

Product datasheet

anti-Denosumab ELISA Kit (Prolia[®]) ab237663

Overview

Product name	anti-Denosumab ELISA Kit (Prolia [®])
Detection method	Colorimetric
Sample type	Serum, Plasma
Assay type	Sandwich (qualitative)
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	anti-Denosumab ELISA Kit (Prolia [®]) (ab237663) is a highly specific and sensitive kit designed for the <i>in vitro</i> determination of the antibody against Denosumab in biological matrices such as human serum and plasma.

Denosumab (Prolia[®]) is a human monoclonal antibody that binds to RANKL, a transmembrane or soluble protein essential for the formation, function, and survival of osteoclasts, the cells responsible for bone resorption. Prolia prevents RANKL from activating its receptor, RANK, on the surface of osteoclasts and their precursors. Prevention of the RANKL/RANK interaction inhibits osteoclast formation, function, and survival, thereby decreasing bone resorption and increasing bone mass and strength in both cortical and trabecular bone. 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.

Platform Microplate

Properties

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

Components	96 tests
Micro ELISA Plate	1 unit
Positive Control	1 x 0.25ml
Negative Control	1 x 0.5ml
Assay Buffer	1 x 12ml

Components	96 tests
Peroxidase Conjugate	1 x 12ml
TMB substrate	1 x 12ml
Stop Solution	1 x 12ml
Wash buffer (20X)	1 x 50ml
Plate sealers	2 units

Relevance

Denosumab (Prolia®) is a human monoclonal antibody that binds to RANKL, a transmembrane or soluble protein essential for the formation, function, and survival of osteoclasts, the cells responsible for bone resorption. Prolia prevents RANKL from activating its receptor, RANK, on the surface of osteoclasts and their precursors. Prevention of the RANKL/RANK interaction inhibits osteoclast formation, function, and survival, thereby decreasing bone resorption and increasing bone mass and strength in both cortical and trabecular bone.

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