

C U S T O M Monoclonal A N T I B O D I E S

MONOCLONAL ANTIBODY PRODUCTION

Consultation on epitope selection

Evaluation of suitable screening methods

Frequent interactions between researcher and AnaSpec in all phases of development

Project segmentation into different phases to provide researcher flexibility for terminating project at the end of each phase

Phase I: Immunogen Design and Immunization (9 wks)

Phase II: Cell Fusion and Parental Clone Screening (3 wks)

Phase III: Monoclonal Selection and Cryogenic Vial Preservation (3 wks)

Phase IV: Large Scale Antibody Production (3-6 wks)

For additional information about our custom antibody production services, please contact ab@anaspec.com. AnaSpec provides both small and large scale monoclonal antibody production.

ADDED VALUES TO ANASPEC'S CUSTOM ANITBODY SERVICE:

1. High purity peptides
2. Expert consultation on epitope selection
3. Proprietary QC methods for KLH and BSA peptide conjugations
4. Large scale antibody production, in-vivo or in-vitro
5. Effective affinity purification of phospho-antibodies
6. Online tracking of antibody projects
7. Literature citations of monoclonal antibodies generated by AnaSpec
8. Customers' references available



ISO 9001:2008 Company

To search for a list of custom antibodies made by AnaSpec, please visit www.highwire.org and type in AnaSpec in the keyword search box.

