

# Monoclonal antibody production



## A comprehensive monoclonal antibody production service.

Starting from in-house or customer's hybridomas, we propose different production schemes. Each strategy is based on the mAb amount and the clone productivity.



### **In vitro production**

All our custom monoclonal antibodies are produced exclusively *in vitro*.



### **Up to 100 g mAb**

We are equipped to produce from mg to grams of purified monoclonal antibodies. Test studies are performed for requests  $\geq 200$  mg.



### **Optimized protocols**

Depending on the requested amount and the hybridoma productivity, we produce monoclonal antibodies in flasks, roller bottles or bioreactors.

## Downstream processes

We offer a variety of manufacturing options, to deliver monoclonal antibodies in a format that will optimally suit the customer's application.



### **Purity**

- **Polishing**  
Mixed-mode chromatography
- **Purification**  
Affinity, ion exchange, gel filtration, IgM purification
- **Clearance**  
DNA and endotoxin removal



### **Format**

- **Concentration/diafiltration**
- **Formatting**  
Conjugation and freeze-drying service

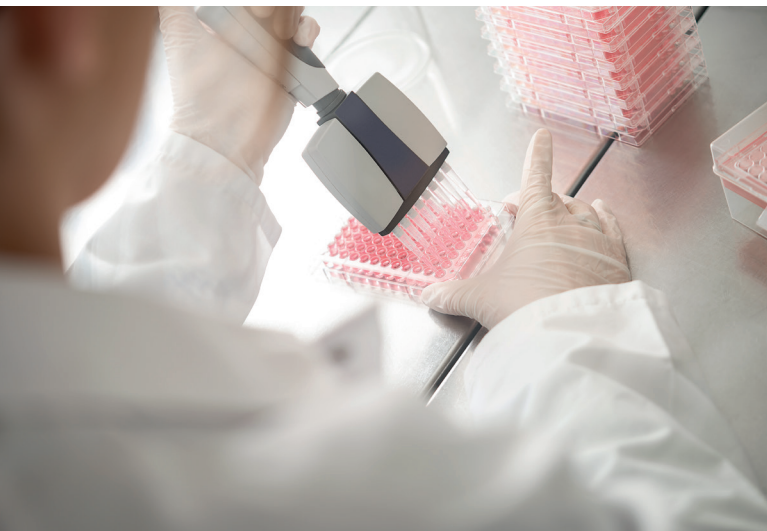
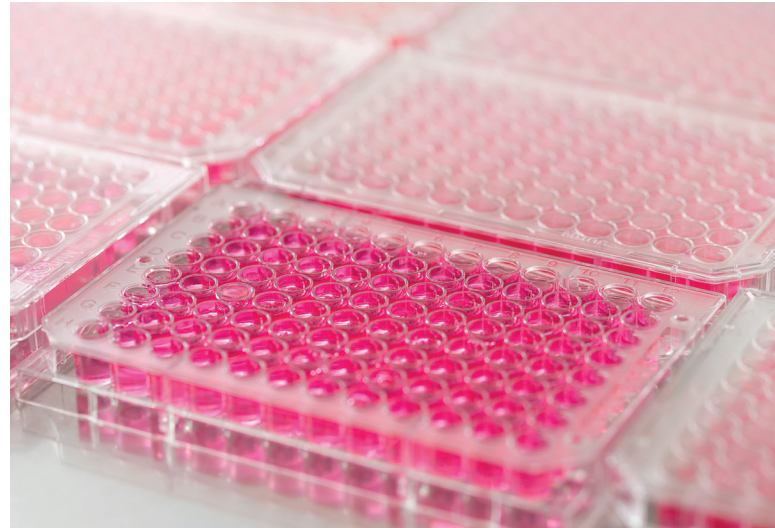


### **Quality controls**

- **Physico-chemical**  
Purity, Quantity, Integrity, Sterility, ...
- **Technical**  
Method validation
- **Functional**  
Binding, cytotoxicity, ADCC, CDC, residual endotoxins, *in vivo* testing...

## Additional services for antibodies

- Culture media selection and serum-free production,
- Isotyping, sequencing, cell-line adaptation, cell banking, productivity assessment, variable chains sequencing, cloning for recombinant antibody generation.
- ELISA assays development.
- Mycoplasma testing and cell line decontamination.



## Hybridoma generation

We provide a comprehensive **murine hybridomas development offer**. We generate custom hybridomas from a large panel of antigens, and provide a personalized follow up at every step, from antigen design up to the cell banking.

## Why work with Eurogentec?

- Belgian company with state-of-the-art facilities.
- Over 25 years of experience in antibody manufacturing.
- A Team of antibody experts, dedicated to find the best strategy for the customer's request.



Learn more on [www.eurogentec.com](http://www.eurogentec.com)