Cat. #MK301)  
TRACP & ALP Double Stain Kit (Cat. #MK300)  
Anti-Rat Bone Specific Alkaline Phosphatase, Polyclonal (Cat. #M190)  
Wash and Stop Solution for ELISA without Sulfuric Acid (Cat. #MK021)

**XI. Precautions**

1. Do not mix-use kits or reagents from different lots.
2. Do not expose (5) Substrate Solution (TMBZ) to strong light during storage or incubation. Avoid contact of Substrate Solution and Stop Solution with skin or mucous membranes. If these reagents come into contact with skin, wash thoroughly with water.
3. Use metal-free pipettes when handling (5) Substrate Solution (TMBZ).
4. Do not use (5) Substrate Solution (TMBZ) that has developed color.
5. Each reaction varies depending on time and temperature. Therefore, a new standard curve must be established for each assay.
6. Handle blood samples with great care.

---

**NOTE:** This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc.

Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from TAKARA BIO INC.

If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at [www.takara-bio.com](http://www.takara-bio.com).

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.

---