

Custom polyclonal antibody programs

Classical programs



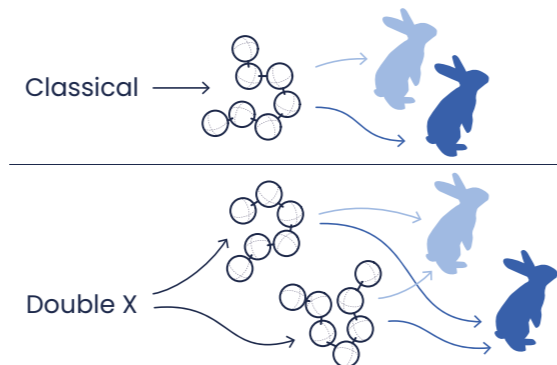
Speedy 28-day programs



		Anti-antigen		Double X (?)		Anti-PTM		Anti-antigen		Double X (?)		Anti-PTM		
		You provide the antigen* (peptide, protein, complex sample)		We design and synthesize two peptides to increase the chances of success		We design and synthesize two peptides: one modified with the PTM and one unmodified		You provide the antigen* (peptide, protein, complex sample)		We design and synthesize two peptides to increase the chances of success		We design and synthesize two peptides: one modified with the PTM and one unmodified		
Peptide	Length	x		10 - 16 aa each		10 - 13 aa each		x		10 - 16 aa each		10 - 13 aa each		
	Coupling to a carrier (KLH, BSA, OVA, THY or MAP carrier)	Optional		5 mg each		5 mg of modified peptide		Optional		5 mg of each		5 mg of modified peptide		
Immunization	Host ^b	Nbr	Cat#		Nbr	Cat#		Nbr	Cat#		Nbr	Cat#		
		Rabbit	2	AS-PNOR-3MORAB		2	AS-DOUB-DXRAB		2	AS-PSPE-PTMRAB		2	AS-SUPR-DXRAB	
		Rat	2	AS-PNOR-3MORAT		2	AS-DOUB-DXRAT		6	AS-PSPE-PTMRAT		2	AS-SUPR-DXRAT	
		Guinea pig	2	AS-PNOR-3MOGPG		2	AS-DOUB-DXGP		6	AS-PSPE-PTMGP		2	AS-SUPR-DXGP	
		Goat	1	AS-PNOR-3MOGOAT		1	AS-DOUB-DXGOAT		x	-		1	AS-SUPR-GOATAG	
		Mouse	3	On request		3	On request		x	-		x	-	
		Chicken	2	AS-PNOR-3MOHEN		2	AS-DOUB-DXHEN		x	-		x	-	
Large animal (pig, sheep, cow...)	1	On request		1	On request		x	-		x	-			
ELISA		x		Optional		Optional		x		Optional		Optional		
Antibody Processing	Affinity Purification	Optional		Optional		Double purification of 50 mL serum		Optional		Optional		Double purification of 50 mL serum		
	QC validation	Optional		Optional		ELISA and SDS-PAGE		Optional		Optional		ELISA and SDS-PAGE		
	Delivery	Serum Purified antibody (Optional)		Serum Remaining peptide Purified antibody (Optional)		Remaining serum Purified antibody		Serum Purified antibody (Optional)		Serum Remaining peptides Purified antibody (Optional)		Remaining serum Purified antibody		

Double X programs

Double X programs increase the chances of success by using two peptides corresponding to distinct regions of the targeted protein. Each host will receive a combination of the two peptides.



Legend

- x Not applicable
- Nbr Number of host
- PTM Post Translational Modification*
*Phospho; Acetyl Lys; DiAcetyl Ser; Oxidized Cys; MonoMethyl Arg; DiMethyl Arg (sym); DiMethyl Arg (asym); MonoMethyl Lys; DiMethyl Lys; TriMethyl Lys. More modifications on request.

Notes

- For more details browse our page www.eurogentec.com/en/how-to-provide-my-antigen
- Other hosts are available on request
- The guarantee only applies for peptides when approved by Eurogentec
- The guarantee does not apply for Goat immunization
- This example is subject to variations (injections and intermediate bleeds) depending on the programme selected.