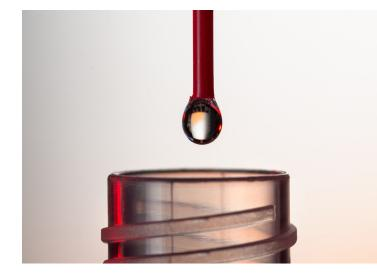


Large scale antibody production

We offer <mark>large</mark> scale antibody solutions for the industry.

We produce and purify both polyclonal and monoclonal custom antibodies with flexibility. Our team of experts is ready to handle your request and provide you with the most suited deliverable.



eurogentec.com

Polyclonal antibody production

We produce and deliver serum or purified polyclonal antibodies. We handle large demanding production projects, according to our own SOP's or we can adapt our processes to be in line with the customer requirements.



🕆 Large volume

We can produce up to several liters of serum from rabbits or goats, in a single batch.



Long-term programs

We can host immunized animals to generate large serum volume over time, delivered at your convenience.



Purified antibodies

We can purify grams of custom antibodies, using a tailored set of purification strategies.

In vitro monoclonal antibody production

We produce monoclonal antibodies *in vitro* from hybridomas. Depending on the customer's request and the production scheme, we can generate up to 100 g of high-quality monoclonal antibodies.



In vitro production

All our custom monoclonal antibodies are produced using an *in vitro* method.



We are equipped to produce from mg to grams purified monoclonal antibodies. Test studies are performed for request from 200 mg.



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



Why partner with Eurogentec?

- A dedicated **project manager** follows up throughout the production process
- We adjust the immunization and purification programs according to the customer's needs with remarkable **flexibility**.
- We have **great expertise** in the development and production of antibodies for over 25 years.
- Large animal **housing capacity** & downstream processes

Quality

- Comprehensive Quality Management System (QMS)
- Belgian state-of-the-art facilities
- 3Rs guidelines for Animal Welfare
- GMP BioManufacturing facilities for the production of antibody fragments by biofermentation.





Res-FLYER-EN-Large-scale-antibody-Oct-2021-VI



Learn more on www.eurogentec.com