

MOUSE MONOCLONAL ANTIBODY DEVELOPMENT



REFERENCES AND PRICES

Phase	Reference
Phase 1 - Immunisation	AS-ACMC-IM (X4) 4 mice Or AS-ACMC-IMFAST (X4) 4 mice-Fast protocol
Phase 2 – Fusion & hybridoma production	AS-ACMC-FU08
Phase 3 – Screening	AS-ACMC-AN08S
Phase 4 – Cloning & Isotyping	AS-ACMC-CIFH (1-2 clones)
Phase 4 – Additional clone	AS-ACMC-CIAH

If no positive hybridoma is obtained after phase 3, only phase 1 will be charged.

Additional quality control	Reference
Western blot	AS-ACMC-QCWB
Antibody Affinity evaluation with Biacore 2000	AS-ACMC-QCAFFI
FACS analysis	AS-ACMC-QCFACS

How to order

To start discussing your project and get a quotation adapted to your needs, please contact us at monoclonals@eurogentec.com

BENEFIT FROM OUR VAST EXPERIENCE, AND CONTINUAL DEVELOPMENTS IN MONOCLONAL PRODUCTION

MOUSE MONOCLONAL ANTIBODY DEVELOPMENT

ANTI-PROTEIN | ANTI-PEPTIDE

PROJECT DEVELOPMENT

Prior to project initiation, you can benefit from our personalised project proposals and expertise.

To qualify and build your project or to receive a quotation adapted to your needs, please contact your local representative or send us an e-mail at monoclonals@eurogentec.com

STARTING MATERIAL

You send us:

- Your **purified protein**
- Or your **peptide sequence**
- Or your **protein's accession number**

6 WEEKS

Standard Programme

- 3 immunisations (Days 0 - 21 - 36)
- 3 boosts (Before fusion)

PHASE 1 IMMUNISATION OF 4 MICE

Mouse lymphocytes
Specific Antibody Secreting Cells
Ig positive (ELISA)

PHASE 2 FUSION/HYBRIDOMA PRODUCTION

Mouse lymphocytes
Myeloma
Immortality Property
Hypoxanthine Phosphoribosyl Transferase (HPRT) negative, Ig negative

Hybridoma

Spleen cell + Myeloma cell
Sp2/OAg 14
(2 fusions per programme)

2 WEEKS

Accelerated Programme

- 1 intrapodal immunisation
- 1 boost (Day 14)
- FUSION (Day 17)

PHASE 3 SCREENING FOR POSITIVE HYBRIDOMAS

Flask contains a **mixture of hybridomas** and non fused cells

Lymphocytes die off on their own over time

Hypoxanthine Aminopterin Thymidine (HAT) treatment kills off excess myeloma cells

Only hybridomas survive

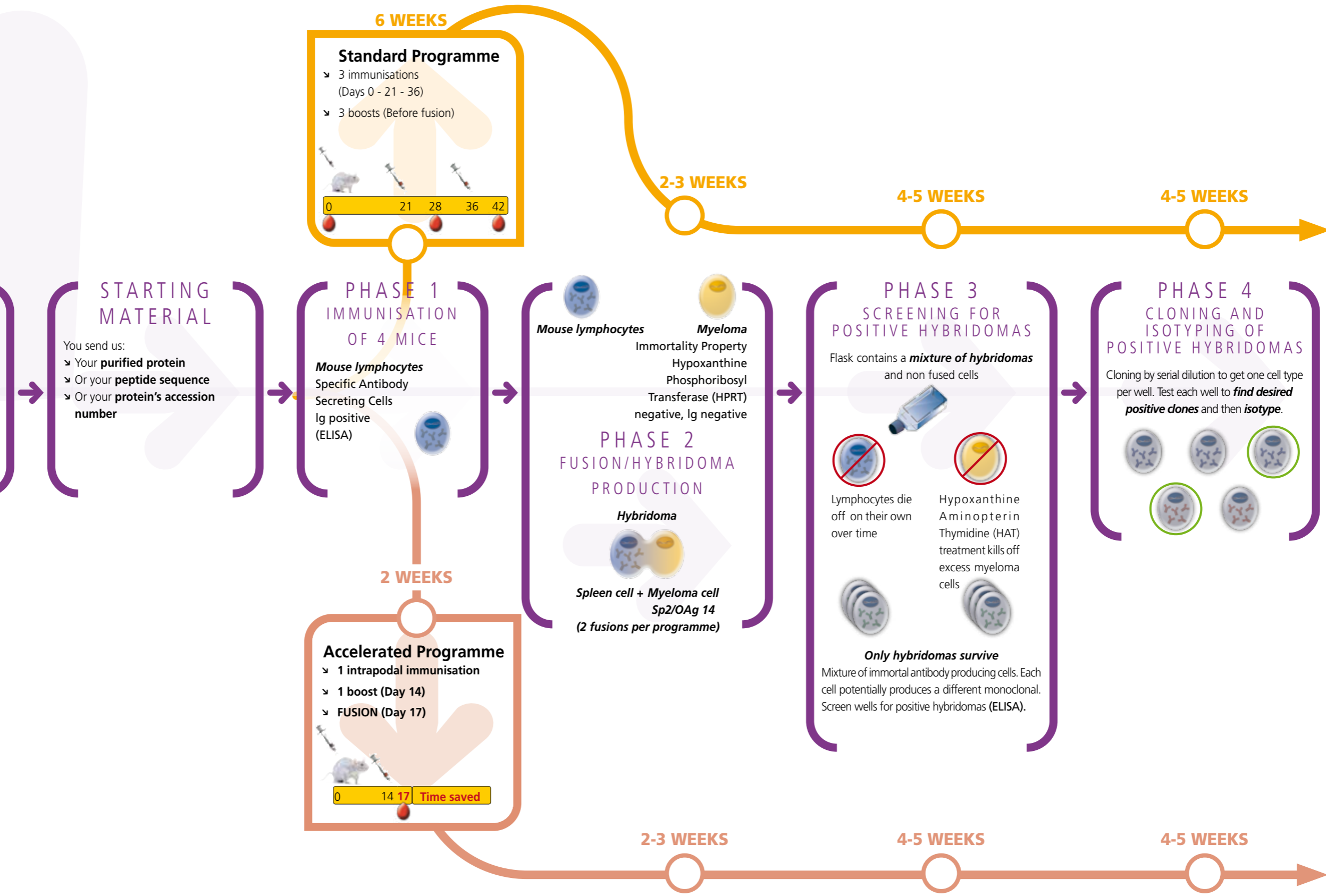
Mixture of immortal antibody producing cells. Each cell potentially produces a different monoclonal. Screen wells for positive hybridomas (ELISA).

PHASE 4 CLONING AND ISOTYPING OF POSITIVE HYBRIDOMAS

Cloning by serial dilution to get one cell type per well. Test each well to **find desired positive clones** and then **isotype**.

- Options**
- Additional clones can be selected during phase 4
 - Selection of the right Mo-Ab couple for sandwich assay development
- Additional QC**
- Specificity → Western Blot
 - Affinity → Surface Plasmon Resonance (Biacore)

- Additional Services**
- *in vitro* production of mAb
 - Antibody purification
 - Fluorescent + Enzyme coupling
 - Coupling to solid supports
 - Cell line storage
 - Luminex® Assay development
 - Flow Cytometry development
 - Peptide libraries
 - Blocking peptides



You receive

- ELISA** against the supplied antigen and serum samples
- Fusion report**
- Screening report**
 - ELISA against your antigen.
 - Delivery of > 1 ml positive supernatants (Maximum 30).
 - Customer selects best group of clones based on internal application results
- Hybridomas ELISA and isotyping report.**
 - Delivery of > 10 ml of supernatant from 1-2 positive clones.
 - Frozen hybridomas.

16-19 WEEKS

12-15 WEEKS