



Therapeutic IgG1 ELISA Assay Kit

Catalog Number:

IG131-K01 (1 x 96 wells)

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

v. 220106 (2.16.23)

Eagle Biosciences, Inc.

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

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

www.EagleBio.com



INTENDED USE

The Eagle Biosciences Therapeutic IgG1 ELISA Assay Kit (enzyme-linked immunoassay kit) is intended for the determination of IgG1 in serum and plasma. The Eagle Biosciences Therapeutic IgG1 ELISA Assay Kit is for research use only and not to be used in diagnostic procedures.

INTRODUCTION

Recombinant human monoclonal antibodies have become a major focus of new drug discovery in the Pharmaceutical and Biotech industries, with hundreds of new human monoclonal antibodies currently in preclinical development. Fully human recombinant antibodies have gradually replaced humanized mouse or rat antibodies, with IgG1 and IgG4 as the preferred engineered isotypes¹. As part of the developmental process, human therapeutic antibody drug candidate molecules must be tested in non-human primate species to assess their pharmacokinetic profile and potential toxicity²³. Primate and human antibodies are highly homologous, making it difficult to distinguish the human antibody drug candidate from the endogenous primate antibodies in a serum or plasma sample by traditional human IgG-specific immunoassays. Therefore, we developed two highly sensitive ELISAs that distinguish human IgG1 or human IgG4 from rhesus or cynomolgus monkey immunoglobulins. This Recombinant (Therapeutic) Human IgG1 ELISA can be used to accurately quantify human IgG1 in the serum or plasma of rhesus or cynomolgus monkeys without cross reacting with antibodies of those species. This same ELISA can also be used to quantify human IgG1 in the plasma of other preclinical test species (mouse, rat, rabbit, guinea pig, etc.) if desired. This assay is supplied with a recombinant human IgG1 kappa allotype G1m17.1 standard. The customer may wish to substitute their own specific drug candidate IgG1 for this standard in the assay.

PRINCIPLE OF THE ASSAY

This is a sandwich enzyme-linked immunosorbent assay (ELISA). The plate is pre-coated with anti-Human IgG Antibody and blocked, ready for the addition of samples and standards. The assay should take approximately 3 hours to run, plus any required sample preparation time.

MATERIALS PROVIDED

Component	Description	Volume	Storage
Anti-Human IgG Plate	96-well plate coated and blocked	1 plate	4°C
Assay Buffer	Buffer to dilute kit components and samples	100 mL	4°C
10x Wash Buffer	Buffer used to wash the plate	30 mL	4°C
Human IgG1 Standard	Lyophilized Recombinant Human IgG1	0.2 µg	4°C
10x Detection Antibody	Anti-Human IgG ¹ -HRP Conjugate	1.5 mL	4°C
TMB Substrate	Stabilized TMB color reagent	1.5 mL	4°C

MATERIALS NEEDED BUT NOT PROVIDED

1. Microplate Reader with 450 nm filter
2. Adjustable micropipettes (10 - 1000 µL) and tips
3. 3N Sulfuric Acid (H₂SO₄)
4. Plate shaker
5. Plate cover or plastic film

STORAGE

1. Store the components of this kit at the temperature specified on the labels.
2. Unopened reagents are stable until the indicated kit expiration date.



3. Reconstituted Standard may be used up to 1 month after reconstitution if stored at 4°C.
4. Diluted Detection Antibody must be used within 2 hours of dilution.

WARNINGS AND PRECAUTIONS

1. Use aseptic technique when opening and dispensing reagents.
2. This kit is designed to work properly as provided and instructed. Additions, deletions or substitutions to the procedure or reagents are not recommended, as they may be detrimental to the assay.

PROCEDURAL NOTES

1. To minimize errors in absorbance measurements due to handling, wipe the exterior bottoms of the microplate wells with a lint-free paper towel prior to inserting into the plate reader.

SAMPLE PREPARATION

If assessing human IgG1 in monkey plasma or serum, dilute the serum at least 1:5 into Assay Buffer before testing. Undiluted monkey plasma or serum contains contaminants that interfere with the accurate functioning of this ELISA.

REAGENT PREPARATION

1. **Human IgG1 Standard:** Resuspend lyophilized standard vial by adding 0.2mL of Assay Buffer directly to vial, gently invert to suspend completely, to create 1µg/mL stock solution.
2. **10x Wash Buffer:** Dilute the wash buffer 1:10 by adding 30mL of 10x Wash Buffer to 270mL of dH₂O.
3. **10x Detection Antibody:** Immediately prior to use, dilute detection antibody 1:10 by adding 1.2mL 10x Detection Antibody to 10.8mL Assay Buffer. Use within 2 hours of diluting.

STANDAR CURVE PREPARATION

Set up for the standard curve by labeling dilution tubes and dispensing the indicated volumes of Assay Buffer and 1µg/mL Standard Stock Solution according to Table 1 below.

Table 1: Standard Curve Preparation

Standard	IgG1 Concentration (ng/mL)	Assay Buffer (µL)	Transfer Volume (µL)	Transfer Source	Final Volume (µL)
S ₇	20	980	20	1µg/mL Stock	500
S ₆	10	500	500	S ₇	500
S ₅	5	500	500	S ₆	500
S ₄	2.5	500	500	S ₅	500
S ₃	1.25	500	500	S ₄	500
S ₂	0.63	500	500	S ₃	500
S ₁	0.31	500	500	S ₂	1000
S ₀	0	1000	-	-	1000



ASSAY PROCEDURE

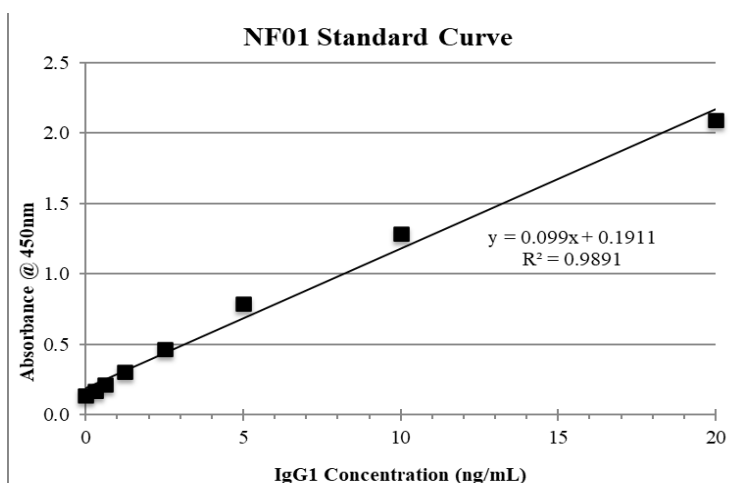
1. Add 100 μ L of Standards and Samples to the corresponding wells on the microplate in duplicate. Incubate at room temperature for 1.5 hours on an orbital plate shaker. See Scheme 1 below for a suggested plate layout.
2. Dump the contents of the plate and wash each well three times with 300 μ L of Wash Buffer. After the final wash, tap the plate upside-down on a lint-free paper towel to make sure there is no solution left in the wells.
3. Add 100 μ L of the Detection Antibody to each well. Incubate at room temperature for 1 hour on orbital plate shaker.
4. Wash the plate as in step 2.
5. Add 100 μ L of TMB Substrate to each well. Allow the color to develop for 10 minutes at room temperature.
6. Stop the reaction by adding 100 μ L per well of 3N Sulfuric Acid (H₂SO₄).
7. Read the plate at 450 nm in a microplate reader within 15 minutes

1	2	3	4	5	6	7	8	9	10	11	12
S0	S0	U1	U1	U9	U9	U17	U17	U25	U25	U33	U33
S1	S1	U2	U2	U10	U10	U18	U18	U26	U26	U34	U34
S2	S2	U3	U3	U11	U11	U19	U19	U27	U27	U35	U35
S3	S3	U4	U4	U12	U12	U20	U20	U28	U28	U36	U36
S4	S4	U5	U5	U13	U13	U21	U21	U29	U29	U37	U37
S5	S5	U6	U6	U14	U14	U22	U22	U30	U30	U38	U38
S6	S6	U7	U7	U15	U15	U23	U23	U31	U31	U39	U39
S7	S7	U8	U8	U16	U16	U24	U24	U32	U32	U40	U40

CALCULATIONS

- a. Average all duplicate well absorbance values.
- b. Plot a standard curve using the absorbance values of each Standard (y-axis) versus the Standard concentration (x-axis).
- c. Determine the concentration of unknown using the equation of the line.

Figure 1: Typical Standard Curve





REFERENCES

1. Buntz, B. (2021, May 14). 50 of 2020's best-selling pharmaceuticals. Drug Discovery and Development. Retrieved January 20, 2022, from <https://www.drugdiscoverytrends.com/50-of-2020s-best-sellingpharmaceuticals/>
2. Salfeld, J. (2007) Isotype selection in antibody engineering. Nat Biotechnol 25, 1369–1372.
3. Stubenrauch K, Wessels U, Lenz H.; (2009) Evaluation of an immunoassay for human-specific quantitation of therapeutic antibodies in serum samples from non-human primates. J Pharm Biomed Anal.; 49(4):1003-8

DISCLAIMER

This information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. Eagle Biosciences, Inc. shall not be held liable for any damage resulting from handling or from contact with the above product. See catalog for additional terms and conditions of sale.



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 or at 866-411-8023.