



EAGLE  
BIOSCIENCES

# Nivolumab mAb-based ELISA Assay Kit

Catalog Number:

**IG-AB113 (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 Nivolumab (Opdivo®) is a fully human IgG4 monoclonal antibody that binds specifically to programmed death-1 (PD-1), a negative regulatory receptor expressed by activated T and B lymphocytes. Binding of Nivolumab to the PD-1 receptor blocks its interaction with the ligands, PDL1 and PD-L2, thereby attenuating PD-1-mediated inhibition of the immune response, including the antitumor immune response.

The specificity of this test system is achieved by using a monoclonal antibody named "NY-11C3" for the coating of the microtiter plate. This antibody is specific for Nivolumab only and does not cross react with other PD-1 catchers.

## **Intended Use**

Enzyme immunoassay for the specific determination of free Nivolumab in serum and plasma. For research use only - not for use in diagnostic procedures.

## **Assay Principle**

This ELISA is based on Nivolumab-specific mouse monoclonal antibody (catcher Ab, *ImmunoGuide* clone NY-11C3). Diluted standards and samples are incubated in the microtiter plate coated with IG-NY-11C3 mAb. After incubation, the wells are washed. A horseradish peroxidase (HRP)-conjugated anti-human IgG antibody is added and binds to the Fc part of Nivolumab. 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 Nivolumab in the sample or standard. Results of samples can be determined by using the standard curve.



### **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-Nivolumab monoclonal antibody.
5 x 0.5 mL	Nivolumab Standards A-E, Concentrate (10X) 6000; 2000; 600; 200; and 0 ng/mL Used for construction of the standard curve. Contains Nivolumab, 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) Contains orange dye, proteins and preservative.
1 x 12 mL	Enzyme Conjugate Red colored. Ready to use. Contains horseradish peroxidase(HRP)-conjugated anti-human IgG antibody, 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 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 100µl of Assay Buffer into each of the wells to be used.
<b>2.</b>	Pipette 50 µL of each 1:10 Diluted Standard, and 1:200 Diluted Samples into the microtiter te. Bubble formation during the pipetting of ust be avoided. standards and samples m Standard A <u>Wells</u> A1: Standard B B1: Standard C C1: Standard D D1: Standard E 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 a bench top for 60 min at room temperature (RT, 2025°C).
<b>4.</b>	Remove adhesive seal. Aspirate or decant the incubation solution. Wash the plate 5 X 350 µL of Diluted Wash Buffer per well. Remove excess solution by tapping the inverted plate on a paper towel.
<b>5.</b>	Pipette 100 µL of Enzyme Conjugate (HRP-anti human IgG) 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 µL of Diluted Wash Buffer per well. Remove excess solution by tapping the inverted plate on a paper towel.
<b>8.</b>	Pipette 100 µL of Ready-to-Use TMB Substrate Solution into each well.
<b>9.</b>	Incubate 10 min at RT. Avoid exposure to direct sunlight.
<b>10.</b>	Stop the substrate reaction by adding 100 µL of Stop Solution into each well. Briefly mix contents by gently shaking the plate. Color changes from blue to yellow.
<b>11.</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 the final result. We are recommending that each laboratory determines the best initial dilution for their samples in order to minimize retesting.

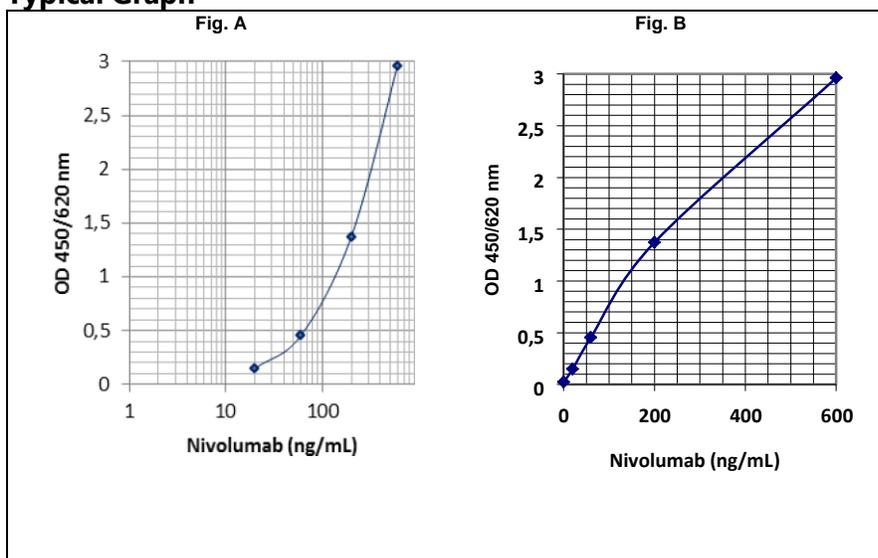
## Typical Calibration Curve

(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)	600	200	60	20	0
Mean OD450/620 nm	2.961	1.373	0.454	0.150	0.023



## 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 48 different native human sera. All produced OD<sub>450/620 nm</sub> values (ranged from 0.043 to 0.100) less than the mean OD (0.150) of standard D (20 ng/mL). In addition, binding of Nivolumab to the solid phase is inhibited by recombinant human PD-1 protein in a concentration dependent manner. Therefore, the *ImmunoGuide* Nivolumab ELISA (mAb-Based) measures the biologically active free form of Nivolumab, i.e. not pre-occupied by human PD-1 antigen. No cross reaction was observed with sera spiked with the other therapeutic antibodies including Cetuximab, Aflibercept, Rituximab, Trastuzumab, Bevacizumab, Vedolizumab, Infliximab and Adalimumab at concentrations up to 200 µg/mL. All produced mean OD<sub>450/620 nm</sub> values (ranged from 0.024 to 0.030) less than standard D.

### Sensitivity:

The lowest detectable level that can be clearly distinguished from the zero standard is 15 ng/mL (zero standard +2SD read from the curve) under the abovedescribed conditions. Analytical sensitivity is 15 ng/mL, and corresponding to the detection limit (limit of quantification) of 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 Nivolumab.



## Automation

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

## References

1. Yamamoto N, Nokihara H, Yamada Y, et al., Phase I study of Nivolumab, an anti-PD-1 antibody, in patients with malignant solid tumors, *Invest New Drugs*, DOI 10.1007/s10637-016-0411-2.
2. Bajaj G, Wang X, Agrawal S, Gupta M, Roy A and Feng Y, Model-Based Population Pharmacokinetic Analysis of Nivolumab in Patients With Solid Tumors, *CPT Pharmacometrics Syst. Pharmacol.* 2016; 00, 00; doi:10.1002/psp4.12143; published online on 0 Month 2016.
3. Agrawal S, Statkevich P, Bajaj G, Feng Y, Saeger S, et al., Evaluation of Immunogenicity of Nivolumab Monotherapy and Its Clinical Relevance in Patients With Metastatic Solid Tumors, *The Journal of Clinical Pharmacology*, 2016; 00(0) 1–7. DOI: 10.1002/jcph.818
4. Wang C, Thudium K.B, Han M, et al, In Vitro Characterization of the Anti-PD-1 Antibody Nivolumab, BMS-936558, and In Vivo Toxicology in Non-Human Primates, *Cancer Immunol Res*, 2014; 2(9); 846–56.
5. Agrawal S, Waxman I, Lambert A, Roy A, Darbenzio R, Evaluation of the potential for QTc prolongation in patients with solid tumors receiving nivolumab, *Cancer Chemother Pharmacol*, 2016; 77: 635–641.
6. Li Y, Fu T, Liu T, Guo H et al. Characterization of alanine to valine sequence variants in the Fc region of nivolumab biosimilar produced in Chinese hamster ovary cells, *mAbs*, 2016; 8:5, 951-960, DOI: 10.1080/19420862.2016.1172150
7. Brahmer JR, Drake CG, Wollner I, Powderly JD, Phase I Study of Single-Agent Anti-Programmed Death-1 (MDX-1106) in Refractory Solid Tumors: Safety, Clinical Activity, Pharmacodynamics, and Immunologic Correlates, *J Clin Oncol*, 2010; 28: 3167-75
8. Wolchok JD, Kluger H, Callahan MK, Postow MA, et al. Nivolumab plus Ipilimumab in Advanced Melanoma, *N Engl J Med* 2013;369:122-33.
9. Scott LJ, Nivolumab: A Review in Advanced Melanoma, *Drugs*, 2015; 75: 1413–1424.
10. Kanda S, Goto K, Shiraishi H, Kubo E, Tanaka A, Utsumi H, Sunami K, Kitazono S, Safety and efficacy of nivolumab and standard chemotherapy drug combination in patients with advanced non-small-cell lung cancer: a four arms phase Ib study, *Annals of Oncology*, 2016; 27: 2242–50.
11. Kimura T, Fukushima S, Miyashita A, Aoi J, Jinnin M, Kosaka T, et al. Myasthenic crisis and polymyositis induced by one dose of nivolumab, *Cancer Sci*, 2016; 107: 1055–58.
12. Nakamura Y, Kitano S, Takahashi A, Tsutsumida A, et al. Nivolumab for advanced melanoma: pretreatment prognostic factors and early outcome markers during therapy, *Oncotarget*, 2016; 7(47): 77404-77415.
13. Keating GM, Nivolumab: A Review in Advanced Squamous Non-Small Cell Lung Cancer, *Drugs*, 2015; 75: 1925–34.
14. Zhou GW, Xiong Y, Chen S, Xia F, Li Q, Hu J, Anti-PD-1/PD-L1 antibody therapy for pretreated advanced non-small-cell lung cancer: A meta-analysis of randomized clinical trials, *Medicine*, 2016; 95: 35.
15. Jung K, Zeng X, Bilusic M, Nivolumab-associated acute glomerulonephritis: a case report and literature review, *BMC Nephrology* 2016; 17: 188.



16. Munakata W, Ohashi K, Yamauchi N, Tobinai K, Fulminant type I diabetes mellitus associated with nivolumab in a patient with relapsed classical Hodgkin lymphoma, *Int J Hematol*, DOI 10.1007/s12185-016-2101-4.
17. Paoluzzi L, Cacavio A, Ghesani M, Karambelkar A, Rapkiewicz A, Weber J, Rosen G, Response to anti-PD1 therapy with nivolumab in metastatic sarcomas, *Clin Sarcoma Res*, 2016; 6: 24.
18. Gardiner D, Lalezari J, Lawitz E, DiMicco M, Ghalib R, A Randomized, Double-Blind, Placebo-Controlled Assessment of BMS-936558, a Fully Human Monoclonal Antibody to Programmed Death-1 (PD-1), in Patients with Chronic Hepatitis C Virus Infection, *PLoS ONE* 8(5): e63818. doi:10.1371/journal.pone.0063818, May 2013 | Volume 8 | Issue 5 | e63818
19. Agrawal S, Feng Y, Roy A, Kollia G, Lestini B, Nivolumab dose selection: challenges, opportunities, and lessons learned for cancer immunotherapy, *Journal for ImmunoTherapy of Cancer*, 2016; 4: 72-83.
20. Costa R, Carneiro BA, Agulnik M, Rademaker AW, et al., Toxicity profile of approved anti-PD-1 monoclonal antibodies in solid tumors: a systematic review and meta-analysis of randomized clinical trials, *Oncotarget*, Advance Publications 2016.



## **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.