

Version 2 Last updated 27 June 2020

# ab237663 anti-Denosumab ELISA Kit (Prolia<sup>®</sup>)

For the measurement of the antibody against Denosumab in human serum and plasma.

This product is for research use only and is not intended for diagnostic use.

## Table of Contents

1. Overview	3
2. Protocol Summary	4
3. Precautions	5
4. Storage and Stability	5
5. Limitations	5
6. Materials Supplied	6
7. Materials Required, Not Supplied	6
8. Technical Hints	6
9. Reagent Preparation	8
10. Sample Preparation	8
11. Assay Procedure	9
12. Calculations	10
13. Cross Reactivity	10
14. Notes	11

# 1. Overview

anti-Denosumab ELISA Kit (Prolia®) (ab237663) is a highly specific and sensitive kit designed for the in vitro determination of the antibody against Denosumab in biological matrices such as human serum and plasma.

Denosumab (Prolia®) is a therapeutic fusion protein specific for Tumor Necrosis Factor-Alpha (TNF- $\alpha$ ) and is used to treat rheumatic arthritis, intestinal disorders, dermatological diseases and cancer. Denosumab specifically binds to TNF  $\alpha$  and blocks its interaction with cell surface TNF receptors and reduces the inflammation and subsequently improves the patient's health. However, some patients develop unwanted immunogenicity, which leads to production of anti-drug-antibodies (ADAs) inactivating the therapeutic effects of the treatment and, in rare cases, inducing adverse effects.

## 2. Protocol Summary

Prepare all reagents, samples, and standards as instructed. Add 100  $\mu\text{L}$  of Assay Buffer to each well.



Add 10  $\mu\text{L}$  of negative control, positive control and samples to appropriate wells. Cover and incubate for 60 minutes at room temperature



Discard incubation solution and wash plate 3 times with 300  $\mu\text{L}$  diluted Wash Buffer



Add 100  $\mu\text{L}$  peroxidase conjugate to each well. Cover and incubate for 60 minutes at room temperature



Discard the solution and wash plate 3 times with 300  $\mu\text{L}$  diluted Wash Buffer



Add 100  $\mu\text{L}$  TMB Substrate and incubate the plate in the dark at room temperature for 20 minutes.



Add 100  $\mu\text{L}$  Stop Solution and read OD at 450 nm within 20 minutes.

### 3. Precautions

Please read these instructions carefully prior to beginning the assay.

- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:  
[www.abcam.com/assaykitguidelines](http://www.abcam.com/assaykitguidelines)
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

### 4. Storage and Stability

Store kit at +4°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components.

### 5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors.

### 6. Materials Supplied

Item	Quantity	Storage
------	----------	---------

		<b>condition</b>
Micro ELISA Plate	1 unit	+4°C
Positive Control	0.3 mL	+4°C
Negative Control	1 mL	+4°C
Assay Buffer	12 mL	+4°C
Peroxidase Conjugate	12 mL	+4°C
TMB Substrate	12 mL	+4°C
Stop Solution	12 mL	+4°C
Wash Buffer (20X)	50 mL	+4°C
Plate sealers	2 units	+4°C

## 7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at OD 450 nm
  - Deionized water.
  - Multi- and single-channel pipettes.
  - Tubes for sample dilution.
  - Plate shaker for all incubation steps.
  - Absorbent paper

## 8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.

- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background.
- All samples should be mixed thoroughly and gently.
- Avoid multiple freeze/thaw of samples.
- Incubate ELISA plates on a plate shaker during all incubation steps.
- When generating positive control samples, it is advisable to change pipette tips after each step.

## 9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. Before using the kit, spin tubes and bring down all components to the bottom of tubes.
- Prepare only as much reagent as is needed on the day of the experiment.

### 9.1 20X Wash Buffer:

Dilute the 20X Wash Buffer to 1X solution in ddH<sub>2</sub>O (10 mL of Wash Buffer stock to 190 mL of ddH<sub>2</sub>O). Mix the 1X solution thoroughly by vortex manually. The working stock can be stable for 2 weeks after preparation at 4°C.

## 10. Sample Preparation

### General sample information:

- We recommend that you use fresh samples for the most reproducible assay.

### 10.1 Serum/plasma:

1. Samples are stable at 4°C for 7 days and -20°C for 6 months. Avoid freeze-and-thaw cycle.

**Δ Note:** The usual precautions for venipuncture should be observed.

## 11. Assay Procedure

- Prepare reagents within 30 minutes before the experiment.
  - Equilibrate all materials and prepared reagents to room temperature 15 minutes prior to use.
  - We recommend that you assay all standards, controls and samples in duplicate.
- 11.1 Pipette 100  $\mu$ l of Assay Buffer into each of the wells to be used.
  - 11.2 Add 10  $\mu$ L negative control (2 wells), positive control, and samples into appropriate wells. Cover wells and incubate for 60 minutes at room temperature (RT).
  - 11.3 Discard incubation solution. Wash plate 3 times each with 300  $\mu$ L of diluted Wash Buffer. Remove excess solution by tapping the inverted plate on a paper towel.
  - 11.4 Add 100  $\mu$ L of Peroxidase Conjugate into each well. Cover wells with adhesive plate sealer and incubate at room temperature for 60 minutes.
  - 11.5 Discard the solution and wash the wells as step 13.3.
  - 11.6 Add 100  $\mu$ L of 1X TMB substrate solution and incubate the plate in the dark at room temperature for 20 minutes.
  - 11.7 Add 100  $\mu$ L of Stop solution to stop the reaction.
  - 11.8 Read the absorbance in a microplate reader set to 450 nm within 20 minutes. (Reference wavelength to 650 nm).

## 12. Calculations

### 12.1 Qualitative Interpretation

- For the run to be valid, the OD 450/650 nm of positive control should be  $>1.500$  and the OD 450/650 nm of each negative control should be  $<0.150$ , if not, improper technique or reagent deterioration may be suspected and the run should be repeated.
- If "Sample OD450/650 / Negative Control OD450/650" is  $< 3$ , the sample is NEGATIVE for Antibody to Denosumab.
- If "Sample OD450/650 / Negative Control OD450/650" is  $\geq 3$ , the sample is POSITIVE for Antibody to Denosumab.

## 13. Cross Reactivity

**Cross Reactivity:** Denosumab (Prolia®) infusion camouflages/masks the presence of antibody to Denosumab (ATD) in serum/plasma samples. Therefore, blood sampling time is critical for detection of ATD. It is convenient to obtain blood sample just before the infusion or at least 2 weeks after the infusion of Denosumab.

## 14. Notes

## Technical Support

Copyright © 2018 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

### **Austria**

wissenschaftlicherdienst@abcam.com | 019-288-259

### **France**

supportscientifique@abcam.com | 01.46.94.62.96

### **Germany**

wissenschaftlicherdienst@abcam.com | 030-896-779-154

### **Spain**

soportecientifico@abcam.com | 91-114-65-60

### **Switzerland**

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

### **UK, EU and ROW**

technical@abcam.com | +44(0)1223-696000

### **Canada**

ca.technical@abcam.com | 877-749-8807

### **US and Latin America**

us.technical@abcam.com | 888-772-2226

### **Asia Pacific**

hk.technical@abcam.com | (852) 2603-6823

### **China**

cn.technical@abcam.com | +86-21-5110-5938 | 400-628-6880

### **Japan**

technical@abcam.co.jp | +81-(0)3-6231-0940

### **Singapore**

sg.technical@abcam.com | 800 188-5244

### **Australia**

au.technical@abcam.com | +61-(0)3-8652-1450

### **New Zealand**

nz.technical@abc.com | +64-(0)9-909-7829