



EAGLE  
BIOSCIENCES

# **Glutamic Acid Decarboxylase ( GAD ) ELISA Assay Kit**

Catalog Number: GDA31-K01 ( 1x 96 Wells )

For Research Use Only

*v. 1.0 (effective 06 SEP 2023)*

---

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)



## **INTENDED USE**

The Eagle Biosciences GAD autoantibody ELISA kit is a highly sensitive and specific ELISA kit for precision detection and quantitative measurement of GAD autoantibody (GADA) titres in human serum (plasma samples are not recommended). This kit is for research use only and not for use in diagnostic procedures.

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

## **ASSAY BACKGROUND**

Glutamic acid decarboxylase (GAD) autoantibodies are found in 70% to 80% of individuals with new-onset type 1 diabetes, making it the most frequent autoantibody in autoimmune diabetes. GAD autoantibodies can be detected in serum for many years post diagnosis, and high concentrations of GAD autoantibodies have been considered as a marker of faster  $\beta$ -cell exhaustion in these patients. Furthermore, GAD autoantibodies in non-diabetic individuals predicts the later development of type 1 diabetes. Besides autoimmune diabetes, GAD autoantibodies also exist in Stiff Man Syndrome, autoimmune poly-endocrinopathies, and some of Grave's Disease patients.

## **ASSAY PRINCIPLES**

In this GADA ELISA, recombinant GAD protein coated onto plate wells can specifically recognize the GAD autoantibodies in human sera and calibrators. After a 1-hour incubation, GAD autoantibodies are captured by immobilized GAD protein while the unbound components were discarded and washed away. Afterwards, biotinylated GAD protein (GAD Biotin) is added for another round of incubation for 1 hour, wherein the GAD-Biotin detects GAD autoantibodies previously bound to GAD protein on the plate. After removal of nonspecific bindings, bound GAD-Biotin is revealed by addition of streptavidin horseradish peroxidase (STV-HRP), which specifically binds with biotin, followed by the substrate 3,3',5,5'-Tetramethylbenzidine (TMB), which results in formation of a blue color. Color reaction will be further stopped by 2M H<sub>2</sub>SO<sub>4</sub>, transforming the blue color to yellow signals. The absorbance of yellow reaction mixture is measured by plate reader at 450nm and 405nm. The higher the reading is, the higher concentration of GAD autoantibodies. Low concentration of GAD autoantibodies (<18 u/ml) is recommended to be read off the 450nm calibration curve, while high value of GAD



autoantibodies to be read off 405nm standard curve. The measuring interval is 5-2000U/ml (units are NIBSC 97/550)

### **STORAGE AND PREPARATION OF TEST SERUM SAMPLES**

Test samples are suggested to be assayed immediately after separation of serum, or preferably stored at -200C or below in aliquots. Duplicate test is recommended therefore 50µl is sufficient for each aliquot (25µl per test). Lipemic or hemolyzed sera, as well as plasma samples are not recommended. When required, vortex test serum samples at room temperature to ensure homogeneity. Then centrifuge samples at 10-15,000 rpm for 5 minutes prior to assay to remove particulates. Please do not omit this centrifugation step if samples are cloudy and containing particles.

### **MATERIALS NEEDED BUT NOT SUPPLIED**

1. Pipettes capable to dispense 1000 µl, 100 µl, and 25 µl
2. Multichannel pipettes
3. Plate shake capable of shaking at 500 rpm.
4. 96-well plate reader capable of absorbance measurement at 450 nm
5. 405 nm Distilled water

### **REAGENTS PROVIDED AND REAGENT PREPARATION**

A	GAD protein coated ELISA plate	12 strips of 8 wells (96 wells in total) in a frame and sealed in a foil bag. Make sure test strips are firmly fitted into the provided frame before use. Equilibrate test strips to room temperature before use. After opening, seal unused strips in the original self-seal foil bag with desiccant. Store the re-sealed foil bag at 2-8C for up to 16 weeks.
B 1-7	Calibrators	1ml x 6 5, 10, 18, 35, 120, 250, 2000 (U/ml) (Units are WHO standard NIBSC 97/550) Ready to use
C	Positive control	1 ml x 1



		Ready to use
D	Negative control	1 ml x 1 Ready to use
E	GAD-Biotin	3 vials Lyophilized
F	Reconstitution buffer	15ml x 1 Pale yellow Ready-to-use for reconstitution of GAD Biotin.
G	20 x Streptavidin horseradish peroxidase (STV-HRP)	1ml x 1 Dilute 1 in 20 with dilution buffer ( H ). For example, 0.5 mL ( G ) + 9.5 mL ( H ).
H	Dilution buffer ( For STV-HRP )	15ml x 1 Ready to use
I	Substrate solution ( 3,3',5,5'-Tetramethylbenzidine, TMB )	Ready to use Equilibrate to room temperature for 15 minutes before use
J	Stop solution	15ml x 1 2M H <sub>2</sub> SO <sub>4</sub> Ready to use
K	10X Wash buffer	50ml x 1 Dilute with distilled water ( 1:10 )

## ASSAY PROCEDURE

### PREPARATION OF REAGENTS

1. 1X 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.

2. 1x detection GAD-Biotin solution

Reconstitute each vial of lyophilized GAD-Biotin with 4.5ml IMD reconstitution buffer (F). Return the reconstituted detection solution to 2-8°C immediately. The detection solution can be stored at 40C for up to 3 days after reconstitution or stored at -800C or -200C for long-term use in aliquots (< 3 freeze/thaw cycles).

3. 1x STV-HRP solution

Dilute 1 in 20 with dilution buffer (H). For example, 0.5 mL (G) + 9.5 mL (H). The 1×STV-HRP buffer may be stored at 2-8°C for up to one month.

**Please pre-balance all the reagents to room temperature (20-25°) for at least 30 minutes before use.**

<b>Step 1</b>	Pipette 25µl of test serum samples, calibrators (B1-7) and controls (C and D) into respective plate wells in duplicate. Leave at least one well for blank.
<b>Step 2</b>	Cover the frame and shake the wells at room temperature at 500 rpm for 1 hour on a plate shaker.
<b>Step 3</b>	Discard the content and tap the plate on a clean paper towel to remove residual solution in each well. Add 300 µl of 1×Wash buffer to each well and incubate for 1 minute. Discard the 1×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.
<b>Step 4</b>	Add 100 µl of reconstituted GAD-Biotin to each well.
<b>Step 5</b>	Cover the frame and shake the wells at room temperature at 500 rpm for 1 hour on a plate shaker.
<b>Step 6</b>	Wash each well 3 times as described in step 3.
<b>Step 7</b>	Pipette 100 µl of 1x STV-HRP solution to each well
<b>Step 8</b>	Cover the frame and shake the wells at room temperature at 500 rpm for 20



	minutes on a plate shaker.
<b>Step 9</b>	Wash each well 4 times as described in step 3.
<b>Step 10</b>	Add 100 $\mu$ l of Substrate solution (I) to each well, incubate at room temperature for 15 minutes. <b>Protect from light.</b>
<b>Step 11</b>	Add 100 $\mu$ l of Stop solution (J) to each well, gently tap the plate frame for a few seconds to ensure thorough mixing.
<b>Step 12</b>	Measure absorbance of each well at 450 nm and/or 405 nm immediately.

## DATA ANALYSIS

1. Subtract absorbance of blank from that of standards and samples.
2. A calibration curve can be established by plotting calibrator concentration on x axis (log-scale) against the absorbance of the calibrators on the y-axis (linear scale). The best fit line can be generated with any curve-fitting software by regression analysis. Any curve of 4-parameter or log-lin curve fitting can be used for calculation.
3. The GAD autoantibody concentrations in sera can be read off the calibration curve. Negative control can be assigned a value of 0.5U/ml to assist the statistical software to process the data analysis. Samples with high GAD autoantibody concentrations can be diluted with GAD autoantibody negative serum or the kit negative control. Some sera will not dilute in a linear way according to the kit calibrators (standardized against NIBSC 97/550). Most test sera will have values below 250 u/mL and the 2000 u/mL calibrator need not always be included.

## TYPICAL RESULTS

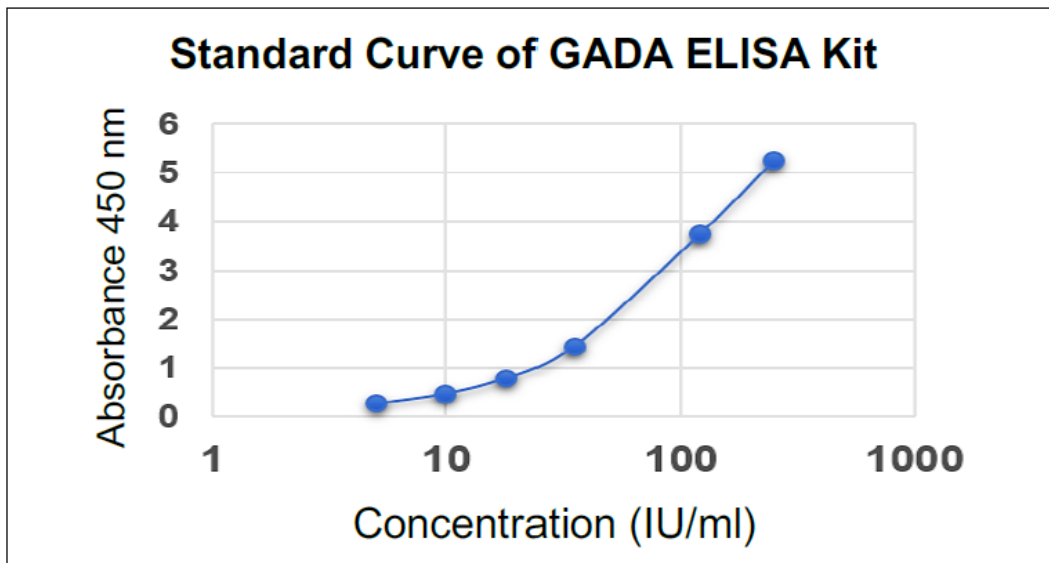
*Example only, not for calculation of actual results.*

Calibrator *	Concentration (IU/ml)	OD450	OD405
1	2000	7.86	2.3425
2	250	5.24	1.56
3	120	3.732	1.0865



4	35	1.445	0.439
5	18	0.794	0.2565
6	10	0.4785	0.168
7	5	0.276	0.1125

<b>Negative Control</b>	0	0.096	0.0655
<b>Positive Control</b>	15.18	0.746	0.235



#### ASSAY CUT-OFF VALUE

< 5 IU/ml	Negative
≥ 5IU/ml	Positive

This cut-off value has been validated at IMD. However, it is recommended that each laboratory should establish its own normal and pathological reference range for GAD autoantibody level.



Furthermore, it is also recommended that each laboratory should include its own panel of control samples in the assay.

## CLINICAL EVALUATION

Clinical Evaluation		
Sensitivity	92.5% (n=80)	
Specificity	95% (n=150)	
Inter Assay Precision		
Sample	IU/ml (n=20)	CV
1	120	5.75%
2	18	6.66%
3	5	5.26%
Intra Assay Precision		
Sample	IU/ml (n=20)	CV
1	120	6.61%
2	18	4.63%
3	5	5.65%

## REFERENCE

1. Stenström, Gunnar, et al. "Latent autoimmune diabetes in adults: definition, prevalence,  $\beta$ -cell function, and treatment." *Diabetes* 54.suppl 2 (2005): S68-S72
2. Schmidli, Robert S., Peter G. Colman, and Ezio Bonifacio. "Disease sensitivity and specificity of 52 assays for glutamic acid decarboxylase antibodies: the Second International GADAB Workshop." *Diabetes* 44.6 (1995): 636-640.
3. Moffett, Ashley, et al. "Variation of maternal KIR and fetal HLA-C genes in reproductive failure: too early for clinical intervention." *Reproductive biomedicine online* 33.6 (2016):





## **PRECAUTIONS**

- This kit is for research use only
- Compare contents and packing list, if there is breakage or shortage, notify Eagle Biosciences immediately
- Do not pipette reagents by mouth
- Do not smoke, eat or drink while performing assay
- Wear disposable gloves and proper lab protection and attire
- Treat all samples as potentially infectious
- Do not mix reagents from other lots
- Avoid contact with TMB and Stop solutions. If contact occurs, rinse thoroughly with water
- Eagle Biosciences is not responsible for outcomes as results of tampering with the reagents or using them unconventionally

## **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](mailto:info@eaglebio.com) or at 866-411-8023.