

Periodic Table of Antibodies

		Number of Specificities																
		1				2				3								
Fc containing antibodies			2		3		4		25		26		27		28		29	
	7		13		14		15		30		31		32		33		34	
	5		19		73		75		37		38		39		40		41	
	8		83		84		85		42		43		44		45		46	
	27		104		106		126		79		90		91		92		93	
	45		151		158		163		125		135		140		149		150	
	64		171		179		180		159		160		161		162		167	
Ab fragments									168		170		173		176		177	
	6		7		8		9		44		45		46		47		48	
	22		123		124		134		49		50		60		81		82	
	53		165		166				99		107		118		119		131	
Fc fusion proteins									132		136		137		138		141	
	0		21		22		23		132		136		137		138		141	
	4		61		62		72		132		136		137		138		141	
	4		80		87		144		132		136		137		138		141	
Fc domains		10		11		12												
		33		77		78		86		103		114		115		157		
IgG-dAb (HC N-term) IgG-dAb (LC N-term) Protein IgG-fusion (HC C-term) Half antibody-scFv (HC C-term) Half antibody-scFv (LC C-term) scFv-Fc-scFv Heterodimeric IgG (common LC) Heterodimeric IgG (common HC) Heterodimeric IgG (HC+LC) Heterodimeric Fab/scFv-Fc Heterodimeric Fab/dAb-Fc Heterodimeric scFv-Fc Tandem scFv Fab-scFv (HC C-term) Fab-scFv (LC C-term) Fab-(scFv)~ (C-term) scFv-Fab-scFv (type 1) scFv-Fab-scFv (type 2) Fab-CH2-scFv IgG-scFv (HC+LC C-term) Heterodimeric IgG-scFv (type 1) Heterodimeric IgG-scFv (type 2) Heterodimeric IgG-scFv (type 3) TCB Heterodimeric Fab-scFv/scFv-Fc Heterodimeric IgG-fusion (type 1) Heterodimeric IgG-fusion (type 2) 5Heterodimeric IgG-fusion (type 3) scFv fusion (C-term) Heterodimeric Fc fusion (N-term) Heterodimeric Fc fusion (C-term) Bispecific Fc fusion (type 1) Bispecific Fc fusion (type 2) Bispecific Fc fusion (type 3) Heterodimeric bispecific Fc fusion (type 1) Heterodimeric bispecific Fc fusion (type 2) Heterodimeric bispecific Fc fusion (type 3) Heterodimeric bispecific Fc fusion (type 4) Heterodimeric Fab/protein Fc fusion (type 1) Trispecific Fc fusion (type 1) Tandem Fc fusion (type 1) Tandembody Triple Fc fusion (type 1) Triplebody Probody Heterodimeric protein Homodimeric protein mAb² Fcab Fab-fusion (type 1) Fab-fusion (type 2) Extended hinge IgG IgA with J-chain IgM with J-chain Protein-protein fusion Tandem Fc fusion (type 2) Heterodimeric Fab-Fc Peptide-IgG-scFv Heterodimeric Fab-Fc-scFv/scFv-Fc Heterodimeric Fab/protein Fc fusion (type 2) OctomAb Heterodimeric VHH-Fc Half antibody-scFv (HC + LC C-term) Heterodimeric IgG-scFv (type 4) scFv-scFv-dAb scFv-dAb-scFv dAb-scFv-scFv scFv fusion (N-term) N-term Fc fusion-scFv Trispecific Fc fusion (type 2) Tribody Single chain homotrimer Fab-IgG Bispecific Fab-IgG IgG-Fab Minibody-scFv Heterodimeric IgG-scFv (type 5) Tandem scFv-protein fusion Fab-fusion (type 3) Fab-fusion (type 4) Protein-dAb-sc homotrimer (type 1) 113. Protein-dAb-sc homotrimer (type 2) 114. Protein-sc homotrimer (type 1) 115. Protein-sc homotrimer (type 2) 116. Heterodimeric BiTE-Fc 117. Trispecific dAb 118. Bispecific Fab2 (type 1) 119. Bispecific Fab2 (type 2) 120. Fab3 121. 121. Fab4 122. 122. Fab5 123. 123. Fab6 124. 124. Fab7 125. 125. Heterodimeric IgG-scFv (type 6) 126. 126. IgY 127. 127. Tetraivalent monospecific IgG-scFv 128. 128. Hexavalent monospecific IgG-scFv 129. 129. scFv-monomeric Fc 130. 130. Fc-scFv 131. 131. Fab-VHH (HC C-term) 132. 132. Fab-VHH (LC C-term) 133. 133. Heterodimeric IgG-scFv (type 7) 134. 134. Fab8 135. 135. Heterodimeric IgG-fusion (type 4) 136. 136. TandAb 137. Tandem Fab (domain swapped) 138. DART 139. Heterodimeric protein-Fab/scFv-Fc 140. Heterodimeric IgG-fusion (type 5) 141. Monospecific tandem scFv 142. Monospecific tandem scFv-fusion 143. IgG-dAb-protein fusion (type 1) 144. Fc fusion protein (N and C-term) 145. Tetraivalent monospecific IgG-scFv (LC fusion) 146. Trimerized protein 147. Tetramerized protein 148. Pentamerized protein 149. Heterodimeric Fab/scFab-Fc 150. Duobody 151. VH only IgG 152. dAb fusion (C-term) 153. Tetrameric VHH 154. Protein-Fc-sc homotrimer 155. Heterodimeric protein-Fc-sc homotrimer (type 1) 156. Heterodimeric protein-Fc-sc homotrimer (type 2) 157. Protein-protein-protein fusion 158. Heterodimeric VHH-Fc (type 2) 159. Heterodimeric Fab/scFv-Fc (type 2) 160. dAb-Fc-dAb 161. dAb-Fc-protein 162. Heterodimeric IgG-scFab 163. Heterodimeric mAb2 164. Heterodimeric mono scFv-Fc 165. ScFv-CL 166. Diabody (monospecific) 167. Heterodimeric scFv-Fc (type 3) 168. TandAb-Fc fusion 169. TandAb-protein fusion 170. Bispecific Fab-IgG (type 2) 171. Fab-scFc 172. Bispecific 2:1 IgG-scFv (Proprietary LC) 173. Bispecific 2:1 scFv KiH-scFv 174. Trispecific 1:1:2 scFv Fab KiH scFv 175. Trispecific 2:2:2 scFv Fab Fc-scFv (LC N-term & HC C-term) 176. Heterodimeric scFv-protein KiH bispecific 177. IgG-protein fusion (LC C-term) 178. Bispecific dAb 179. IgY - His 180. IgG - Cys </																		