

# MOUSE MONOCLONAL ANTIBODY APPLICATIONS



## REFERENCES AND PRICES

Application	Service	References	Price
In vitro production	Pilot phase (including protein A purification and QC)	AS-ACMC-IVPPILOT	On request
	Production of 100 mg Ig (including protein A purification and QC)	AS-ACMC-IVPPROD100	On request
	Production of 200 mg Ig (including protein A purification and QC)	AS-ACMC-IVPPROD200	On request
	0.22 µm sterile filtration	AS-ACMC-IVPSTE200	On request
	Sterility test (bacteria and fungi)	AS-ACMC-IVPSTE210	On request
	Monoclonal antibody production roller bottles, 4 weeks	AS-ACMC-ROLLB	On request
	Monoclonal antibody production in CL1000 Bioreactor, 4 weeks	AS-ACMC-CL1000	On request
	-IgG purification - maximum 100 mg antibody from 500 ml supernatant, including QC by SDS-PAGE	AS-PURI-MCIGG500	On request
Antibody fragmentation	Test phase (1 mg) including purification	AS-ACMC-FRAGTEST	On request
	10 mg including purification and QC	AS-ACMC-FRAG10	On request
Optional quality control	Western blot	AS-ACMC-QCWB	On request
	Antibody Affinity evaluation with Biacore 2000	AS-ACMC-QCAFFI	On request
	FACS analysis	AS-ACMC-QCFACS	On request
Antibody labelling	Phycoerythrine	AS-ACMC-LABPE	On request
	Rhodamine	AS-ACMC-LABRHO	On request
	FITC	AS-ACMC-LABFITC	On request
	Alexa®	AS-ACMC-LABALEXA	On request
	BODIPY®	AS-ACMC-LABBODI	On request
	Cy3®	AS-ACMC-LABCY3	On request
	Cy®5	AS-ACMC-LABCY5	On request
	Texas Red®	AS-ACMC-LABTEXAS	On request
	Horse Radish Peroxydase (HRP)	AS-ACMC-LABHRP	On request
	Alkaline Phosphatase (AP)	AS-ACMC-LABAP	On request
	Biotin	AS-ACMC-LABBIOT	On request
Digoxigenin	AS-ACMC-LABDIG	On request	

PRO-MONOCLONARP-0409-V1

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

## How to order

To start discussing your project and get a quotation adapted to your needs, please contact us at [monoclonals@eurogentec.com](mailto:monoclonals@eurogentec.com)

# MOUSE MONOCLONAL ANTIBODY APPLICATIONS

## OUR ADDITIONAL SERVICES

### Antibody labelling

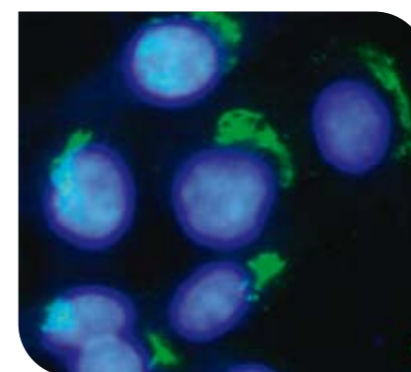
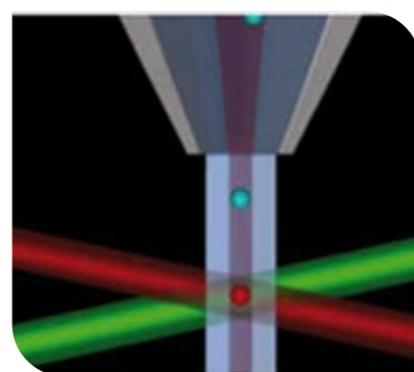
Depending on your application (DuoLink®, *in vivo/in vitro* localisation, FACS ...), your antibodies can be labelled with:

- Phycoerythrine
- Rhodamine
- FITC
- Alexa®
- BODIPY®
- Cy®3, Cy5®
- Texas Red®
- Horse Radish Peroxydase (HRP)
- Alkaline Phosphatase
- Biotin
- Digoxigenin...



### Fragmentation

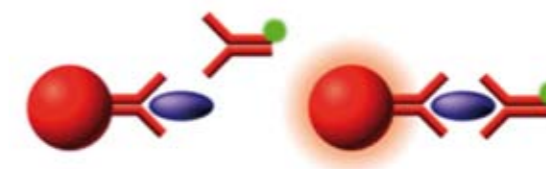
Your antibodies are selectively cleaved into Fab, F(ab')<sub>2</sub> and Fc fragments.



### Assay development

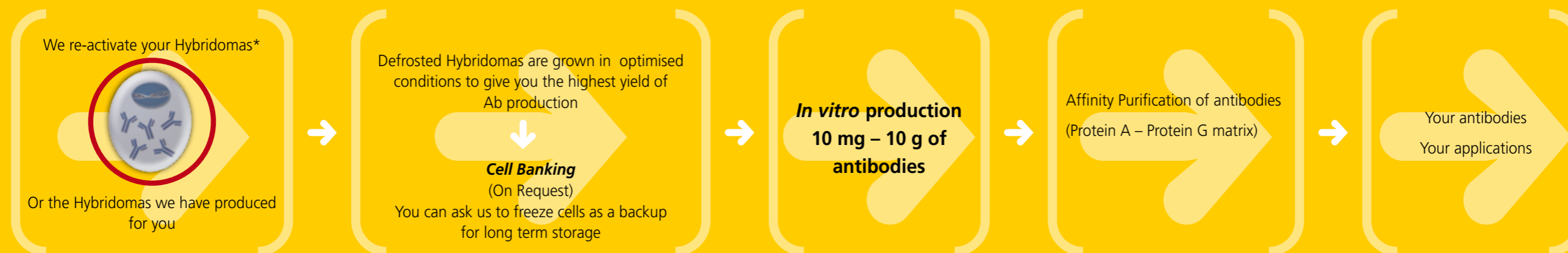
**ELISA:** Your antibodies can be coated on plates.

**Luminex:** Your antibodies can be coated on microspheres.



**Purification:** Your antibodies can be coated on beads for affinity purification, co-immunoprecipitation...

## OUR FLOWCHART



\* We can reactivate your rat Hybridomas (On request)