

# Bovine IFN $\gamma$ ELISA Reagent Kit

## ESS0026B

| Number   | Description   |
|----------|---|
| ESS0026B | <p><b>Bovine IFN<math>\gamma</math> ELISA Reagent Kit</b>, pre-titered coating and detection antibodies, recommended buffers and specific assay protocol optimized for the quantitative measurement of bovine IFN<math>\gamma</math> in cell culture supernatants</p> <p>Kit provides sufficient reagents for approximately five 96-well plates, provided the Bovine IFN<math>\gamma</math> ELISA Reagent Kit Protocol is followed.</p> |

| Kit Contents   | Size    | Assay Dilution |
|--|---------|----------------|
| Anti-Bovine IFN $\gamma$ Coating Antibody            | 0.625mL | 1:100          |
| Lyophilized Recombinant Bovine IFN $\gamma$ Standard | 5 vials | See vial label |
| Anti-Bovine IFN $\gamma$ Detection Antibody          | 0.625mL | 1:100          |
| Streptavidin-HRP                                     | 0.25mL  | 1:400          |
| Substrate Solution                                   | 55mL    | Ready to use   |
| Stop Solution, 0.16M Sulfuric Acid                   | 55mL    | Ready to use   |

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

**Storage:** Upon receipt store the kit at 2-8°C.

### Introduction

The Invitrogen™ Bovine IFN $\gamma$  pre-titered coating and detection antibodies, recommended buffers, and specific assay protocol have been optimized for the quantitative measurement of bovine IFN $\gamma$  in cell culture supernates.

### Materials Required

- 8-well strip plates, clear, corner notched (Product No. 15031)
- Plate sealers for 96-well plates (Product No. 15036)
- Reagent reservoir, sterile, 50mL capacity, 40pk (Product No. 15075)
- Microtiter plate reader with software capable of measurement at or near 450nm
- Plate washer—automated or manual (squirt bottle, manifold dispenser, or equivalent)
- Distilled or deionized water
- Calibrated adjustable precision pipettes and glass or plastic tubes for diluting solutions; beakers, flask and cylinders for preparation of reagents

---

## ELISA Reagent Kit Buffers

- D-PBS: 0.008M sodium phosphate, 0.002M potassium phosphate, 0.14M sodium chloride, 0.01M potassium chloride, pH 7.4, 0.2µm filtered (e.g., Thermo Scientific™ BupH™ Modified Dulbecco's Phosphate Buffered Saline Packs, Product No. 28374)
- Carbonate-bicarbonate Buffer: 0.2M sodium carbonate-bicarbonate buffer, pH 9.4, 0.2µm filtered (e.g., BupH™ Carbonate/Bicarbonate Buffer, Product No. 28382)
- Blocking Buffer: 4% bovine serum albumin (BSA), 5% sucrose in D-PBS, 0.2µm filtered OR ELISA Blocker Blocking Buffer, Product No. N502
- Reagent Diluent: 4% BSA in D-PBS (pH 7.4), 0.2µm filtered
- Wash Buffer: 0.05% Tween™-20 Detergent (e.g., 0.5% Thermo Scientific™ Surfact-Amps™ 20 Detergent Solution, Product No. 28320) in D-PBS, pH 7.4 OR ELISA Wash Buffer (30X), Product No. N503

**Note:** Mix new solution daily.

## Assay Protocol

Kit components are titrated to give optimal results using the Bovine IFN $\gamma$  ELISA Reagent Kit Protocol for cell culture supernatants. Any change, including component concentration, volumes, incubation times or temperatures, buffer content or number of wash steps may significantly affect the ELISA results and require optimization to give the best results.

**Note:** Allow all reagents and buffers to equilibrate to room temperature (22-25°C) before use. Thaw one aliquot of coating and detecting antibody for each plate. Do not use a water bath.

### A. Plate Preparation

1. Dilute the Coating Antibody 1:100 in carbonate-bicarbonate buffer by adding 110µL Coating Antibody to 10.89mL of carbonate-bicarbonate buffer.
2. Add 100µL of diluted Coating Antibody to each well. Cover plate with plate sealer and incubate overnight at room temperature.
3. Aspirate Coating Antibody solution and add 300µL of Blocking Buffer to each well. Cover plate with plate sealer and incubate for 1 hour at room temperature.
4. Aspirate Blocking Buffer and proceed to assay or allow to dry overnight at room temperature. When sealed with desiccant, plates can be stored at 2-8°C for 6 months.

### B. Assay Procedure

1. Reconstitute standard with Reagent Diluent with volume stated on vial label. The concentration of the reconstituted standard is 4000pg/mL.
2. Dilute reconstituted standard 1:2 in Reagent Diluent to prepare top Standard (2000pg/mL). Using Reagent Diluent, prepare 1:2 serial dilutions of top Standard and dilute any supernatant expected to read above the top standard. Add 100µL of sample or Standard to each well. Cover plate with plate sealer and incubate for 1 hour at room temperature.
3. Aspirate and wash three times with Wash Buffer using 300µL per well.
4. Dilute the Detection Antibody 1:100 in Reagent Diluent by adding 110µL of Detection Antibody to 10.89mL of Reagent Diluent.
5. Add 100µL of Detection Antibody to each well. Cover plate with plate sealer and incubate for 1 hour at room temperature.
6. Aspirate and wash three times with Wash Buffer, using 300µL per well.
7. Dilute Streptavidin-HRP 1:400 in Reagent Diluent by adding 30µL of Streptavidin-HRP to 12mL of Reagent Diluent.

8. Add 100µL of diluted Streptavidin-HRP reagent to each well. Cover plate with plate sealer and incubate for 30 minutes at room temperature.
9. Aspirate and wash three times with Wash Buffer, using 300µL per well.
10. Add 100µL of Substrate Solution to each well. Cover plate with plate sealer and incubate in the dark for 20 minutes at room temperature.
11. Stop the reaction by adding 100µL of Stop Solution to each well.
12. Measure the absorbance at  $A_{450}$  minus  $A_{550}$ .

### C. Absorbance Measurement

Measure absorbance on an ELISA plate reader set at 450nm and 550nm. Subtract 550nm values from 450nm values to correct for optical imperfections in the microplate. If an absorbance at 550nm is not available, measure the absorbance at 450nm only.

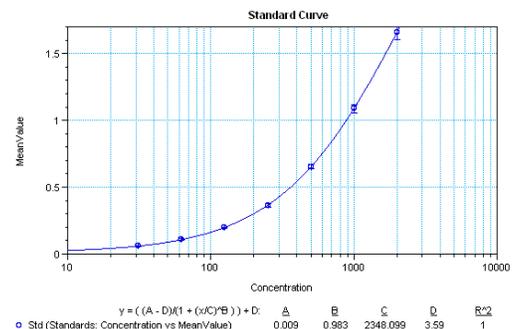
**Note:** When the 550nm measurement is omitted, absorbance values will be higher.

**Note:** Evaluate the plate within 30 minutes of stopping the reaction.

### D. Calculation of Results

- The standard curve is used to determine bovine IFN $\gamma$  amount in an unknown sample. Generate the standard curve by plotting the average absorbance obtained for each Standard concentration on the vertical (Y) axis vs. the corresponding bovine IFN $\gamma$  concentration (pg/mL) on the horizontal (X) axis.
- Calculate results using graph paper or curve-fitting statistical software. Determine the bovine IFN $\gamma$  amount in each sample by interpolating from the absorbance value (Y-axis) to bovine IFN $\gamma$  concentration (X-axis) using the standard curve.
- If the test sample was diluted, multiply the interpolated value obtained from the standard curve by the dilution factor to calculate pg/mL of bovine IFN $\gamma$  in the sample.
- Absorbance values obtained for duplicates should be within 10% of the mean value. Carefully consider duplicate values that differ from the mean by greater than 10%.

### Standard Curve Example



**Standard curve based on data obtained using the Bovine IFN $\gamma$  ELISA Reagent Kit Protocol.**

**NOTE: This standard curve is for demonstration only. A standard curve must be run with each assay.**

### Performance Characteristics

**Specificity:** This ELISA is specific for the measurement of natural and recombinant bovine IFN $\gamma$ . It does not cross-react with recombinant human, mouse, or rat IFN $\gamma$ . They do not cross-react with recombinant bovine IL-1 $\beta$ , IL-2, IL-4, IL-6, IL-8, or TNF $\alpha$ .

---

## General Reference

*Immunoassay: A Practical Guide*. Chan and Perlstein, Eds. (1987). Academic Press: New York. p.71.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

---

### Product label explanation of symbols and warnings

|   |                |   |            |   |                        |   |        |   |              |   |                              |   |   |
|---|----------------|---|------------|---|------------------------|---|--------|---|--------------|---|------------------------------|---|---|
|  | Catalog Number |  | Batch code |  | Temperature limitation |  | Use by |  | Manufacturer |  | Consult instructions for use |  | Caution, consult accompanying documents |
|---|----------------|---|------------|---|------------------------|---|--------|---|--------------|---|------------------------------|---|---|

Manufacturer's address: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.