

Anti-hCD20-hIgG4 (S228P)

Engineered monoclonal human IgG4 antibody against human CD20

Catalog code: hcd20-mab14, hcd20-mab14-03

<https://www.invivogen.com/anti-human-cd20-igg4s228p-rituximab>

For research use only, not for diagnostic or therapeutic use

Version 23L21-MM

PRODUCT INFORMATION

Contents: Anti-hCD20-hIgG4 (S228P) purified monoclonal antibody (mAb) is provided azide-free and lyophilized. It is available in two quantities:

hcd20-mab14: 100 µg Anti-hCD20-hIgG4 (S228P)

hcd20-mab14-03: 3 x 100 µg Anti-hCD20-hIgG4 (S228P)

Target: Human CD20

Clonality: Monoclonal antibody

Isotype: Human IgG4 (S228P), kappa

Source: CHO cells

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

Purity: Purified by affinity chromatography with protein G

Storage and stability

- Product is shipped at room temperature. Upon receipt, store at -20°C.
- 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

- Binding to human CD20 has been tested using flow cytometry.
- The complete sequence of this antibody has been verified.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Anti-hCD20-hIgG4 (S228P) features the variable region of rituximab and a mutated constant region of the human IgG4 isotype reported to reduce Fab-arm exchange¹. Rituximab is a mouse/human chimeric monoclonal antibody that targets the CD20 antigen found on the surface of malignant and normal B lymphocytes. Binding of rituximab to CD20 results in cell destruction through different mechanisms including direct signaling of apoptosis, complement activation and cell-mediated cytotoxicity. Rituximab has been approved by the FDA for the treatment of various lymphoid malignancies, including B-cell non-Hodgkin's lymphoma and B-cell chronic lymphocytic leukemia.

Anti-hCD20-hIgG4 (S228P) is an engineered human IgG4 isotype. IgG4 antibodies display low antibody-dependent cytotoxicity (CDC). IgG4s are dynamic molecules that exchange Fab arms by swapping a heavy chain and attached light chain (half molecule) with a heavy-light chain pair from another molecule, resulting in bispecific antibodies^{1, 2}. IgG4 molecules thereby lose their ability to cross-link antigen and to form immune complexes under most conditions². Thus, mutations that prevent Fab-arm exchange in vivo should be considered when designing therapeutic IgG4. The constant region of Anti-hCD20-hIgG4 (S228P) contains the S228P mutation which reduces Fab-arm exchange by stabilizing the disulfides in the core-hinge of the IgG4 molecules¹. Anti-hCD20-hIgG4 (S228P) was generated by recombinant DNA technology. It has been produced in CHO cells and purified by affinity chromatography with protein G.

1. Labrijn AF. et al., 2009. Therapeutic IgG4 antibodies engage in Fab-arm exchange with endogenous human IgG4 in vivo. Nat Biotechnol. 27(8):767-71.

2. van der Neut Kolfsochten M. et al., 2007. Anti-inflammatory activity of human IgG4 antibodies by dynamic Fab arm exchange. Science. 317(5844):1554-7.

APPLICATIONS

Anti-hCD20-hIgG4 (S228P) targets the human CD20 antigen found on the surface of malignant and normal B lymphocytes.

ANTIBODY ISOTYPE COLLECTION

For your research, InvivoGen provides an anti-hCD20 isotype collection. This collection consists of monoclonal antibodies comprising the variable region of rituximab, and the constant region of human (h) and mouse (m) isotypes; hIgG1, hIgG1fut, hIgG1NQ, hIgG2, hIgG3, hIgG4, hIgG4(S228P), hIgA2, mIgG1, and mIgG2a. The isotypes differ in their functional locations and effector functions, such as complement-dependent cytotoxicity (CDC) and antibody-dependent cell-mediated cytotoxicity (ADCC), as presented in the table below.

Effector functions	Human isotypes							Mouse isotypes	
	IgG1	IgG1 fut	IgG1 NQ	IgG2	IgG3	IgG4 and IgG4 (S228P)	IgA2	IgG1	IgG2a
ADCC	++	++++	-	+/-	++	+/-	+	+/-	++
ADCP	+++	+++	-	+/-	++	+	+	+	+++
CDC	++	++	+/-	+	+++	-	-	-	++

METHODS

Antibody resuspension (100 µg/ml)

Note: Ensure you see the lyophilized pellet before resuspension.

- Add 1 ml of sterile water to 100 µg and gently pipette until completely resuspended.

RELATED PRODUCTS

Product	Catalog Code
Anti-β-Gal-hIgG4 (S228P)	bgal-mab114
Anti-hCD20-hIgG1	hcd20-mab1
Anti-hCD20-hIgG1fut	hcd20-mab13
Anti-hCD20-hIgG1NQ	hcd20-mab12
Anti-hCD20-hIgG2	hcd20-mab2
Anti-hCD20-hIgG3	hcd20-mab3
Anti-hCD20-hIgG4	hcd20-mab4

Other antibody isotype families are available, such as Anti-hTNF-α, Anti-hPD1 and Anti-βGal (control). For more information, please visit www.invivogen.com/biosimilar-antibody-isotypes.

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

