



**EAGLE**  
BIOSCIENCES

# **Ustekinumab mAb-based ELISA Assay Kit**

Catalog Number:

**IG-AB121 (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](http://www.EagleBio.com)



## Introduction

The drug Ustekinumab (trade name Stelara®) is a humanized immunoglobulin G1 · monoclonal antibody that binds with high specificity to the p40 protein subunit used by both the IL-12 and IL-23 cytokines.

The *ImmunoGuide* Ustekinumab ELISA (mAb-based) kit can be efficiently used for measuring free Ustekinumab pegol levels in serum and plasma.

## Intended Use

Enzyme immunoassay for the specific and quantitative determination of free Ustekinumab in serum and plasma.

## Assay Principle

This ELISA is based on Ustekinumab-specific monoclonal antibody (catcher Ab, *ImmunoGuide* clone 9C7). Diluted standards and samples are incubated in the microtiter plate coated with IG-9C7 mAb. After incubation, the wells are washed. A biotinylated anti-human IgG monoclonal antibody (clone IG-1B5; specific for the Fc part of all human IgG i.e. IgG1, IgG2, IgG3 and IgG4) is added and binds to the Fc part of Ustekinumab. Following incubation, wells are washed and the horseradish peroxidase (HRP)-conjugated streptavidin is added and binds to the biotinylated 1B5 mAb. Following incubation, wells are washed and the bound enzymatic activity is detected by addition of chromogen-substrate. The color developed is proportional to the amount of Ustekinumab in the sample or standard. Results of samples can be determined by using the standard curve. Preincubation of Ustekinumab with recombinant human interleukin-12 (IL-12), contains p40 protein subunit as that of IL-23, inhibited the reaction. Therefore, the *ImmunoGuide* Ustekinumab ELISA (mAb-based) measures the free form of Ustekinumab

## Warnings and Precautions

1. Before starting the assay, read the instructions completely and carefully. Use the valid version of the package insert provided with the kit. Be sure that everything is understood. For further information (clinical background, test performance, automation protocols, alternative applications, literature, etc.) please refer to the local distributor.
2. In case of severe damage of the kit package, please contact *Eagle Biosciences* or your supplier in writing, latest one week after receiving the kit. Do not use damaged components in test runs but keep safe for complaint related issues.
3. Obey lot number and expiry date. Do not mix reagents of different lots. Do not use expired reagents.
4. Follow good laboratory practice and safety guidelines. Wear lab coats, disposable latex gloves and protective glasses where necessary.
5. Reagents of this kit containing hazardous material may cause eye and skin irritations. See MATERIALS SUPPLIED and labels for details.
6. Chemicals and prepared or used reagents have to be treated as hazardous waste according to the national biohazard safety guidelines or regulations.
7. Avoid contact with Stop solution. It may cause skin irritations and burns.
8. If any component of this kit contains human serum or plasma it is indicated and if so, it has been tested and were found to be negative for HIV I/II, HBsAg and HCV. However, the presence of these or other infectious agents cannot be excluded absolutely and therefore reagents should be treated as potential biohazards in use and for disposal.
9. Some reagents contain preservatives. In case of contact with eyes or skin, flush immediately with water.



### Storage and Stability

The kit is shipped at ambient temperature and should be stored at 2-8°C. Keep away from heat or direct sun light. The storage and stability of specimen and prepared reagents is stated in the corresponding chapters. The microtiter strips are stable up to the expiry date of the kit in the broken, but tightly closed bag when stored at 2-8°C.

### Kit Components/Materials Provided

Quantity	Component
1 x 12 x 8	Microtiter ELISA Plate Break apart strips coated with anti-Ustekinumab monoclonal antibody.
5 x 0.5 mL	Ustekinumab Standards A-E, Concentrate (10X) 600; 200; 60; 20; and 0 ng/mL Used for construction of the standard curve. Contains Ustekinumab, proteins, preservative and stabilizer.
1 x 12 mL	Assay Buffer Blue colored. Ready to use. Contains proteins and preservative.
1 x 60 mL	Dilution Buffer, Concentrate (5X) Orange colored. Contains proteins and preservative.
1 x 12 mL	Biotinylated $\gamma$ -hIgG Green colored. Ready to use. Contains biotinylated anti-human IgG monoclonal antibody, proteins, stabilizers and preservative.
1 x 12 mL	Enzyme Conjugate Red colored. Ready to use. Contains horseradish peroxidase(HRP)-conjugated streptavidin (HRP-Streptavidin), Proclin <sup>®</sup> and stabilizers.
1 x 12 mL	TMB Substrate Solution Ready to use. Contains 3,3',5,5'-Tetramethylbenzidine (TMB).
1 x 12 mL	Stop Solution Ready to use. 1 N Hydrochloric acid (HCl).
1 x 50 mL	Wash Buffer, Concentrate (20x) Contains buffer, Tween <sup>®</sup> 20 and Kathon <sup>™</sup> .
2 x 1	Adhesive Seal For sealing the microtiter plate during incubation.

### Required Materials that are not supplied

1. Micropipettes (< 3% CV) and tips to deliver 5-1000  $\mu$ L.
2. Bidistilled or deionised water and calibrated glasswares (e.g. flasks or cylinders).
3. Wash bottle, automated or semi-automated microtiter plate washing system.
4. Microtiter plate reader capable of reading absorbance at 450 nm (reference wavelength at 600-650 nm is optional).
5. Absorbent paper towels, standard laboratory glass or plastic vials, and a timer.



### Handling/Storage

The usual precautions for venipuncture should be observed. It is important to preserve the chemical integrity of a blood specimen from the moment it is collected until it is assayed. Do not use grossly hemolytic, icteric or grossly lipemic specimens. Samples appearing turbid should be centrifuged before testing to remove any particulate material.

Storage: 2-8°C, Stability: 3d

Keep away from heat or direct sun light Avoid repeated freeze-thaw cycles

### Procedural Notes

1. Any improper handling of samples or modification of the test procedure may influence the results. The indicated pipetting volumes, incubation times, temperatures and pre-treatment steps have to be performed strictly according to the instructions. Use calibrated pipettes and devices only.
2. Once the test has been started, all steps should be completed without interruption. Make sure that required reagents, materials and devices are prepared readily at the appropriate time. Allow all reagents and specimens to reach room temperature (20-25 °C) and gently swirl each vial of liquid reagent and sample before use. Mix reagents without foaming.
3. Avoid contamination of reagents, pipettes and wells/tubes. Use new disposable plastic pipette tips for each reagent, standard or specimen. Do not interchange the caps of vials. Always cap not used vials. Do not reuse wells or reagents.
4. Use a pipetting scheme to verify an appropriate plate layout.
5. Incubation time affects results. All wells should be handled in the same order and time sequences. It is recommended to use an 8-channel Micropipettor for pipetting of solutions in all wells.
6. Microplate washing is important. Improperly washed wells will give erroneous results. It is recommended to use a multichannel pipette or an automatic microplate washing system. Do not allow the wells to dry between incubations. Do not scratch coated wells during rinsing and aspiration. Rinse and fill all reagents with care. While rinsing, check that all wells are filled precisely with Wash Buffer, and that there are no residues in the wells.
7. Humidity affects the coated wells. Do not open the pouch until it reaches room temperature. Unused wells should be returned immediately to the resealed pouch including the desiccant.

### Preparation of Components

Dilute/dissolve	Component		Diluent	Relation	Remarks	Storage	Stability
10 mL	Wash Buffer	up to 200 mL	Distilled Water	1:20	Warm up at 37°C to dissolve crystals. Mix vigorously.	2-8 °C	4 w
10 mL	Dilution Buffer	up to 50 mL	Distilled Water	1:5		2-8 °C	4 w



### **Dilutions of Standards and Samples**

The dilutions depicted below are examples of how to obtain final dilutions for standards and samples. Standards and samples should be properly diluted as homogenous mixture before starting the assay procedure. It is recommended mixing the standards and samples well to homogenous solution at each dilution step. We are recommending that each laboratory determines the best initial dilution for their samples in order to minimize retesting.

1. 10  $\mu\text{L}$  of standard added to 90  $\mu\text{L}$  of 1X dilution buffer, giving the total volume of 100  $\mu\text{L}$  and a dilution of 1:10.
2. 10  $\mu\text{L}$  of sample added to 1990  $\mu\text{L}$  of 1X dilution buffer, giving the total volume of 2000  $\mu\text{L}$  and a dilution of 1:200.
3. Samples with a drug concentration above the measuring range should be rated as ">highest standard". The result should not be extrapolated. The sample in question should be further diluted with 1X Dilution Buffer and then retested.

### **Test Procedure**

- Before performing the assay, samples and assay kit should be brought to room temperature (about 30 minutes beforehand) and ensure the homogeneity of the solution.
- All Standards should be run with each series of unknown samples.
- Standards should be subject to the same manipulations and incubation times as the samples being tested.
- All steps of the test should be completed without interruption.
- Use new disposable plastic pipette tips for each reagent, standard or specimen in order to avoid cross contamination.
- The total pipetting time needed for dispensing all samples into the wells should not exceed 5 minutes. If this is difficult to achieve the samples should be pre-dispensed in a separate neutral polypropylene microplate and then transferred into the reaction ELISA plate by a multi channel pipette.



## Assay Procedure

<b>1.</b>	Pipette 100 $\mu$ L of Assay Buffer into each of the wells to be used.
<b>2.</b>	Pipette 75 $\mu$ L of each 1:10 Diluted Standard, and 1:200 Diluted Samples into the respective wells of the microtiter plate. Bubble formation during the pipetting of standards and samples must be avoided. Wells A1: Standard A B1: Standard B C1: Standard C D1: Standard D E1: Standard E F1 and so on: Samples (Serum/Plasma)
<b>3.</b>	Cover the plate with adhesive seal. Shake plate carefully by tapping several times. Incubate the plate on bench top for 60 min at room temperature (RT, 20-25°C).
<b>4.</b>	Remove adhesive seal. Aspirate or decant the incubation solution. Wash the plate 5 X 350 $\mu$ L of Diluted Wash Buffer per well. Remove excess solution by tapping the inverted plate on a paper towel.
<b>5.</b>	Pipette 100 $\mu$ L of Biotinylated $\alpha$ -hIgG into each well.
<b>6.</b>	Cover plate with adhesive seal. Shake plate carefully by tapping several times. Incubate the plate on a bench top for 30 min at RT.
<b>7.</b>	Remove adhesive seal. Aspirate or decant the incubation solution. Wash the plate 5 X 350 $\mu$ L of Diluted Wash Buffer per well. Remove excess solution by tapping the inverted plate on a paper towel.
<b>8.</b>	Pipette 100 $\mu$ L of Enzyme Conjugate (HRP-Streptavidin) into each well.
<b>9.</b>	Cover plate with adhesive seal. Shake plate carefully by tapping several times. Incubate the plate on a bench top for 30 min at RT.
<b>10.</b>	Remove adhesive seal. Aspirate or decant the incubation solution. Wash the plate 5 X 350 $\mu$ L of Diluted Wash Buffer per well. Remove excess solution by tapping the inverted plate on a paper towel.
<b>11.</b>	Pipette 100 $\mu$ L of Ready-to-Use TMB Substrate Solution into each well.
<b>12.</b>	Incubate 20 min at RT. Avoid exposure to direct sunlight.
<b>13.</b>	Stop the substrate reaction by adding 100 $\mu$ L of Stop Solution into each well. Color changes from blue to yellow. Briefly mix contents by gently shaking the plate.
<b>14.</b>	Measure optical density (OD) with a photometer at 450 nm (Reference at OD620 nm is optional) within 15 min after pipetting the Stop Solution.



### Quality Control

The test results are only valid if the test has been performed following the instructions. Moreover the user must strictly adhere to the rules of GLP (Good Laboratory Practice) or other applicable standards/laws. All standards/controls must be found within the acceptable ranges as stated above and/or label. If the criteria are not met, the run is not valid and should be repeated. In case of any deviation, the following technical issues should be reviewed: Expiration dates of (prepared) reagents, storage conditions, pipettes, devices, incubation conditions and washing methods.

### Calculations of Results

A standard curve should be constructed using the standard concentration (X-axis) versus the OD450 (or OD450/620) values (Y-axis). This can be done manually using graph paper or with a computer program. Concerning the data regression by computer, it is recommended to primarily use the "4 Parameter Logistic (4PL)" or alternatively the "point-to-point calculation". In case of manual plot there are 2 options: Semilog graph (see Fig. A) or linear graph (see Fig. B). Semilog graph paper is available at <http://www.papersnake.com/logarithmic/semilogarithmic/>. The concentration of the samples can be read from this standard curve as follows. Using the absorbance value for each sample, determine the corresponding concentration of the drug from the standard curve. This value always has to be multiplied by the individual dilution factor (usually 200). If any diluted sample is reading greater than the highest standard, it should be further diluted appropriately with 1X Dilution Buffer and retested. Also, this second dilution has to be used for calculation of the final result. We are recommending that each laboratory determines the best initial dilution for their samples in order to minimize retesting.

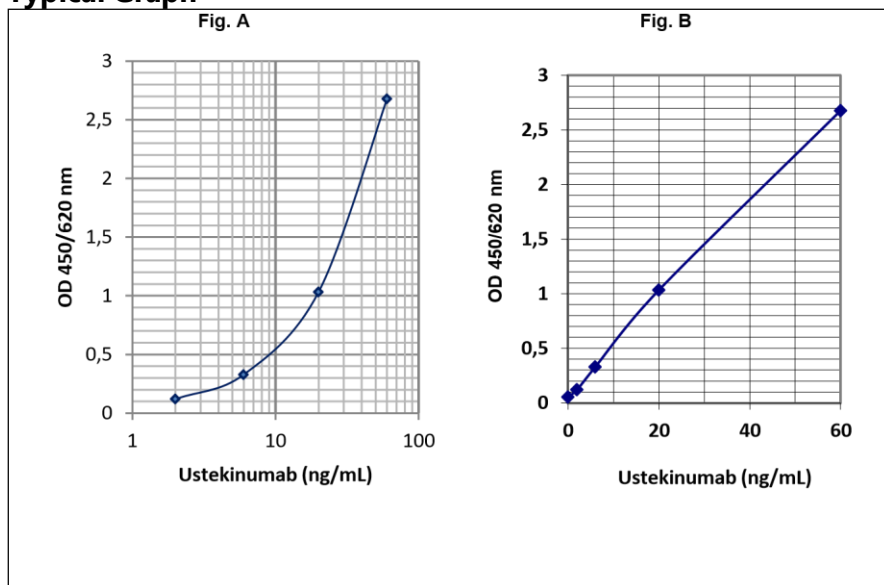
### Typical Calibration Curve

(All steps were performed at 23°C. Just an example. Do not use it for calculation!)

1:10 Diluted Standard	A	B	C	D	E
Concentration (ng/mL)	60	20	6	2	0
Mean OD450/620 nm	2.678	1.033	0.329	0.120	0.054



## Typical Graph



## Assay Characteristics

### Specificity:

There is no cross reaction with any other proteins present in native human serum. A screening test was performed with 21 different native human sera. All produced OD<sub>450/620 nm</sub> values (ranged from 0.036 to 0.068) less than the mean OD (0.120) of standard D (2 ng/mL). In addition, binding of Ustekinumab to the solid phase is inhibited by p40-containing recombinant human interleukin-12 (hIL-12) protein. Therefore, the *ImmunoGuide* Ustekinumab ELISA (mAb-Based) measures the biologically active free form of Ustekinumab, i.e. not pre-occupied by human IL-12 or IL-23 antigen. No cross reaction was observed with sera spiked with the other therapeutic antibodies including Infliximab, Rituximab, Cetuximab, Vedolizumab, Tocilizumab, Trastuzumab, Nivolumab and Bevacizumab at concentrations tested up to 40 µg/mL. All produced mean OD<sub>450/620 nm</sub> values (ranging from 0.035 to 0.058) less than standard D.

### Sensitivity:

The lowest detectable level that can be clearly distinguished from the zero standard is 1.5 ng/mL (zero standard +2SD read from the curve) under the abovedescribed conditions. Analytical sensitivity is 1.5 ng/mL, and corresponding to the detection limit (limit of quantification) of 0.3 µg/mL for undiluted clinical samples because the serum or plasma samples are instructed to be diluted at 1:200 before starting the assay.

### Precision:

Intra-assay CV: <10%.  
Inter-assay CV: <10%.

### Recovery:

Recovery rate was found to be >95% with native serum and plasma samples when spiked with exogenous Ustekinumab.





## Automation

The *ImmunoGuide* Ustekinumab ELISA (mAb-based) is suitable also for being used by an automated ELISA processor.

## References

1. Feagan BG, Sandborn WJ, Gasink C, Jacobstein D, Lang Y, Friedman JR, Blank MA, Johans J, Gao LL, Miao Y, Adedokun OJ, Sands BE, Hanauer SB, Vermeire S, Targan S, Ghosh S, de Villiers WJ, Colombel JF, Tulassay Z, Seidler U, Salzberg BA, Desreumaux P, Lee SD, Loftus EV Jr, Dieleman LA, Katz S, Rutgeerts P; UNITI-IM-UNITI Study Group, Ustekinumab as Induction and Maintenance Therapy for Crohn's Disease. *N Engl J Med.* 2016;375(20):1946-1960.
2. Zhu Y, Wang Q, Frederick B, Bouman-Thio E, Marini JC, Keen M, Petty KJ, Davis HM, Zhou H., Comparison of the pharmacokinetics of subcutaneous ustekinumab between Chinese and non-Chinese healthy male subjects across two Phase 1 studies. *Clin Drug Investig.* 2013;33(4):291-301.
3. Kavanaugh A, Puig L, Gottlieb AB, Ritchlin C, Li S, Wang Y, Mendelsohn AM, Song M, Zhu Y, Rahman P, McInnes IB; PSUMMIT 1 Study Group., Maintenance of Clinical Efficacy and Radiographic Benefit Through Two Years of Ustekinumab Therapy in Patients With Active Psoriatic Arthritis: Results From a Randomized, Placebo-Controlled Phase III Trial. *Arthritis Care Res (Hoboken).* 2015;67(12):1739-49.
4. Lamb YN, Duggan ST., Ustekinumab: A Review in Moderate to Severe Crohn's Disease. *Drugs.* 2017; 77(10):1105-1114.
5. Smolen JS, Agarwal SK, Ilivanova E, Xu XL, Miao Y, Zhuang Y, Nnane I, Radziszewski W, Greenspan A, Beutler A, Baker D., A randomised phase II study evaluating the efficacy and safety of subcutaneously administered ustekinumab and guselkumab in patients with active rheumatoid arthritis despite treatment with methotrexate. *Ann Rheum Dis.* 2017; 76(5): 831-839.
6. Deepak P, Loftus EV Jr., Ustekinumab in treatment of Crohn's disease: design, development, and potential place in therapy. *Drug Des Devel Ther.* 2016;10:3685-3698. eCollection 2016.
7. Lebwohl M, Yeilding N, Szapary P, Wang Y, Li S, Zhu Y, Reich K, Langley RG, Papp KA., Impact of weight on the efficacy and safety of ustekinumab in patients with moderate to severe psoriasis: rationale for dosing recommendations. *J Am Acad Dermatol.* 2010; 63(4): 571-9.
8. Menting SP, van den Reek JM, Baerveldt EM, de Jong EM, Prens EP, Lecluse LL, Wolbink GJ, Van der Kleij D, Spuls PI, Rispens T. The correlation of clinical efficacy, serum trough levels and antidrug antibodies in ustekinumab-treated patients with psoriasis in a clinical-practice setting. *Br J Dermatol.* 2015; 173(3): 855-7
9. van Bezooijen JS, van Doorn MBA, Schreurs MWJ, Koch BCP, Te Velthuis H, Prens EP, van Gelder T., Prolongation of Biologic Dosing Intervals in Patients With Stable Psoriasis: A Feasibility Study. *Ther Drug Monit.* 2017; 39(4): 379-386.
10. Chiu HY, Chu TW, Cheng YP, Tsai TF., The Association between Clinical Response to Ustekinumab and Immunogenicity to Ustekinumab and Prior Adalimumab. *PLoS One.* 2015; 10(11):e0142930. doi: 10.1371/journal.pone.0142930. eCollection 2015.
11. Hu C, Wasfi Y, Zhuang Y, Zhou H., Information contributed by meta-analysis in exposure-response modeling: application to phase 2 dose selection of guselkumab in patients with moderate-to-severe psoriasis. *J Pharmacokinet Pharmacodyn.* 2014; 41(3): 239-50.
12. Gottlieb A, Narang K., Ustekinumab in the treatment of psoriatic arthritis: latest findings and clinical potential. *Ther Adv Musculoskelet Dis.* 2013; 5(5): 277-85.



13. Martin PL, Bugelski PJ., Concordance of preclinical and clinical pharmacology and toxicology of monoclonal antibodies and fusion proteins: soluble targets. *Br J Pharmacol*. 2012; 166(3): 806-22.
14. Zhu Y, Hu C, Lu M, Liao S, Marini JC, Yohrling J, Yeilding N, Davis HM, Zhou H., Population pharmacokinetic modeling of ustekinumab, a human monoclonal antibody targeting IL-12/23p40, in patients with moderate to severe plaque psoriasis. *J Clin Pharmacol*. 2009; 49(2): 162-75.
15. Segal BM, Constantinescu CS, Raychaudhuri A, Kim L, Fidelus-Gort R, Kasper LH; Ustekinumab MS Investigators. Repeated subcutaneous injections of IL12/23 p40 neutralising antibody, ustekinumab, in patients with relapsing-remitting multiple sclerosis: a phase II, double-blind, placebo-controlled, randomised, dose-ranging study. *Lancet Neurol*. 2008; 7(9): 796-804.
16. Scanlon JV, Exter BP, Steinberg M, Jarvis CI. Ustekinumab: treatment of adult moderate-to-severe chronic plaque psoriasis. *Ann Pharmacother*. 2009; 43(9):1456-65.
17. Merola JF, Lockshin B, Mody EA. Switching biologics in the treatment of psoriatic arthritis. *Semin Arthritis Rheum*. 2017 Feb 8. pii: S0049-0172(16)30157-3. doi: 10.1016/j.semarthrit.2017.02.001. [Epub ahead of print]
18. Yeilding N, Szapary P, Brodmerkel C, Benson J, Plotnick M, Zhou H, Goyal K, Schenkel B, Giles-Komar J, Mascelli MA, Guzzo C. Development of the IL-12/23 antagonist ustekinumab in psoriasis: past, present, and future perspectives. *Ann N Y Acad Sci*. 2011; 1222: 30-9.
19. Farhangian ME, Feldman SR. Immunogenicity of biologic treatments for psoriasis: therapeutic consequences and the potential value of concomitant methotrexate. *Am J Clin Dermatol*. 2015; 16(4): 285-94.  
Constantin MM, Poenaru E, Constantin T, Poenaru C, Purcarea VL, Mateescu BR. Biological therapies in moderate and severe psoriasis: perspectives and certainties. *J Med Life*. 2014; 7 Spec No. 2:15-7.



## 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.