

Product datasheet

Nivolumab ELISA Kit (Opdivo®) ab237651

1 Image

Overview

Product name Nivolumab ELISA Kit (Opdivo®)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 15%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 15%

Sample type Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 10 ng/ml

Range 30 ng/ml - 3000 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum		85% - 115%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Nivolumab ELISA Kit (Opdivo®) (ab237651) has been developed for specific quantification of Nivolumab concentration in human serum or plasma with high sensitivity and reproducibility. The density of color is proportional to the amount of Nivolumab captured from the samples and can be quantified when compared with standard curve.

Nivolumab (Opdivo®) is a human immunoglobulin G4 (IgG4) monoclonal antibody that binds to the programmed cell death 1 (PD-1) receptor and selectively blocks interaction with its programmed death ligands PD-L1 and PD-L2. Upregulation of PD-1 ligands occurs in some tumors and

signalling through this pathway can contribute to inhibition of active T-cell immune surveillance of tumour tissue. The inhibitory effect of PD-1 and its ligands occurs through the promotion of apoptosis in antigen specific T cells while simultaneously blocking apoptosis in suppressor T cells. Blocking PD-1 activity has been shown to lead to decreased tumour growth in mouse tumour models.

Tested applications

Suitable for: Sandwich ELISA

Platform

Microplate (12 x 8 well strips)

Properties

Storage instructions

Store at +4°C. Please refer to protocols.

Components	96 tests
Micro ELISA Plate	1 unit
Nivolumab Standard S1	1 x 0.3ml
Nivolumab Standard S2	1 x 0.3ml
Nivolumab Standard S3	1 x 0.3ml
Nivolumab Standard S4	1 x 0.3ml
Nivolumab Standard S5	1 x 0.3ml
Nivolumab Standard S6	1 x 0.3ml
Nivolumab Standard S7	1 x 0.3ml
Nivolumab Standard S8	1 x 0.3ml
Assay Buffer	2 x 50ml
HRP-conjugate Probe	1 x 12ml
TMB substrate	1 x 12ml
Stop Solution	1 x 12ml
Wash buffer (20X)	1 x 50ml
Plate sealers	2 units

Relevance

Nivolumab (Opdivo®) is a human immunoglobulin G4 (IgG4) monoclonal antibody that binds to the programmed cell death 1 (PD-1) receptor and selectively blocks interaction with its programmed death ligands PD-L1 and PD-L2. Upregulation of PD-1 ligands occurs in some tumors and signalling through this pathway can contribute to inhibition of active T-cell immune surveillance of tumour tissue. The inhibitory effect of PD-1 and its ligands occurs through the promotion of apoptosis in antigen specific T cells while simultaneously blocking apoptosis in suppressor T cells. Blocking PD-1 activity has been shown to lead to decreased tumour growth in mouse tumour models.

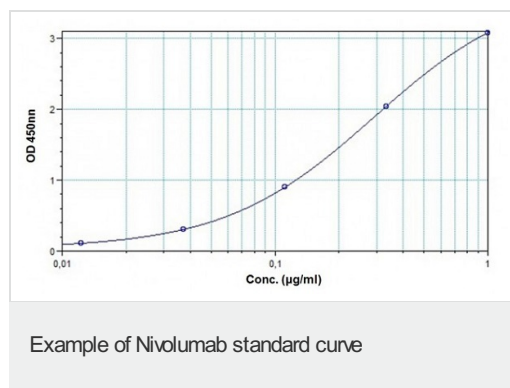
Applications

Our [Abpromise guarantee](#) covers the use of **ab237651** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Typical Standard Curve: This standard curve is for demonstration only. A standard curve must be run with each assay.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors