

ENGINEERING FC REGIONS FOR ALTERED PROPERTIES

The Fc region of an antibody mediates its serum half-life and effector functions, such as complement-dependent cytotoxicity (CDC), antibody-dependent cellular cytotoxicity (ADCC) and antibody-dependent cell phagocytosis (ADCP).

Engineering the Fc region of a therapeutic monoclonal antibody or Fc fusion protein allows the generation of molecules that are better suited to the pharmacology activity required of them^[1].

ENGINEERED FC REGIONS FOR INCREASED HALF-LIFE

One approach to improve the efficacy of a therapeutic antibody is to increase its serum persistence, thereby allowing higher circulating levels, less frequent administration and reduced doses. The half-life of an IgG depends on its pH-dependent binding to the neonatal receptor FcRn. FcRn, which is expressed on the surface of endothelial cells, binds the IgG in a pH-dependent manner and protects it from degradation.

Some antibodies that selectively bind the FcRn at pH 6.0, but not pH 7.4, exhibit a higher half-life in a variety of animal models.

Several mutations located at the interface between the CH2 and CH3 domains, such as T250Q/M428L^[2] and M252Y/S254T/T256E + H433K/N434F^[3], have been shown to increase the binding affinity to FcRn and the half-life of IgG1 in vivo.

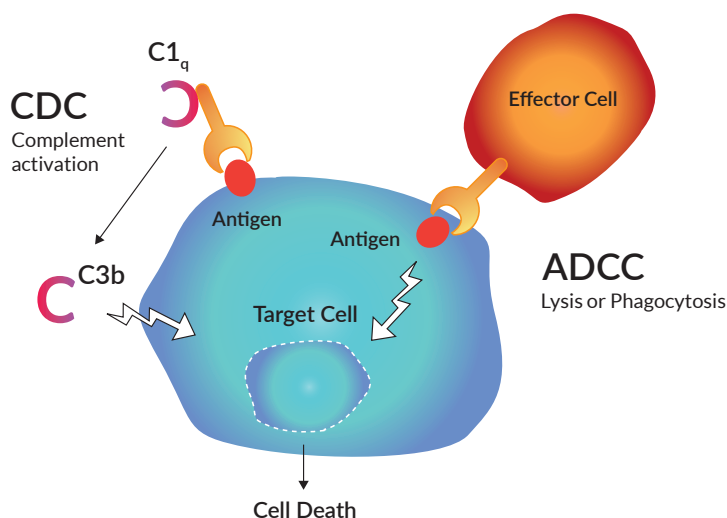
However, there is not always a direct relationship between increased FcRn binding and improved half-life^[4].

ENGINEERED FC REGIONS FOR ALTERED EFFECTOR FUNCTION

Depending on the therapeutic antibody or Fc fusion protein application, it may be desired to either reduce or increase the effector function. For antibodies that target cell-surface molecules, especially those on immune cells, abrogating effector functions is required. Conversely, for antibodies intended for oncology use, increasing effector functions may improve the therapeutic activity.

The four human IgG isotypes bind the activating Fcγ receptors (FcγRI, FcγRIIa, FcγRIIIa), the inhibitory FcγRIIb receptor, and the first component of complement (C1q) with different affinities, yielding very different effector functions^[5].

Binding of IgG to the FcγRs or C1q depends on residues located in the hinge region and the CH2 domain. Two regions of the CH2 domain are critical for FcγRs and C1q binding, and have unique



sequences in IgG2 and IgG4. Substitutions into human IgG1 of IgG2 residues at positions 233-236 and IgG4 residues at positions 327, 330 and 331 were shown to greatly reduce ADCC and CDC^[6,7]. Furthermore, Idusogie et al. demonstrated that alanine substitution at different positions, including K322, significantly reduced complement activation^[8]. Similarly, mutations in the CH2 domain of murine IgG2A were shown to reduce the binding to FcγRI, and C1q^[9].

Numerous mutations have been made in the CH2 domain of human IgG1 and their effect on ADCC and CDC tested in vitro^[7-10]. Notably, alanine substitution at position 333 was reported to increase both ADCC and CDC^[7,9]. Lazar et al. described a triple mutant (S239D/I332E/A330L) with a higher affinity for FcγRIIIa and a lower affinity for FcγRIIb resulting in enhanced ADCC^[11]. The same mutations were used to generate an antibody with increased ADCC^[12]. Richards et al. studied a slightly different triple mutant (S239D/I332E/G236A) with improved FcγRIIIa affinity and FcγRIIa/FcγRIIb ratio that mediates enhanced phagocytosis of target cells by macrophages^[13].

Due to their lack of effector functions, IgG4 antibodies represent the preferred IgG subclass for receptor blocking without cell depletion. IgG4 molecules can exchange half-molecules in a dynamic process termed Fab-arm exchange. This phenomenon can occur between therapeutic antibodies and endogenous IgG4.

The S228P mutation has been shown to prevent this recombination process allowing the design of less unpredictable therapeutic IgG4 antibodies^[14].

ENGINEERED FC	ISOTYPE	SPECIES	MUTATIONS	FCR/C1Q BINDING	EFFECTOR FUNCTION	REFS
hlgG1e1-Fc	IgG1	Human	T250Q/M428L	Increased binding to FcRn	Increased half-life	2
hlgG1e2-Fc	IgG1	Human	1M252Y/S254T/ T256E + H433K/ N434F	Increased binding to FcRn	Increased half-life	3
hlgG1e3-Fc	IgG1	Human	E233P/L234V/ L235A/?G236 + A327G/A330S/ P331S	Reduced binding to Fc γ RI	Reduced ADCC and CDC	6,7
hlgG1e4-Fc	IgG1	Human	E333A	Increased binding to Fc γ RIIIa	Increased ADCC and CDC	8,10
hlgG1e5-Fc	IgG1	Human	S239D/A330L/ I332E	Increased binding to Fc γ RIIIa	Increased ADCC	11,12
hlgG1e6-Fc	IgG1	Human	P257I/Q311	Increased binding to FcRn	Unchanged half-life	4
hlgG1e7-Fc	IgG1	Human	K326W/E333S	Increased binding to C1q	Increased CDC	9
hlgG1e9-Fc	IgG1	Human	S239D/I332E/ G236A	Increased Fc γ RIIIa/Fc γ RIIIb ratio	Increased macrophage phagocytosis	13
hlgG2e1-Fc	IgG1	Human	K322A	Reduced binding to C1q	Reduced CDC	8
hlgG4e1-Fc	IgG4	Human	S228P	-	Reduced Fab-arm exchange	14
mlgG2Ae1-Fc	IgG2a	Mouse	L235E + E318A/ K320A/K322A	Reduced binding to Fc γ RI and C1q	Reduced ADCC and CDC	9

REFERENCES

1. Strohl WR., 2009. Optimization of Fc-mediated effector functions of monoclonal antibodies. *Curr Opin Biotechnol.* 20(6):685-91. Review.
2. Hinton PR. et al., 2004. Engineered human IgG antibodies with longer serum half-lives in primates. *J Biol Chem.* 279(8):6213-6.
3. Vaccaro C. et al., 2005. Engineering the Fc region of immunoglobulin G to modulate in vivo antibody levels. *Nat Biotechnol.* 23(10):1283-8.
4. Datta-Mannan A. et al., 2007. Humanized IgG1 Variants with Differential Binding Properties to the Neonatal Fc Receptor : Relationship to Pharmacokinetics in Mice and Primates. *Drug Metab. Dispos.* 35: 86 - 94.
5. Bruhns P. et al., 2009. Specificity and affinity of human Fc γ receptors and their polymorphic variants for human IgG subclasses. *Blood.* 113(16):3716-25.
6. Armour KL. et al., 1999. Recombinant human IgG molecules lacking Fc γ receptor I binding and monocyte triggering activities. *Eur J Immunol.* 29(8):2613-24.
7. Shields RL. et al., 2001. High resolution mapping of the binding site on human IgG1 for Fc gamma RI, Fc gamma RII, Fc gamma RIII, and FcRn and design of IgG1 variants with improved binding to the Fc gamma R. *J Biol Chem.* 276(9):6591-604.
8. Idusogie EE. et al., 2000. Mapping of the C1q binding site on rituxan, a chimeric antibody with a human IgG1 Fc. *J Immunol.* 164(8):4178-84.
9. Steurer W. et al., 1995. Ex vivo coating of islet cell allografts with murine CTLA4/Fc promotes graft tolerance. *J Immunol.* 155(3):1165- 74.
10. Idusogie EE. et al., 2001. Engineered antibodies with increased activity to recruit complement. *J Immunol.* 166(4):2571-5.
11. Lazar GA. et al., 2006. Engineered antibody Fc variants with enhanced effector function. *PNAS* 103(11): 4005-4010.
12. Ryan MC. et al., 2007. Antibody targeting of B-cell maturation antigen on malignant plasma cells. *Mol. Cancer Ther.* 6: 3009 - 3018.
13. Richards JO., et al., 2008. Optimization of antibody binding to Fc γ RIIIa enhances macrophage phagocytosis of tumor cells. *Mol Cancer Ther.* 7(8):2517-27.
14. 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.

For more information, please visit : www.invivogen.com