

## Product datasheet

# Bovine IL-2 ELISA Kit ab193682

[2 Images](#)

### Overview

**Product name** Bovine IL-2 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

**Sample type** Cell culture supernatant, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 0.2 ng/ml

**Range** 0.2 ng/ml - 200 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	133.7	126% - 142%
Serum	194.8	125% - 145%
Plasma	102.1	90% - 111%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Cow

**Product overview**

Abcam's IL-2 Bovine ELISA Kit (ab193682) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of bovine IL-2 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for bovine IL-2 coated on a 96-well plate. Standards and

samples are pipetted into the wells and the immobilized antibody captures IL-2 present in the samples. The wells are washed and biotinylated anti-bovine IL-2 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of IL-2 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Tested applications**

**Suitable for:** Sandwich ELISA

**Platform**

Pre-coated microplate (12 x 8 well strips)

**Properties**

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated Bovine IL-2 Detection Antibody	2 vials
Bovine IL-2 Standards (lyophilized)	2 vials
Pre-coated bovine IL-2 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function**

Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

**Involvement in disease**

Note=A chromosomal aberration involving IL2 is found in a form of T-cell acute lymphoblastic leukemia (T-ALL). Translocation t(4;16)(q26;p13) with involves TNFRSF17.

**Sequence similarities**

Belongs to the IL-2 family.

**Cellular localization**

Secreted.

**Applications**

Our [Abpromise guarantee](#) covers the use of **ab193682** in the following tested applications.

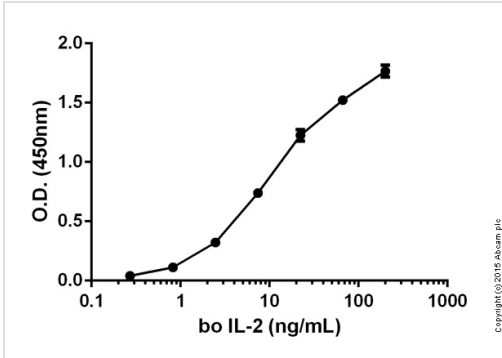
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
-------------	-----------	-------

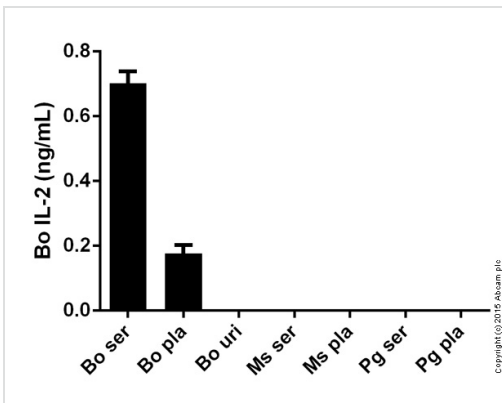
Sandwich ELISA

Use at an assay dependent concentration.

## Images



Sandwich ELISA -ab193682 Bo IL-2



Sandwich ELISA - ab193682 Bovine IL-2

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors