

Classical programmes



Speedy 28-day programmes



	Classical programmes				Speedy 28-day programmes																	
	Anti-antigen	Anti-peptide	Double X	Anti-PTM**	Anti-antigen	Speedy-Mini	Anti-peptide	Double X	Anti-PTM**													
	You provide the antigena (peptide, protein, complex sample)	Eurogentec designs and synthesizes one peptide	Eurogentec designs and synthesizes two peptides to increase the chances of success	Eurogentec designs and synthesizes two peptides: one modified with the PTM and one unmodified	You provide the antigena (peptide, protein, complex sample)	Eurogentec designs and synthesizes one peptide	"Eurogentec designs and synthesizes one peptide"	Eurogentec designs and synthesizes two peptides to increase the chances of success	"Eurogentec designs and synthesizes two peptides: one modified with the PTM and one unmodified"													
Peptide	Length	10 - 16 aa	10 - 16 aa each	10 - 13 aa each	×	10 - 16 aa	10 - 16 aa	10 - 16 aa each	10 - 13 aa each													
	Quantity	×	15-25 mg	15-25 mg each	15-20 mg each	×	8-10 mg	15-25 mg	15-20 mg each													
	Purity		> 70 %	> 70 %	> 70 %		> 70 %	> 70 %	> 70 %													
	Coupling to a carrier (KLH, BSA, OVA, THY or MAP carrier)	+	5 mg	5 mg each	5 mg of modified peptide	+	2-3 mg to KLH	5 mg	5 mg of each	5 mg of modified peptide												
Immunization	Host ^b	Nbr	Cat#	Nbr	Cat#	Nbr	Cat#	Nbr	PTM	Cat#	Nbr	Cat#	Nbr	Cat#	Nbr	Cat#	Nbr	PTM	Cat#			
		2	AS-PNOR-3MORAB	2	AS-PCAP-RABBIT	2	AS-DOUB-LX	2	Phospho	AS-PSPE-CIFIC	2	AS-SUPR-ANTIGEN	1	AS-SMAF-SINGLE	2	AS-SUPR-SINGLE	2	AS-SUPR-DX	2	Phospho	AS-SUPR-PSPEC	
		2	-	2	-	2	-	2	Acetyl Lys	AS-PTMA-LYSACET	2	-	2	-	2	-	2	-	2	Acetyl Lys	AS-SUPR-LYSACET	
		2	-	2	-	2	-	2	DiAcetyl Ser	AS-PTMA-SERACET2	2	-	2	-	2	-	2	-	2	Citruiline	AS-SUPR-CITSPEC	
		2	-	2	-	2	-	2	Oxidized Cys	AS-PTMA-OXCYS	2	-	2	-	2	-	2	-	2	Oxidized Cys	AS-SUPR-OXCYS	
		2	-	2	-	2	-	2	MonoMethyl Arg	AS-PTMA-ARGMEI	2	-	2	-	2	-	2	-	2	MonoMethyl Arg	AS-SUPR-ARGMEI	
		2	-	2	-	2	-	2	DiMethyl Arg (sym)	AS-PTMA-ARGME2SYM	2	-	2	-	2	-	2	-	2	DiMethyl Arg (sym)	AS-SUPR-ARGME2SYM	
		2	-	2	-	2	-	2	DiMethyl Arg (asym)	AS-PTMA-ARGME2ASY	2	-	2	-	2	-	2	-	2	DiMethyl Arg (asym)	AS-SUPR-ARGME2ASY	
		2	-	2	-	2	-	2	MonoMethyl Lys	AS-PTMA-LYSMEI	2	-	2	-	2	-	2	-	2	MonoMethyl Lys	AS-SUPR-LYSMEI	
		2	-	2	-	2	-	2	DiMethyl Lys	AS-PTMA-LYSME2	2	-	2	-	2	-	2	-	2	DiMethyl Lys	AS-SUPR-LYSME2	
		2	-	2	-	2	-	2	TriMethyl Lys	AS-PTMA-LYSME3	2	-	2	-	2	-	2	-	2	-	-	
		Rat	2	AS-PNOR-3MORAT	2	AS-PCAP-RAT	2	AS-DOUB-LXRAT	6	Phospho	AS-PSPE-CIFICRAT	2	AS-SUPR-RATANTIGEN	×	-	2	AS-SUPR-RATSINGLE	2	AS-SUPR-RATDX	6	on request	-
		Guinea pig	2	AS-PNOR-3MOGPG	2	AS-PCAP-GUIPG	2	AS-DOUB-LXGUI	6	Phospho	AS-PSPE-CIFICGP	2	AS-SUPR-GPANTIGEN	×	-	2	AS-SUPR-GPSINGLE	2	AS-SUPR-GPDX	6	Phospho	AS-SUPR-PSPECGP
		Goat	1	AS-PNOR-3MOGOAT	1	AS-PCAP-GOAT	1	AS-DOUB-LXGOAT	×	-	-	1	AS-SUPR-GOATAG	×	-	1	AS-SUPR-GOATSING	1	AS-SUPR-GOATDX	×	-	-
		Mouse	3	on request	3	AS-PCAP-MOUSE	3	on request	×	-	-	×	-	×	-	×	-	×	-	×	-	-
		Chicken	2	AS-PNOR-3MOHEN	2	AS-PCAP-HEN	2	AS-DOUB-LXHEN	2	Acetyl Lys	AS-PTMA-LYSACETHEN	×	-	×	-	×	-	×	-	×	-	-
		Large animal (pig, sheep, cow,...)	1	on request	1	on request	1	on request	×	-	-	×	-	×	-	×	-	×	-	×	-	-
			-	-	-	-	-	-		More modifications on request		-	-	-	-	-	-	-	-	-	-	More modifications on request
		Antibody Processing	ELISA		×	+	+	+		+		×		+		+		+		+		+
											Double purification of 50 mL serum				Specific response at 1/20000 guarantee ^c No Result, No Serum, No Invoice ^d			Specific response at 1/20000 guarantee ^c No Result, No Serum, No Invoice ^d			Specific response at 1/20000 guarantee ^c (at least for one of the 2 peptides) ^d	
Antibody Processing	Affinity Purification	+	+	+	+	+		+		+		+		+		+		+		+		
	QC validation	+	+	+	+	+		+		+		+		+		+		+		+		
	Delivery	Serum Purified antibody (Optional)	Serum Purified antibody (Optional) Remaining peptide	Serum Purified antibody (Optional) Remaining peptide	Serum Purified antibody (Optional) Remaining peptide	Serum Purified antibody (Optional) Remaining peptide		Purified antibody Remaining serum Remaining peptides			Serum Purified antibody (Optional)	Serum Purified antibody	Serum Purified antibody (Optional) Remaining peptide	Serum Purified antibody (Optional) Remaining peptides	Serum Purified antibody (Optional) Remaining peptides		Purified antibody Remaining serum Remaining peptides					

× = not applicable + = Optional

a. See the section "How to provide my antigen" on p.25 for more details

b. Other hosts are available on request

c. The guarantee only applies for peptides when approved by Eurogentec

d. The guarantee does not apply for Goat immunization

* This example is subject to variations (injections and intermediate bleeds) depending on the programme selected.

**Post Translational Modification

Double X programmes

Double X programmes 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.

