

SPEEDY 28-DAY POLYCLONALS



TITRE guaranteed

THE NEW REFERENCE
FOR CUSTOM ANTIBODIES

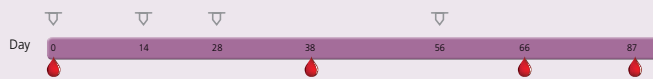
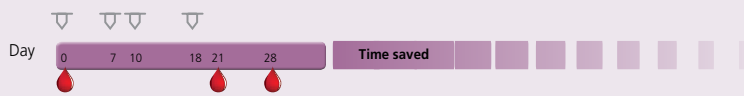


Eurogentec
Experience true partnership

GET RESULTS SOONER!

Eurogentec Speedy 28-day uses a proprietary mixture of immunostimulatory compounds (Freund's free)

▼ Speedy 28-day protocol

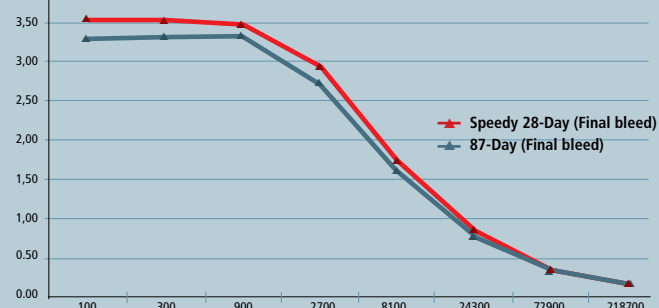


▲ Standard 87-day protocol

ANTIBODY TITRE

Our Speedy 28-day offers as good or better titre than our 87-day immunisation programme.

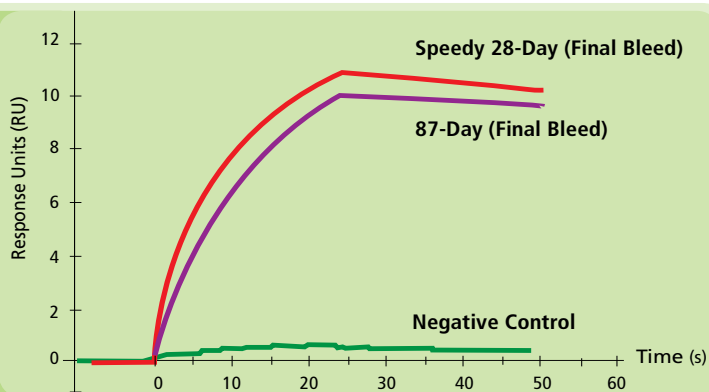
► ELISA results of the final bleed serum against a good responding peptide*



ANTIBODY AFFINITY

Our Speedy 28-day offers as good or better affinity than our 87-day protocol.

► Biacore Affinity Test*



TITRE GUARANTEE

NEW

- Free ELISA at day 21 for each **anti-peptide programme** (Epiproteomics programme excluded)
- We guarantee a significant response at 1/20 000 for at least one animal, or your money back
- For DoubleX programmes, we guarantee the response for at least one of the peptides
- The guarantee is only valid for peptides approved by Eurogentec
- We cannot guarantee the functionality of your antibody in your application

AVAILABLE HOSTS

- Rabbit, guinea pig, rat and goat are compatible with the Speedy 28-day and can be immunised with:
- Your protein
 - A single peptide
 - Two peptides (DoubleX)

* The same KLH coupled peptide was used to immunise rabbits according to our 87-day and 28-day protocols.

What is included?

- > Peptide design from your antigen sequence
 - > Peptide synthesis (> 70 %, 15-25 mg, 16 aa)
 - > Peptide coupling (KLH, BSA, OVA)
 - > 2 SPF Rabbits, Guinea pigs, Rats or one non-SPF Goat
 - > 28-day immunisation protocol (not including peptide synthesis)
 - > All immunisations, blood collections and animal care
 - > Delivery of the free remaining peptide
 - > A single shipping on dry ice
- Modification specific programmes (Epiroteomics) are also available in Rabbits.

How to start?

Anti-peptide antibodies

- > Send us your antigen sequence in single amino acid format at **peptide.design@eurogentec.com**
Please specify your preferred approach: single peptide or DoubleX
- > We will design your peptide(s) for free
- > Within 3 days, you will receive our recommended sequence(s)
- > Send us your complete anti-peptide order form**
- > You will receive an immunisation calendar
- > We will start the Speedy 28-day immunisation

Anti-protein antibodies

- > Send us your complete anti-protein order form**
- > Send us your Protein lyophilised or in solution (PAGE slices are not compatible with the Speedy 28-day)
- > You will receive an immunisation calendar
- > We will start the Speedy 28-day immunisation

**Order form can be downloaded on www.eurogentec.com/antibody-order-form.html

ORDERING INFORMATION

SPEEDY PACKAGES

Antigen	Host	REFERENCE
Your protein	Rabbit	AS-SUPR-ANTIGEN
	Guinea pig	AS-SUPR-GPANTIGEN
	Rat	AS-SUPR-RATANTIGEN
	Goat	AS-SUPR-GOATAG
1 peptide	Rabbit	AS-SUPR-SINGLE
	Guinea pig	AS-SUPR-GPSINGLE
	Rat	AS-SUPR-RATSINGLE
	Goat	AS-SUPR-GOATSING
DoubleX - 2 peptides	Rabbit	AS-SUPR-DX
	Guinea pig	AS-SUPR-GPDX
	Rat	AS-SUPR-RATDX
	Goat	AS-SUPR-GOATDX
DoubleXP - 2 peptides + affinity purification	Rabbit	AS-SUPR-DXP
	Guinea pig	AS-SUPR-GPDXP
	Rat	AS-SUPR-RATDXP
Phospho-specific***	Rabbit	AS-SUPR-PSPEC

ADDITIONAL SERVICES

Description	REFERENCE
Pre-immune sera from 5 animals for testing	AS-PREI-05
Pre-immune sera from 10 animals for testing	AS-PREI-10
Pre-immune sera from 20 animals for testing	AS-PREI-20
5 ml affinity purification against peptide	AS-PURI-SMALL
20 ml affinity purification against peptide	AS-PURI-MED
Second 20 ml affinity purification against same peptide	AS-PURI-MED02
50 ml affinity purification against peptide	AS-PURI-01
Second 50 ml affinity purification against same peptide	AS-PURI-02

*** Other Epiroteomic modifications are available on request.





EUROPEAN OFFICES

BELGIUM

EUROGENTEC HEADQUARTERS
TEL.: +32 4 372 74 00
FAX: +32 4 372 75 00
info@eurogentec.com

EUROGENTEC BIOLOGICS
TEL.: +32 4 372 74 00
FAX: +32 4 365 16 04
info.biologics@eurogentec.com

UNITED KINGDOM

EUROGENTEC LTD.
TEL.: +44 1794 511 411
FAX: +44 1794 522 417
info.uk@eurogentec.com

FRANCE

EUROGENTEC FRANCE SASU
TEL.: +33 2 41 73 33 73
FAX: +33 2 41 73 10 26
info@eurogentec.com

GERMANY & AUSTRIA

EUROGENTEC DEUTSCHLAND GMBH
TEL.: +49 221 258 94 55
FAX: +49 221 258 94 54
info@eurogentec.com

THE NETHERLANDS

EUROGENTEC NEDERLAND B.V.
TEL.: +31 43 352 06 98
FAX: +31 43 354 19 65
info@eurogentec.com

SWITZERLAND

EUROGENTEC S.A.,
SUCCURSALE DE GENÈVE
FRANÇAIS:
TEL.: +32 4 372 74 00
FAX: +32 4 372 75 00

DEUTSCH:
TEL.: +49 221 258 94 55
FAX: +49 221 258 94 54
info@eurogentec.com

CENTRAL EUROPE

REPRESENTED BY
EUROGENTEC HEADQUARTERS
TEL.: +32 4 372 74 00
FAX: +32 4 372 75 00
info@eurogentec.com

AMERICAN OFFICES

ANASPEC, INC.

TEL.: +1 510 791 9560
TOLL-FREE: +1 800 452 5530
FAX: +1 510 791 9572
info.anaspec@eurogentec.com
www.anaspec.com

EGT PRODUCTION SITE
TEL.: +1 858 793 2661
TOLL-FREE: +1 877 387 6436
FAX: +1 858 793 2666
info.usa@eurogentec.com
usa.eurogentec.com

ASIAN OFFICES

JAPAN

NIPPON EGT CO., LTD.
TEL.: +81 76 411 02 77
FAX: +81 76 452 03 99
info@n-egt.com

SINGAPORE

EUROGENTEC AIT PTE. LTD.
TEL.: +65 6778 9388
+65 6778 6822
FAX: +65 6778 5177



Eurogentec
Experience true partnership