

Bovine IL-6 ELISA Reagent Kit

ESS0029

Number

Description

ESS0029

Bovine IL-6 ELISA Reagent Kit, pre-titered coating and detection antibodies, recommended buffers and specific assay protocol optimized for the quantitative measurement of bovine IL-6 in cell culture supernatants

Kit provides sufficient reagents for approximately five 96-well plates, provided the Bovine IL-6 ELISA Reagent Kit Protocol is followed.

Kit Contents	Size	Assay Dilution
Anti-Bovine IL-6 Coating Antibody	0.625mL	1:100
Lyophilized Recombinant Bovine IL-6 Standard	5 vials	See vial label
Anti-Bovine IL-6 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 at 2-8°C.

Table of Contents

Introduction 1

ELISA Reagent Kit Buffers..... 1

Additional Materials Required..... 1

Assay Protocol 2

Performance Characteristics 3

General References 3

Introduction

The Invitrogen™ Bovine IL-6 ELISA Reagent Kit contains pre-titered coating and detection antibodies, recommended buffers and a specific assay protocol that have been optimized for the quantitative measurement of bovine IL-6 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)

For Research Use Only. Not for use in diagnostic procedures.

MAN0014275 Rev A.00 (30) Effective Date: 26 Apr 2018

Manufacturing site: Bender MedSystems GmbH | Campus Vienna Biocenter 2 1030 | Vienna, Austria | Toll Free in USA 800.955.6288
thermofisher.com

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% 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 IL-6 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 10,000pg/mL.
2. Dilute reconstituted standard 1:2 in Reagent Diluent to prepare top Standard (5000pg/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 with moderate shaking.
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 with moderate shaking.
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.

For Research Use Only. Not for use in diagnostic procedures.

MAN0014275 Rev A.00 (30) Effective Date: 26 Apr 2018

Manufacturing site: Bender MedSystems GmbH | Campus Vienna Biocenter 2 1030 | Vienna, Austria | Toll Free in USA 800.955.6288
thermofisher.com

8. Add 100µL of diluted Streptavidin-HRP reagent to each well. Cover plate with plate sealer and incubate for 30 minutes at room temperature with moderate shaking.
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 A₄₅₀ minus A₅₅₀.

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.

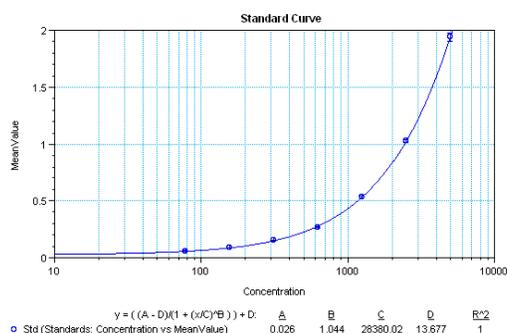
Note: When the 550nm measurement is omitted, absorbance values will be higher.

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

D. Calculation of Results

- The standard curve is used to determine bovine IL-6 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 IL-6 concentration (pg/mL) on the horizontal (X) axis.
- Calculate results using graph paper or curve-fitting statistical software. Determine the bovine IL-6 amount in each sample by interpolating from the absorbance value (Y-axis) to Bovine IL-6 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 IL-6 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 IL-6 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 IL-6. It does not cross-react with recombinant bovine IL-1β, IL-2, IL-4, IL-8, IFNγ, or TNFα (≤ 0.5%).

General Reference

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

For Research Use Only. Not for use in diagnostic procedures.

MAN0014275 Rev A.00 (30) Effective Date: 26 Apr 2018

Manufacturing site: Bender MedSystems GmbH | Campus Vienna Biocenter 2 1030 | Vienna, Austria | Toll Free in USA 800.955.6288
thermofisher.com

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

Product label explanation of symbols and warnings

 REF	Catalog Number	 LOT	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.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For Research Use Only. Not for use in diagnostic procedures.

MAN0014275 Rev A.00 (30) Effective Date: 26 Apr 2018

Manufacturing site: Bender MedSystems GmbH | Campus Vienna Biocenter 2 1030 | Vienna, Austria | Toll Free in USA 800.955.6288
thermofisher.com