

# Anti-mIL-28b-mIgG1 InvivoFit™

Neutralizing monoclonal mouse antibody against murine interleukin 28B

Catalog code: mil28b-mab9-1

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

For research use only

Version 22J24-AK

## PRODUCT INFORMATION

**Contents:** 1 mg purified Anti-mIL-28b-mIgG1 InvivoFit™ monoclonal antibody (mAb) is provided azide-free and lyophilized.

**Target:** Murine 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

**Source:** Chinese hamster ovary (CHO) cells

**Isotype:** Mouse IgG1, kappa

**Purification:** Affinity chromatography with protein A

**Formulation:** 0.2 µm filtered solution in a sodium phosphate buffer with 5% saccharose

**Tested applications:** Neutralization & blocking *in vitro*

### Antibody resuspension (2 mg/ml)

*Note:* Ensure you see the lyophilized pellet before resuspension.

Resuspend Anti-mIL-28b-mIgG1 InvivoFit™ with sterile water:

Add 500 µl to 1 mg.

### 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 complete sequence of this antibody has been verified.

- <5% aggregates (confirmed by size exclusion chromatography).

- Anti-mIL-28b-mIgG1 InvivoFit™ is filter-sterilized (0.2 µm) and its endotoxin level is <1 EU/mg (determined by the LAL assay).

## BACKGROUND

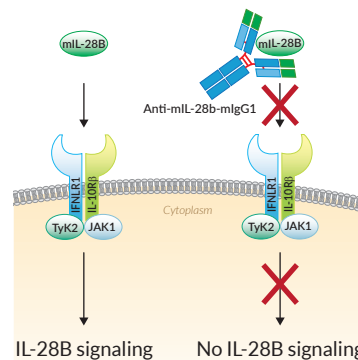
Interleukin-28 (IL-28) is a member of the type III interferon (IFN-λ) cytokine family that exhibits several common features with type I IFNs: antiviral and antitumor activity<sup>1, 2</sup>. IL-28 exerts its action following binding to a heterodimeric protein complex composed of two subunits, IFN lambda receptor 1 (IFNLR1) and IL-10 receptor beta (IL10Rβ), leading to signaling through the Jak/Stat pathway and inducing the expression of IFN-stimulated genes. IL-28 comes in two isoforms, IL-28A (IFN-λ2) and IL-28B (IFN-λ3)<sup>1</sup>.

## DESCRIPTION

Anti-mIL-28b-mIgG1 InvivoFit™ is a recombinant monoclonal antibody (mAb) specific against mIL-28B. Its sequence is 100% murine (constant and variable regions), as the original clone (3C11) was raised in mice using a proprietary method. This feature ensures high antibody performance and overcomes immunogenic events, especially upon repeated injections *in vivo*. Anti-mIL-28b-mIgG1 InvivoFit™ has been selected for its ability to efficiently neutralize the biological activity of mIL-28B. It is generated by recombinant DNA technology, produced in CHO cells and purified by affinity chromatography.

## APPLICATIONS

Anti-mIL-28b-mIgG1 InvivoFit™ is a neutralizing antibody, which blocks mIL-28B-induced cellular activation. InvivoFit™ is a high-quality standard specifically adapted for *in vivo* studies. This product is filter-sterilized (0.2 µm) and filled under strict aseptic conditions in a clean room. This mAb is also available as a standard grade for *in vitro* experiments.



Neutralization activity of Anti-mIL-28b-mIgG1 InvivoFit™

## RELATED PRODUCTS

Product	Cat. Code
Anti-mIL-28b-mIgG1	mil28b-mab9-02
Anti-β-Gal-mIgG1 InvivoFit™	bgal-mab9-1
HEK-Blue™ IFN-α/β Cells	hkb-ifnab
HEK-Blue™ IFN-λ Cells	hkb-ifnl
QUANTI-Blue™ Solution	rep-qbs

1. Donnelly RP, & Kolenko 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.

### 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](mailto:info@invivogen.com)