

Anti-CoV2RBD-imd-mlgG2a

Monoclonal mouse IgG2a antibody against SARS-CoV-2 Spike (REGN10987; Imdevimab)

Catalog code: srbdc4-mab10

<https://www.invivogen.com/sars2-spike-regn-mab>

For research use only, not for diagnostic or therapeutic use

Version 21F11-NJ

PRODUCT INFORMATION

Contents:

• 100 µg of Anti-CoV2RBD-imd-mlgG2a, provided azide-free and lyophilized

Target: SARS-CoV-2 Spike receptor binding domain (S-RBD)

Source: CHO cells

Isotype: Mouse IgG2a

Light chain type: Kappa

Clonality: Monoclonal

Purification: By affinity chromatography with protein A

Formulation: 0.2 µm filtered solution in a sodium phosphate buffer with glycine, saccharose, and stabilizing agents

Storage

- Product is shipped at room temperature. Store lyophilized antibody at -20 °C. Lyophilized product is stable for at least 1 year.

- Reconstituted antibody is stable for 1 month when stored at 4°C and for 1 year when aliquoted and stored at -20 °C. Avoid repeated freeze-thaw cycles.

Quality control

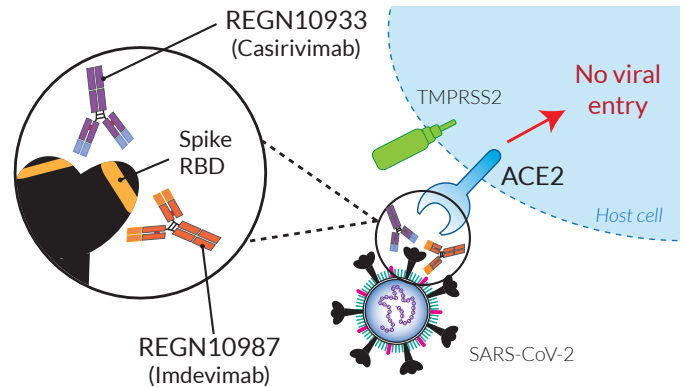
- The complete sequence of the antibody construct has been verified.
- Anti-CoV2RBD-imd-mlgG2a has been functionally validated by ELISA using Spike-RBD proteins derived from SARS-CoV-2 variants.
- Absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and TLR4 cellular assays.

PRODUCT DESCRIPTION

Anti-CoV2RBD-imd-mlgG2a is derived from 'REGN10987', a recombinant SARS-CoV-2 neutralizing monoclonal antibody (mAb)¹. Anti-CoV2RBD-imd-mlgG2a features a variable region that is reactive against the receptor binding domain (RBD) of the Spike protein of SARS-CoV-2, and the constant region of the mouse IgG2a (mlgG2a) isotype. Anti-CoV2RBD-cas-mlgG2a was generated by recombinant DNA technology, produced in CHO cells, and purified by affinity chromatography with protein A.

SARS-CoV-2 Spike RBD mAb (REGN10987; Imdevimab)

The SARS-CoV-2 Spike receptor-binding domain (S-RBD) is an important candidate for both treatment and vaccination strategies in the context of COVID-19. REGN10987 (aka Imdevimab) and REGN10933 (aka Casirivimab) mAbs were generated using both humanized mice and convalescent patients¹. These mAbs bind two distinct and non-overlapping regions of the SARS-CoV2 S-RBD, and potently block its binding to the ACE2 receptor on target cells¹. Treatment using a combination of these mAbs has been authorized under emergency use for COVID-19 outpatients^{2,4}. Prophylactic and therapeutic administration of this mAb cocktail almost completely blocks the establishment of virus infection in non-human primates⁵



and enhances clearance of the virus in patients whose endogenous immune response has not yet been initiated².

APPLICATIONS

Anti-CoV2RBD-imd-mlgG2a can be used either alone or in antibody cocktails to assess its/their neutralization efficiency against SARS-CoV-2 variants. This tool can allow the monitoring of SARS-CoV-2 variant escape.

1. Hansen J. et al. 2020. Studies in humanized mice and convalescent humans yield a SARS-CoV-2 antibody cocktail. Science 369:1010-1014. 2. Weinreich D.M. et al., 2021. REGN-COV2, a Neutralizing Antibody Cocktail, in Outpatients with Covid-19. N Engl J Med. 384(3):238-251. 3. <https://www.fda.gov/media/145610/download>. 4. <https://www.ema.europa.eu/en/news/ema-issues-advice-use-regn-cov2-antibody-combination-casirivimab-imdevimab>. 5. Baum A. et al., 2020. REGN-COV2 antibodies prevent and treat SARS-CoV-2 infection in rhesus macaques and hamsters. Science 370:1110-1115.

METHODS

Anti-CoV2RBD-imd-mlgG2a resuspension (100 µg/ml)

Note: Ensure you see the lyophilized pellet before resuspension.

- Add 1 ml of sterile water to the vial and gently pipette until completely resuspended.

- Prepare aliquots and store at 4°C or -20°C until required.

RELATED PRODUCTS

Product	Catalog Code
Anti-CoV2RBD-cas-mlgG2a	srbdc3-mab10
Anti-CoV2RBD-bam-mlgG2a	srbdc5-mab10
Anti-CoV2RBD-ete-mlgG2a	srbdc6-mab10
Spike-S1-Fc (D614G)	fc-sars2-s1g
A549-hACE2-TMPRSS2	a549-hace2tpsa

Note: For more products related to COVID-19 research, please visit our website <https://www.invivogen.com/covid-19>

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com