



EAGLE  
BIOSCIENCES

# Mouse IL-33 ELISA Assay Kit

Catalog number: ILT11-K01

For the quantitative determination of IL-33  
in mouse serum, plasma and cell culture supernate samples

This package insert must be read in its entirety before using this product  
Use only the current version of product data sheet enclosed with the kit

Catalog Number:

**ILT11-K01 (1 x 96 Wells)**

*For Research Use Only (RUO). Not for use in clinical, diagnostic or therapeutic procedures.  
v. 2.0*

---

EAGLE BIOSCIENCES, INC.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com)



## TABLE OF CONTENTS

Contents	Page
ASSAY PRINCIPLE	1
INTENDED USE	1
REAGENTS SUPPLIED	2
OTHER MATERIALS REQUIRED, BUT NOT PROVIDED	2
STORAGE	2
PREPARATION OF REAGENTS	3
PREPARATION OF STANDARDS AND SAMPLES	4
ASSAY PROCEDURE	5
CALCULATION	5
TYPICAL STANDARD CURVE	6
ASSAY CHARACTERISTICS	7
SUMMARY OF ASSAY PROCEDURE	10



## **ASSAY PRINCIPLE**

This assay is a quantitative sandwich enzyme-linked immunosorbent assay (ELISA). The microtiter plate is pre-coated with a polyclonal antibody specific for mouse IL-33. Standards and samples are pipetted into the wells and any mouse IL-33 present is bound by the immobilized antibody. After washing away any unbound substances, a biotin labelled polyclonal antibody specific for mouse IL-33 is added to the wells. After wash step to remove any unbound reagents, streptavidin-horseradish peroxidase conjugate (STP-HRP) is added. After the last wash step, an HRP substrate solution is added and color develops in proportion to the amount of mouse IL33 bound initially. The assay is stopped, and the optical density of the wells is determined using a microplate reader. Since the increases in absorbance are directly proportional to the amount of captured mouse IL-33, the unknown sample concentration can be interpolated from a reference curve included in each assay.

## **INTENDED USE**

The Eagle Biosciences Mouse IL-33 ELISA kit is designed for quantification of mouse IL-33 in serum, plasma samples and cell culture supernatant samples. The Mouse IL-33 ELISA Assay Kit is for research use only and not to be used for diagnostic purposes.



## **REAGENTS SUPPLIED**

*Each kit is sufficient for one 96-well plate and contains the following components:*

1. Microtiter Strips (96 wells), coated with a polyclonal antibody against mouse IL-33, sealed
2. 10×Wash buffer, 50 mL
3. 5×Assay buffer, 20 mL
4. 100×Detection antibody solution, a biotin labelled polyclonal antibody against mouse IL-33, 0.12 mL
5. Mouse IL-33 standard, 2000 pg of recombinant mouse IL-33, lyophilized
6. 200×STP-HRP solution, 0.06 mL
7. Substrate solution, 12 mL, ready for use
8. Stop solution, 12 mL, ready for use

## **OTHER MATERIALS REQUIRED, BUT NOT PROVIDED**

1. Pipettes and pipette tips
2. 96-well plate or manual strip washer
3. Buffer and reagent reservoirs
4. Paper towels or absorbent paper
5. Plate reader capable of reading absorbency at 450 nm
6. Distilled water or deionized water

## **STORAGE**

The kit should be stored at 2-8°C upon receipt, and all reagents should be equilibrated to room temperature before use. Remove any unused antibody-coated strips from the mouse IL-33 microtiter plate, return them to the foil pouch and re-seal. Once opened, the strips may be stored at 2-8°C for up to one month.



## PREPARATION OF REAGENTS

*Bring all reagents and materials to room temperature before assay.*

### **A. 1×Assay buffer**

Prepare 1×Assay buffer by mixing the 5×Assay buffer (20 mL) with 80 mL of distilled water or deionized water. If precipitates are observed in the 5×Assay buffer bottle, warm the bottle in a 37°C water bath until the precipitates disappear. The 1×Assay buffer may be stored at 2-8°C for up to one month.

### **B. 1×Wash buffer**

Prepare 1×Wash buffer by mixing the 10×Wash buffer (50 mL) with 450 mL of distilled water or deionized water. If precipitates are observed in the 10×Wash buffer bottle, warm the bottle in a 37°C water bath until the precipitates disappear. The 1×Wash buffer may be stored at 2-8°C for up to one month.

### **C. 1×Detection antibody solution**

Spin down the 100×Detection antibody solution briefly and dilute the desired amount of the antibody 1:100 with 1×Assay buffer, 100 µL of the 1×Detection antibody solution is required per well. Prepare only as much 1×Detection antibody solution as needed. Return the 100×Detection antibody solution to 2-8°C immediately after the necessary volume is removed.

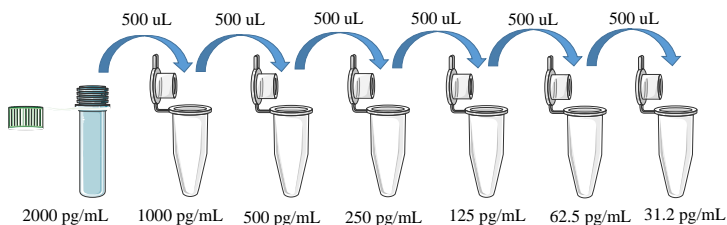
### **D. 1×STP-HRP solution**

Spin down the 200×STP-HRP solution briefly and dilute the desired amount of the 200×STP-HRP solution 1:200 with 1×Assay buffer, 100 µL of the 1×STP-HRP solution is required per well. Prepare only as much 1×STP-HRP solution as needed. Return the 200×STP-HRP solution to 2-8°C immediately after the necessary volume is removed.



## PREPARATION OF STANDARDS AND SAMPLES

**Mouse IL-33 Standards:** Reconstitute the lyophilized standard with 1 mL of 1×Assay buffer to generate a standard stock solution of 2000 pg/mL. Allow the standard to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare serially diluted standards using 1×Assay buffer as shown below.



1×Assay buffer serves as the zero standard (0 pg/mL). The reconstituted standard stock should be aliquoted and stored at -20°C for one month. Avoid repeating freezing/thawing cycles. Please do not store the diluted standard solutions.

### Sample preparation

Serum, plasma or cell culture supernate sample generally requires appropriate dilution in the 1×Assay buffer. It is recommended that the users establish their own dilution factors based on the concentration range of their samples.



## ASSAY PROCEDURE

*It is recommended that all standards and samples should be assayed in duplicate.*

1. Add 100  $\mu$ L of standard or sample per well, incubate at room temperature for 2 hours.
2. Discard the content and tap the plate on a clean paper towel to remove residual solution in each well. Add 300  $\mu$ L of 1 $\times$ Wash buffer to each well and incubate for 1 minute. Discard the 1 $\times$ Wash buffer and tap the plate on a clean paper towel to remove residual wash buffer. Repeat the wash step for a total 3 washes.
3. Add 100  $\mu$ L of 1 $\times$ Detection antibody solution to each well, incubate at room temperature for 1 hour.
4. Wash each well 3 times as in step 2.
5. Add 100  $\mu$ L of 1 $\times$ STP-HRP solution to each well, incubate at room temperature for 20 minutes.
6. Wash each well 4 times as described in step 2.
7. Add 100  $\mu$ L of Substrate solution to each well, incubate at room temperature for 15 minutes. **Protect from light.**
8. Add 100  $\mu$ L of Stop solution to each well, gently tap the plate frame for a few seconds to ensure thorough mixing.
9. Measure absorbance of each well at 450 nm immediately.

## CALCULATION

1. Subtract the absorbance of the blank from that of standards and samples.
2. Generate a standard curve by plotting the absorbance obtained (y-axis) against mouse IL-33 concentrations (x-axis). The best fit line can be generated with any curvefitting software by regression analysis. Any curve of 4-parameter or log-log curve fitting can be used for calculation.
3. Determine mouse IL-33 concentration of samples from standard curve and multiply the value by the dilution factor.

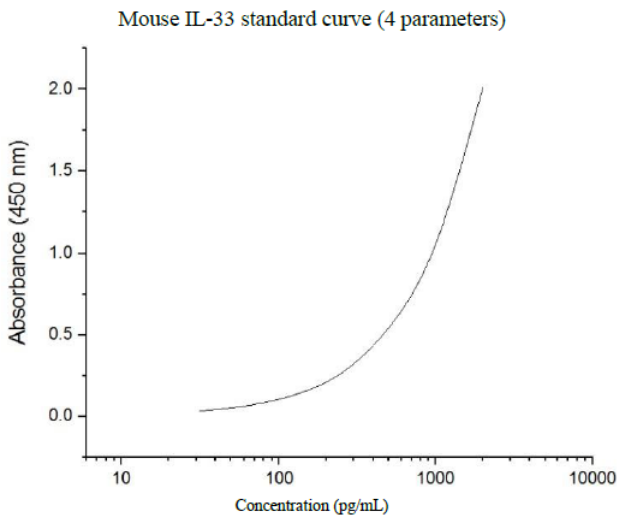


### Typical standard curve

The following standard curve is provided for demonstration only. A standard curve should be generated for each set of sample assay.

Mouse IL-33 (pg/ml)	Absorbance (450 nm)	Blanked Absorbance
0	0.11	0
31.25	0.145	0.027
62.5	0.179	0.062
125	0.236	0.118
250	0.319	0.202
500	0.65	0.532
1000	1.2	1.083
2000	2.35	2.233

Mouse IL-33 standard curve (4 parameters)





## ASSAY CHARACTERISTICS

### A. Sensitivity:

The lowest level of mouse IL-33 that can be detected by this assay is 6.5 pg/mL.

### B. Precision:

Intra-Assay Precision (Precision within an assay)

Three samples of known concentration were tested 8 times on one plate.

Sample	Mean (pg/ml)	SD (pg/ml)	CV (%)
1	1387.6	56.4	4.1
2	700.1	20.7	6.5
3	56.4	3.0	5.7

Intra-Assay Precision (Precision within an assay)

Three samples of known concentration were tested 8 times on separate plates.

Sample	Mean (pg/ml)	SD (pg/ml)	CV (%)
1	1425.7	33.5	2.4
2	713.8	29.6	4.2
3	146.3	18.3	12.5

### C. Spiking:

Cell culture supernate samples were assayed by adding 90  $\mu$ L of sample and 10  $\mu$ L of spike stock solution calculated to yield the intended 0, 150, 750 or 1500 pg/mL spike concentration.

Sample	Spiked level	Expected (pg/ml)	Observed (pg/ml)	Recovery %
Pooled cell culture medium (4X)	Low spike (150 pg/ml)	96.4	111.4	115.6
	Medium spike (750 pg/ml)	696.4	719.4	103.3
	High spike (1500 pg/ml)	1319.4	1343.4	101.8



#### D Linearity:

To assess the linearity of the assay, samples containing and/or spiked with high concentrations of mouse IL-33 were serially diluted with the 1×Assay buffer to produce samples with values within the dynamic range of the assay.

Serial dilution	Measured (pg/ml)	Expected (pg/ml)	Recovery %
Neat	1427.7	1427.7	100
1:2	703.7	713.9	98.6
1:4	329.7	356.9	92.4
1:8	171.2	178.5	95.9

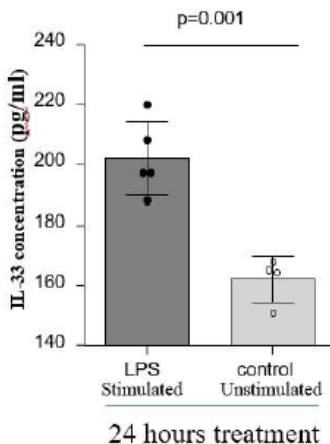
#### F. Validation:

##### Cell culture supernates:

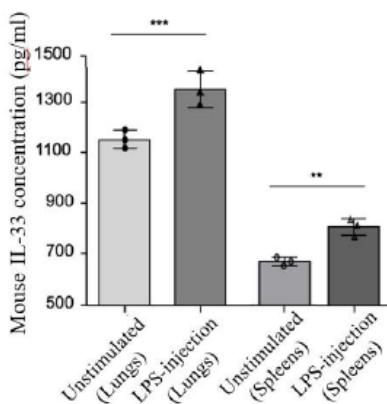
Lungs from mice were chopped into 1-2 mm pieces and cultured in 15 mL RPMI supplemented with 10% FBS, 50  $\mu$ M  $\beta$ -ME, 100 U/mL penicillin, and 100  $\mu$ g/mL streptomycin sulfate or stimulated with 1.0  $\mu$ g/mL lipopolysaccharide (LPS) for 24 hours. Cell culture medium was removed and assayed for levels of mouse IL-33. In addition, we also collected the cell lysates and measured IL-33 levels and compared with the total protein levels in the cell lysates.

Sample (cell lysates)	IL-33 levels/ total protein (pg/mg)
LPS-stimulated	381.7
Unstimulated	190.9

A per-body-weight dose (100 ug/10g) LPS were injected into mice (i.p.) and after 12 hours mice was sacrificed. Lungs and spleens from mice were chopped into 1-2 mm pieces and cultured in 6-well plate. After 24 hours incubation, cell culture medium were removed and assayed for levels of mouse IL-33.

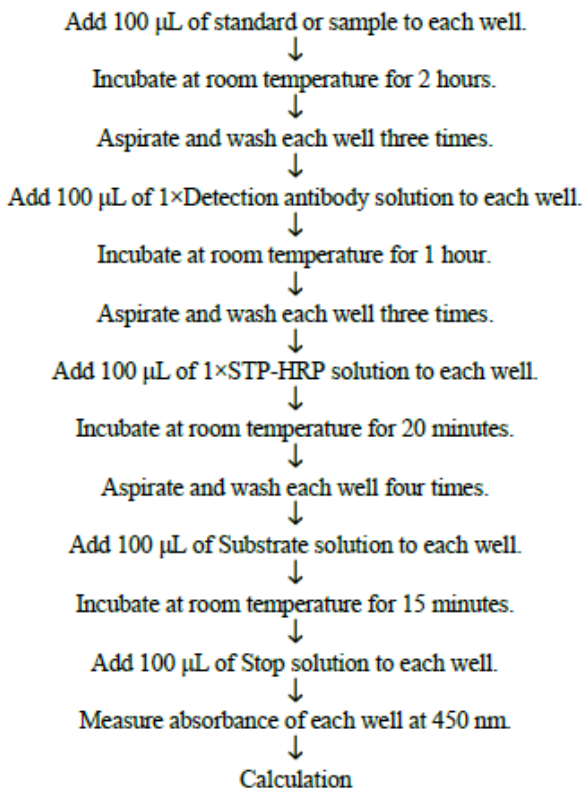


A per-body-weight dose (100 ug/10g) LPS were injected into mice (i.p.) and after 12 hours mice was sacrificed. Lungs and spleens from mice were chopped into 1-2 mm pieces and cultured in 6-well plate. After 24 hours incubation, cell culture medium were removed and assayed for levels of mouse IL-33.





### SUMMARY OF ASSAY PROCEDURE



## Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

*For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.*

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
BIOSCIENCES