



EAGLE
BIOSCIENCES

Human Anti-Human Antibody (HAHA) ELISA Assay Kit

Catalog Number:

HAH31-K01 (1 x 96 wells)

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

v. 1.0

EAGLE BIOSCIENCES, INC.
20A Northwest Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



INTENDED USE

Eagle Biosciences *Human Anti-Human Antibody* (HAHA) ELISA Assay Kit is intended for use in the quantitative determination of human anti-human (IgG) antibody (HAHA) levels in patient serum or plasma samples. The *HAHA* ELISA Assay Kit is for research use only and not intended for diagnostic procedures.

BACKGROUND

Clinically, humanized monoclonal antibodies (IgG) and their fragments are used in vivo diagnosis procedure (radionuclides) and treatment for patients with various diseases. In patients, even a single dose injection of a humanized monoclonal antibody or its fragment may induce immune response directed against this foreign protein (immunogen). Also, people with autoimmune diseases, such as rheumatoid arthritis, lupus, etc. produce autoantibody against human IgG. In the circulation, the presence of human antibody against human IgG would bind to the injected humanized antibody therapeutics or diagnosis and, therefore, diminish the efficacy of either in-vivo diagnosis or treatment. Especially, the HAHA would increase the risk of anaphylactic complications to subsequent administration of the humanized monoclonal antibody-based therapy.

The presence of HAHA in patient serum or plasma specimens may cause both false positive and false negative immunoassay test results depending on assay principles and monoclonal antibodies used in the assay system.

This Eagle Biosciences HAHA ELISA Assay Kit is a ready-to-use test kit with a breakable well microtiter plate and simple test procedures. It also provides a wide measurement range without high dose "hook" effect.

PRINCIPLE OF THE ASSAY

The Eagle Biosciences HAHA ELISA Assay Kit is designed, developed and produced for the quantitative measurement of HAHA in serum and plasma samples. The assay utilizes the two-site "sandwich" technique with two selected antibodies that bind to HAHA.

Assay standards, controls and patient samples are directly added to wells of a microplate that is coated with highly purified human IgG. After the first incubation period, the HAHA binds to the human IgG on the wall of microtiter well and unbound proteins in each microtiter well are washed away. Then a horseradish peroxidase (HRP)-labeled human antibody is added to each microtiter well and a "sandwich" of "well-coated human IgG – HAHA – HRP-conjugated human antibody" is formed. The unbound HRP-conjugated human antibody is removed in the subsequent washing step. For the detection of this immunocomplex, the well is then incubated with a substrate solution in a timed reaction and then measured in a spectrophotometric microplate reader. The enzymatic activity of the immunocomplex bound to HAHA on the wall of the microtiter well is directly proportional to the amount of HAHA in the sample. A standard curve is generated by plotting the absorbance versus the respective HAHA concentration for each standard on point-to-point or 4 parameter curve fit. The concentration of HAHA in test samples is determined directly from this standard curve.

REAGENTS: Preparation and Storage

The HAHA ELISA Assay Kit must be stored at 2 – 8 °C upon receipt. For the expiration date of the kit refer to the label on the kit box. All components are stable until this expiration date.



Prior to use, allow all reagents to equalize to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.

1. Human IgG Coated Microplate

One well breakable microplate with 12 x 8 strips (96 wells total) coated with highly purified human IgG. The plate is framed and sealed in a foil zipper bag with a desiccant. This reagent should be stored at 2 – 8 °C and is stable until the expiration date on the HAHA ELISA Assay Kit box.

2. HRP Conjugated Human Antibody

One vial containing **0.6 mL** HRP-labeled human antibody in a stabilized protein matrix. This reagent must be diluted with Tracer Antibody Diluent before use. This reagent should be stored at 2 – 8 °C and is stable until the expiration date on the HAHA ELISA Assay Kit box.

3. Tracer Antibody Diluent

One vial containing **12 mL** ready-to-use buffer. It should be used only for tracer antibody dilution according to the assay procedures. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the HAHA ELISA Assay Kit box.

4. Assay Buffer

One bottle containing 30 mL of ready-to-use phosphate buffered saline based assay buffer with bovine serum albumin added. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the HAHA ELISA Assay Kit box

5. ELISA Wash Concentrate

One bottle contains 30 mL of 30-fold concentrate. Before use the contents must be diluted with 870 mL of distilled water and mixed well. Upon dilution, this yields a working wash solution containing a surfactant in phosphate-buffered saline with a non-azide and non-mercury-based preservative. The diluted wash buffer should be stored at room temperature and is stable until the expiration date on the HAHA ELISA Assay Kit box.

6. ELISA HRP Substrate

One bottle contains 12 mL of tetramethylbenzidine (TMB) with stabilized hydrogen peroxide. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the HAHA ELISA Assay Kit box.

7. ELISA Stop Solution

One bottle contains 12 mL of 0.5 M sulfuric acid. This reagent should be stored at 2 – 8°C or room temperature and is stable until the expiration date on HAHA ELISA Assay Kit box.

8. HAHA Standards

Five vials containing different levels of HAHA in a liquid protein matrix with a non-azide based preservative. **Refer to vials for exact concentration for each standard.** These reagents should be stored at 2 – 8°C and are stable until the expiration date on the HAHA ELISA Assay Kit box



9. HAHA Controls

Two vials containing different levels of HAHA in a liquid protein matrix with a non-azide based preservative. **Refer to vials for exact concentration range for each control.** Both controls should be stored at 2 – 8°C and are stable until the expiration date on the HAHA ELISA Assay Kit box

SAFETY PRECAUTIONS

The reagents of the HAHA ELISA Assay Kit must be used in a professional laboratory environment and are for research use only. Source material (e.g. highly purified bovine serum albumin) of bovine serum was derived in the contiguous 48 United States. It was obtained only from healthy donor animals maintained under veterinary supervision and found free of contagious diseases. Wear gloves while performing this assay and handle these reagents as if they are potentially infectious. Avoid contact with reagents containing TMB, hydrogen peroxide, or sulfuric acid. TMB may cause irritation to skin and mucous membranes and cause an allergic skin reaction. TMB is a suspected carcinogen. Sulfuric acid may cause severe irritation on contact with skin. Do not get in eyes, on skin, or on clothing. Do not ingest or inhale fumes. On contact, flush with copious amounts of water for at least 15 minutes. Use Good Laboratory Practices.

MATERIALS REQUIRED BUT NOT PROVIDED

- Precision single channel pipettes capable of delivering 25 µL, 50 µL, 100 µL, and 1000 µL etc.
- Repeating dispenser suitable for delivering 100 µL.
- Disposable pipette tips suitable for above volume dispensing.
- Disposable 12 x 75 mm or 13 x 100 glass tubes.
- Disposable plastic 100 mL and 1000 mL bottle with caps.
- Aluminum foil.
- Deionized or distilled water.
- Plastic microtiter well cover or polyethylene film.
- ELISA multichannel wash bottle or automatic (semi-automatic) washing system.
- Spectrophotometric microplate reader capable of reading absorbance at 450 nm.

SPECIMEN COLLECTION

Only 50 µL of human serum or plasma is required for HAHA measurement in duplicate. No special preparation of individual is necessary prior to specimen collection. In the case of serum, whole blood should be collected and must be allowed to clot for a minimum of 30 minutes at room temperature before the serum is separated by centrifugation (850 – 1500xg for 10 minutes). The serum should be separated from the clot within three hours of blood collection and transferred to a clean test tube. Serum or plasma samples should be stored at 2 - 8°C if the assay is to be performed within 72 hours. Otherwise, patient samples should be stored at -20°C or below until measurement. Avoid repeated (more than three times) freezing and thawing of specimen.

ASSAY PROCEDURE

1. Reagent Preparation



- (1) Prior to use allow all reagents of the HAHA ELISA Assay Kit to come to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.
- (2) ELISA Wash Concentrate must be diluted to working solution prior to use. Please see REAGENTS section for details)

2. Assay Procedure

1. Place a sufficient number of human IgG-coated microwell strips/wells in a holder to run HAHA.
2. Test Configuration

ROW	STRIP 1	STRIP 2	STRIP 3
A	STD 1	STD 5	SAMPLE 2
B	STD 1	STD 5	SAMPLE 2
C	STD 2	C 1	SAMPLE 3
D	STD 2	C 1	SAMPLE 3
E	STD 3	C 2	SAMPLE 4
F	STD 3	C 2	SAMPLE 4
G	STD 4	SAMPLE 1	
H	STD 4	SAMPLE 1	

3. Add **25 µL** of standards, controls and patient samples into the designated microwell.
4. Add **100 µL** of assay buffer to each well.
5. Cover the plate with one plate sealer and incubate plate at room temperature, shaking for **45 minutes**.
6. Prepare HAHA Tracer antibody working solution by **1:21** fold dilution of the HRP-conjugated human antibody with the tracer Antibody Diluent. For each strip, it is required to mix 1 mL of the tracer antibody diluent with 50 µL of the tracer antibody in a clean test tube.
7. Remove plate sealer. Aspirate the contents of each well. Wash each well 5 times by dispensing 350 µL of working wash solution into each well and then completely aspirating the contents. Alternatively, an automated microplate washer can be used.
8. Add **100 µL** of above diluted HRP-conjugated Human Antibody working solution to each of the wells.
9. Cover the plate with a plate sealer and an aluminum foil and incubate plate at room temperature, shaking for **45 minutes**.
10. Remove plate sealer. Aspirate the contents of each well. Wash each well 5 times by dispensing 350 µL of working wash solution into each well and then completely aspirating the contents. Alternatively, an automated microplate washer can be used.
11. Add **100 µL** of ELISA HRP Substrate into each of the wells.
12. Cover the plate with a plate sealer and also with an aluminum foil to avoid exposure to light.
13. Incubate plate at room temperature for **20 minutes**.
14. Remove the aluminum foil and plate sealer. Add 100 µL of ELISA Stop Solution into each of the wells. Mix gently.
15. Read the absorbance at **450 nm** within 10 minutes in a microplate reader.



NOTE: to reduce the background, one can set the instrument to dual wavelength measurement at 450 nm with background wavelength correction set at 595 nm, 620 nm or 630 nm.

PROCEDURAL NOTES

1. It is recommended that all controls and unknown samples be assayed in duplicate in the HAHA ELISA Assay Kit. The average absorbance reading of each duplicate should be used for data reduction and the calculation of results.
2. For patient samples with concentration higher than level 5 standard, it is recommended to dilute the specimen with assay buffer at 1:10, 1:100, etc. for a more accurate report.
3. Keep light-sensitive reagents in the original amber bottles.
4. Store any unused human IgG-coated strips in the foil Ziploc bag with desiccant to protect from moisture.
5. Careful technique and use of properly calibrated pipetting devices are necessary to ensure reproducibility of the test.
6. Incubation times or temperatures other than those stated in this insert may affect the results.
7. Avoid air bubbles in the microwell as this could result in lower binding efficiency and higher CV% of duplicate reading.
8. All reagents should be mixed gently and thoroughly prior to use. Avoid foaming.

INTERPRETATION OF RESULTS

1. Calculate the average absorbance for each pair of duplicate test results.
2. Subtract the average absorbance of the STD 1 (0 $\mu\text{g/mL}$) from the average absorbance of all other readings to obtain corrected absorbance.
3. The standard curve is generated by the corrected absorbances of all standard levels on the ordinate against the standard concentration on the abscissa using point-to-point or log-log paper. Appropriate computer assisted data reduction programs may also be used for the calculation of results. We recommend using **4-parameter** or **point-to-point** curve fit.

The HAHA concentrations for the controls and patient samples are read directly from the standard curve using their respective corrected absorbance.

EXAMPLE DATA AND STANDARD CURVE

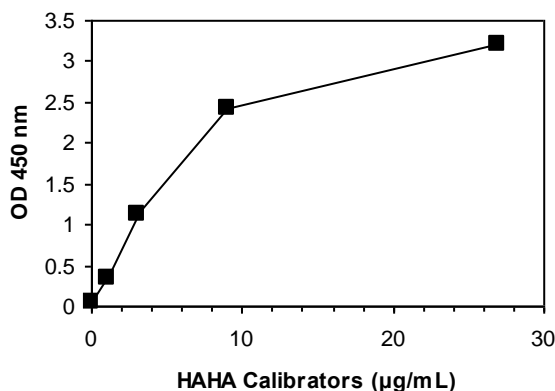
A typical absorbance data and the resulting standard curve from this HAHA ELISA are represented. **This curve should not be used in lieu of standard curve run with each assay.**

Well I.D.	OD 450 nm Absorbance			Results $\mu\text{g/mL}$
	Readings	Average	Corrected	
0 $\mu\text{g/mL}$	0.047 0.046	0.047	0.000	
1 $\mu\text{g/mL}$	0.342 0.338	0.340	0.293	
3	1.178	1.136	1.089	



µg/mL	1.094			
9	2.501	2.418	2.371	
µg/mL	2.336			
27	3.224	3.214	3.167	
µg/mL	3.205			
Control 1	0.598	0.620	0.573	1.71
	0.643			µg/mL
Control 2	1.988	2.000	1.953	7.04
	2.013			µg/mL

HAHA ELISA



EXPECTED VALUES

Eighty-six normal adult sera were measured with this HAHA ELISA Assay Kit. One hundred sixty sera showed the OD reading very close to the zero calibrator. The 99% confidence normal cut-off is **0.5 µg/mL**.

It is highly recommended that each laboratory establishes its own normal cut off level.

LIMITATION OF THE PROCEDURE

1. Since this is the first commercial assay of this kind and there is no Gold Standard concentration or international standard available for HAHA measurement, the values of assay standards were established and validated by Epitope Diagnostics. Results obtained with different assay methods or kits cannot be used interchangeably.
2. For unknown sample value read directly from the assay that is greater than the highest assay standard, it is recommended to measure a further diluted sample for more accurate measurement.
3. Bacterial or fungal contamination of serum specimens or reagents, or cross-contamination between reagents may cause erroneous results.
4. Water deionized with polyester resins may inactivate the horseradish peroxidase enzyme.

QUALITY CONTROL

To assure the validity of the results each assay should include adequate controls with known HAHA levels. We recommend that all assays include the laboratory's own HAHA controls in addition to those provided with this kit.



PERFORMANCE CHARACTERISTICS

Sensitivity

The sensitivity of this HAHA ELISA as determined by the 95% confidence limit on 20 duplicate determination of zero standard is about 0.1 µg/mL.

High Dose “hook” effect

This assay has showed that it did not have any high dose “hook” effect up to 3,000 µg/mL .

Precision

The intra-assay precision was validated by measuring one control sample in a single assay with eight replicate determinations.

Mean HAHA Value (µg/mL)	CV (%)
1.62	5.1

The inter-assay precision is validated by measuring one control sample in duplicate in 6 individual assays.

Mean HAHA Value (µg/mL)	CV (%)
6.93	6.7

REFERENCES

1. Szolar OH, Stranner S, Zinoecker I, Mudde GC, Himmler G, Waxenecker G, Nechansky A. Qualification and application of a surface plasmon resonance-based assay for monitoring potential HAHA responses induced after passive administration of a humanized anti Lewis-Y antibody. *J Pharm Biomed Anal.* 2006 Jun 16;41(4):1347-53. Epub 2006 Apr 27
2. Ritter G, Cohen LS, Williams C Jr, Richards EC, Old LJ, Welt S. Serological analysis of human anti-human antibody responses in colon cancer patients treated with repeated doses of humanized monoclonal antibody A33. *Cancer Res.* 2001 Sep 15;61(18):6851-9
3. Lofgren JA, Dhandapani S, Pennucci JJ, Abbott CM, Mytych DT, Kaliyaperumal A, Swanson SJ, Mullenix MC. Comparing ELISA and surface plasmon resonance for assessing clinical immunogenicity of panitumumab. *J Immunol.* 2007 Jun 1;178(11):7467-72



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 617-419-2019.