



Human HMGB1 ELISA Kit

Enzyme Immunoassay kit for the quantification of Human HMGB1 in Human plasma and cell culture supernatants samples.

Catalog number: ARG81185

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 • info@eaglebio.com



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

TABLE OF CONTENTS

SECTION	Page
INTRODUCTION	3
PRINCIPLE OF THE ASSAY	3
MATERIALS PROVIDED & STORAGE INFORMATION	4
MATERIALS REQUIRED BUT NOT PROVIDED	5
TECHNICAL HINTS AND PRECAUTIONS	5
SAMPLE COLLECTION & STORAGE INFORMATION	6
REAGENT PREPARATION.....	7
ASSAY PROCEDURE	8
CALCULATION OF RESULTS	9
EXAMPLE OF TYPICAL STANDARD CURVE	10
QUALITY ASSURANCE.....	10

MANUFACTURED BY:

Arigo Biolaboratories Corporation

Address: No. 22, Ln. 227, Gongyuan Rd., Hsinchu City 300, Taiwan

Phone: +886 (3) 562 1738

Fax: +886 (3) 561 3008

Email: info@arigobio.com

INTRODUCTION

Multifunctional redox sensitive protein with various roles in different cellular compartments. In the nucleus is one of the major chromatin-associated non-histone proteins and acts as a DNA chaperone involved in replication, transcription, chromatin remodeling, V(D)J recombination, DNA repair and genome stability. Proposed to be a universal biosensor for nucleic acids. Promotes host inflammatory response to sterile and infectious signals and is involved in the coordination and integration of innate and adaptive immune responses. In the cytoplasm functions as sensor and/or chaperone for immunogenic nucleic acids implicating the activation of TLR9-mediated immune responses, and mediates autophagy. Acts as danger associated molecular pattern (DAMP) molecule that amplifies immune responses during tissue injury. Released to the extracellular environment can bind DNA, nucleosomes, IL-1 beta, CXCL12, AGER isoform 2/sRAGE, lipopolysaccharide (LPS) and lipoteichoic acid (LTA), and activates cells through engagement of multiple surface receptors. In the extracellular compartment fully reduced HMGB1 (released by necrosis) acts as a chemokine, disulfide HMGB1 (actively secreted) as a cytokine, and sulfonyl HMGB1 (released from apoptotic cells) promotes immunological tolerance. [provided by Uniprot]

PRINCIPLE OF THE ASSAY

This assay employs the sandwich enzyme immunoassay technique for the detection of Human HMGB1 in plasma and cell culture supernatants samples. An antibody specific for HMGB1 has been pre-coated onto a microtiter plate. Standards or samples are pipetted into the wells and then a Horseradish

Human HMGB1 ELISA Kit ARG81185

Peroxidase (HRP) conjugated primary antibody binds to HMGB1 is added to each well and incubates. After washing away any unbound antibody-enzyme reagent, a substrate solution (TMB) is added to the wells and color develops in proportion to the amount of total HMGB1 bound in the initial step. The color development is stopped by the addition of acid and the intensity of the color is measured at a wavelength of 450nm \pm 2nm. The concentration of total HMGB1 in the sample is then determined by comparing the O.D of samples to the standard curve.

MATERIALS PROVIDED & STORAGE INFORMATION

Store unopened kit at 2-8 °C. Use the kit before expiration date.

Component	Quantity	Storage information
antibody-coated microplate	12 X 8 strips	4°C
Standard (100 ng)	1 vial	4°C, store at -20 °C after reconstitution
Standard reconstitution buffer	1.1 ml	4°C
HRP-antibody Conjugate (200X)	115 μ l	4°C
Sample & Antibody diluent buffer	50 ml	4°C
10X Wash buffer	20 ml	4°C
TMB substrate	12 ml (ready-to-use)	4°C (Protect from light)
STOP solution	6 ml	4°C
Plate sealer	1 strips	Room temperature

MATERIALS REQUIRED BUT NOT PROVIDED

- Microplate reader capable of measuring absorbance at 450nm
- Pipettes and pipette tips
- Deionized or distilled water
- Automated microplate washer (optional)

TECHNICAL HINTS AND PRECAUTIONS

- Wear protective gloves, clothing, eye, and face protection especially while handling blood or body fluid samples.
- Store the kit at 4°C at all times.
- Briefly spin down the HRP-antibody conjugate before use.
- If crystals are observed in the 10X Wash buffer, warm to RT (not more than 50°C) until the crystals are completely dissolved. 1X Wash buffer should be prepared and stored at 4°C before use.
- Ensure complete reconstitution and dilution of reagents prior to use.
- It is highly recommended that the standards, samples and controls be assayed in duplicates.
- Change pipette tips between the addition of different reagent or samples.
- Samples contain azide cannot be assayed.
- Cell culture supernatants samples contain animal serum (E.g. fetal bovine serum (FBS) or calf serum (CS)) might contain HMGB1 from the animal serum and it might affect the accuracy of detection. Therefore, it is recommended to use BSA to replace serum in cell culture medium or chemical modified serum free medium (e.g. OPTI-MEM) for culture cells.

Human HMGB1 ELISA Kit ARG81185

- Human serum/plasma from various autoimmune diseases might contain HMGB1 autoantibodies which might result in interference of this assay.

SAMPLE COLLECTION & STORAGE INFORMATION

The sample collection and storage conditions listed below are intended as general guidelines. Sample stability has not been evaluated.

Plasma: Collect plasma using EDTA or heparin as an anticoagulant. Centrifuge for 15 minutes at 1000xg within 30 minutes of collection. It is important to ensure a platelet free preparation since platelets can release HMGB1. Assay immediately or aliquot and store at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles. Avoid using haemolytic, icteric or lipaemic samples and samples contain azide cannot be assayed.

Cell Culture Supernatants: It is recommended to use serum-free medium or use BSA to replace serum in cell culture medium or use chemical modified serum free medium for cell culture. Users should choose suitable culture medium based on the cell types. Remove particulates by centrifugation and aliquot & store samples at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles.

Note: Due to the expression/releasing levels of different cell lines or conditions are variant, samples from Cell Culture Supernatant might need to be concentrated by speedvac or protein concentrator spin columns before assay. (It is recommended to do pre-test to determine the suitable concentration factor. For our experience, we concentrated 5X for the Cell Culture Supernatant from LPS treated Huh7 cells before assay).

Human HMGB1 ELISA Kit ARG81185

REAGENT PREPARATION

- **Standard:** Reconstitute the standard with **1.0 ml** Standard reconstitution buffer to yield a stock concentration of 100 ng/ml. **Vortex** the tube for **1 min** and **stand for 10 minutes** to completely dissolve contents. Aliquot and store at $\leq -20^{\circ}\text{C}$ for long term storage.

Make sure the standard is dissolved completely before making serial dilutions. The Sample & Antibody diluent buffer serves as zero standard (0 ng/ml), and the standard stock can be diluted with Sample & Antibody diluent buffer as according to the suggested concentration below: 20 ng/ml, 10 ng/ml, 5 ng/ml, 2.5 ng/ml, 1.25 ng/ml, 0.625 ng/ml and 0.3125 ng/ml.

Dilute HMGB1 standard as according to the table below:

Standard	HMGB1 Conc. (ng/ml)	μl of Sample & Antibody diluent buffer	μl of standard
S7	20	400	100 (100 ng/ml Stock)
S6	10	250	250 (S7)
S5	5	250	250 (S6)
S4	2.5	250	250 (S5)
S3	1.25	250	250 (S4)
S2	0.625	250	250 (S3)
S1	0.3125	250	250 (S2)
S0	0	250	0

- **1X Wash buffer:** Dilute 10X wash buffer with distilled water to yield 1X wash buffer. The diluted Wash buffer should be stored at 4°C .
- **HRP-antibody Conjugate:** Diluent 200X HRP-antibody Conjugate into Sample & Antibody diluent buffer to yield 1X Detection antibody

Human HMGB1 ELISA Kit ARG81185

solution. The 1X HRP-antibody Conjugate solution should be used immediately.

- **Sample:** If the initial assay found samples contain HMGB1 higher than the highest standard, the samples can be diluted with Sample & Antibody diluent buffer and re-assay the samples. For the calculation of the concentrations this dilution factor has to be taken into account.

Example:

- a) Dilution 1:10: 10 µl sample + 90 µl Sample & Antibody diluent buffer (mix thoroughly).
- b) Dilution 1:100: 10 µl 1:10 diluted a) + 90 µl Sample & Antibody diluent buffer (mix thoroughly).

ASSAY PROCEDURE

All materials do not need equilibrated to room temperature (RT) before use. Standards, samples and controls should be assayed in duplicates.

1. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, and reseal it.
2. Add 100 µl of standards, samples and zero controls into appropriate wells.
3. Add 200 µl 1XHRP-antibody conjugate into each well. Mix thoroughly by gently shaking the plate. Cover wells and incubate at 4 °C overnight.
4. Aspirate each well and wash, repeating the process 4 times for a total 5 washes. Wash by filling each well with **1X cold wash buffer** (350 µl) using a squirt bottle, manifold dispenser, or autowasher. Complete removal of liquid at each is essential to good performance. After the last wash, remove any remaining buffer by aspirating, decanting or blotting against

Human HMGB1 ELISA Kit ARG81185

clean paper towels.

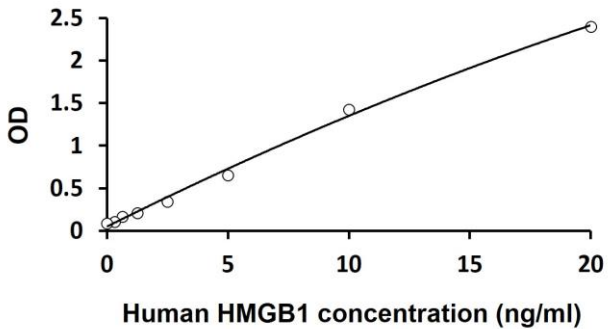
5. Add 100 μ l of TMB substrate to each well. Incubate for 10 minutes at RT in dark. Substrate will change from colorless to different strengths of blue.
6. Add 50 μ l of Stop solution to each well. The color of the solution should change from blue to yellow. Mix thoroughly by gently shaking the plate.
7. Read the OD with a microplate reader at 450 nm immediately.

CALCULATION OF RESULTS

1. Calculate the average absorbance values for each set of standards, controls and patient samples.
2. Using linear graph paper, construct a standard curve by plotting the mean absorbance obtained from each standard against its concentration with absorbance value on the vertical (Y) axis and concentration on the horizontal (X) axis.
3. Using the mean absorbance value for each sample determine the corresponding concentration from the standard curve.
4. Automated method: The results in the IFU have been calculated automatically using a 4 PL (4 Parameter Logistics) curve fit. 4 Parameter Logistics is the preferred method. Other data reduction functions may give slightly different results.
5. If the samples have been diluted or concentrated, the concentration read from the standard curve must be further converted by the appropriate dilution or concentration factor according to the sample preparation procedure as described above.

EXAMPLE OF TYPICAL STANDARD CURVE

The following data is for demonstration only and cannot be used in place of data generations at the time of assay.



QUALITY ASSURANCE

Sensitivity

Standard range: 0.3125 – 20 ng/ml

Minimum Detectable Concentration: 0.3 ng/ml

Precision:

Intra-assay: 4.3%

Inter-assay: 5.2%

Recovery:

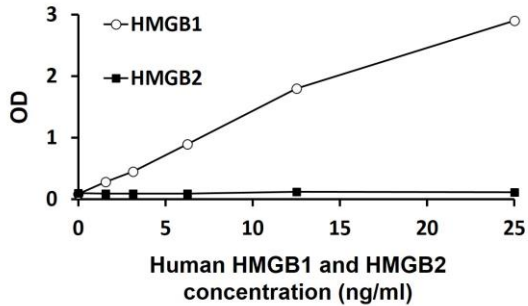
94.8-103%

Human HMGB1 ELISA Kit ARG81185

Specificity

This assay recognizes natural and recombinant total human HMGB1.

No significant cross-reactivity or interference with 25 ng/mL of recombinant human HMGB2 was observed.



The ELISA Kit might react to Mouse HMGB1, Rat HMGB1 based on sequence homology.

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 • info@eaglebio.com



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 or at 866-411-8023.

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, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.