

Anti-mIL-28b-IgG

Neutralizing monoclonal mouse antibody against mouse interleukin 28B

Catalog code: mabg-mil28b, mabg-mil28b-5

<https://www.invivogen.com/anti-mil28b-igg>

For research use only, not for diagnostic or therapeutic use

Version 22H11-MM

PRODUCT INFORMATION

Contents: Anti-mIL-28b-IgG purified antibody is provided azide-free and lyophilized. It is available in two pack sizes:

- 100 µg: mabg-mil28b
- 5 x 100 µg: mabg-mil28b-5

Target: Natural and recombinant mouse interleukin-28B (IL-28B)

Specificity: Reacts with mouse IL-28A and mouse IL-28B. No cross-reactivity with human IL-28A or human IL-28B.

Clone: 3C11

Isotype: Mouse IgG1

Light chain type: Kappa

Immunogen: Mouse IL-28B protein expressed in Swiss mice following DNA immunization

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

Applications: Block/neutralize

Antibody resuspension (0.1 mg/ml)

Add 1 ml of sterile water per 100 µg vial.

Storage and stability

- Product is shipped at room temperature. Upon receipt, store lyophilized antibody at -20 °C.
- Reconstituted antibody is stable for 1 month at 4 °C and for 1 year at -20 °C. Avoid repeated freeze-thaw cycles.

Quality control

- This product has been validated for neutralization using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

BACKGROUND

Interleukin-28 (IL-28) is a member of the type III interferon (IFN lambda) cytokine family that exhibits several common features with type I IFNs: antiviral activity and antitumor activity^{1,2}. IL-28 exerts its action following binding to a heterodimeric protein complex composed of two subunits, IFN lambda receptor 1 (IFNLR1) and IL-10 receptor beta (IL10RB), leading to signaling through the Jak/Stat pathway and inducing the expression of IFN-stimulated genes. IL-28 comes in two isoforms, IL-28A (also known as interferon lambda 2) and IL-28B (also known as interferon lambda 3)¹. The IL-28A isoform has been implicated in the promotion of Th1 cell differentiation and the inhibition of Th2-mediated allergic inflammation³.

1. Donnelly RP. & Kotenko SV., 2010. Interferon-lambda: a new addition to an old family. *J Interferon Cytokine Res.* 30(8):555-64. 2. Li M. et al., 2009. Interferon-lambda: the modulators of antiviral, antitumor, and immune responses. *J. Leukoc. Biol.* 86:23-32. 3. Koltsida O. et al., 2011. IL-28A (IFN-λ2) modulates lung DC function to promote Th1 immune skewing and suppress allergic airway disease. *EMBO Mol Med.* 3(6):348-61.

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

DESCRIPTION

Anti-mIL-28b-IgG is a fully mouse monoclonal antibody against murine IL-28B (mIL-28B). This autoantibody was raised in mice by a proprietary method designed to induce the production of anti-cytokine antibodies directly in the animal. Anti-mIL-28b-IgG has been selected for its ability to efficiently neutralize the biological activity of mIL-28B. This antibody is produced in hybridomas and purified by affinity chromatography.

APPLICATIONS

Anti-mIL-28b-IgG is a neutralizing antibody, it blocks mIL-28-induced cellular activation *in vitro*, as described below. Furthermore, as anti-mIL-28b-IgG is a mouse anti-mouse antibody, it could be used for neutralization assays *in vivo*.

Neutralization

The exact concentration of antibody required to neutralize mIL-28B activity is dependent on the cytokine concentration, cell type and growth conditions. InvivoGen has determined the neutralization dose for this antibody using recombinant murine IL-28B and HEK-Blue™ IFN-α/β cells. These cells are HEK293 cells stably expressing the human STAT2 and IRF9 genes, and an IFN-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene. These cells respond to type I IFNs (IFN-α/β) and to a lesser extent to type III IFNs (IFN-λ).

Procedure for neutralization using HEK-Blue™ IFN-α/β cells

1. Prepare a cell suspension at ~300,000 cells/ml.
2. Add 20 µl of Anti-mIL-28b-IgG or control antibody (10 ng-1 µg/ml final concentration) per well of a 96-well plate.
Note: We recommend using Mouse IgG1 Control (which targets E. coli β-galactosidase) as a negative control.
3. Add 20 µl of recombinant mIL-28B (10 ng/ml final concentration).
4. Incubate 30 minutes at 37 °C.
5. Add 160 µl of cell suspension (~50,000 cells) per well.
6. Incubate overnight at 37 °C.
7. Add 20 µl of supernatant to 180 µl QUANTI-Blue™ Solution in a 96-well plate.
8. Incubate 18-24 hours at 37 °C.
9. Assess SEAP levels with the naked eye or spectrophotometrically by reading the optical density at 655 nm. QUANTI-Blue™ Solution turns blue following cytokine stimulation but remains pink if neutralization occurs.

RELATED PRODUCTS

Product	Description	Cat. Code
HEK-Blue™ IFN-α/β Cells	IFN-α/β reporter cells	hkb-ifnab
Mouse IgG1 Control	Isotype control antibody	mabg1-ctrlm
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs

