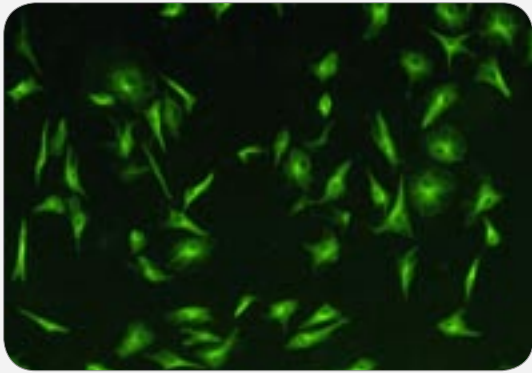


Polyclonal

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



Regular or Phospho-Polyclonal Antibodies (2 rabbits)

56-day standard protocol

5 mL pre-immune serum per rabbit

50 mL anti-serum per rabbit

Additional Services:

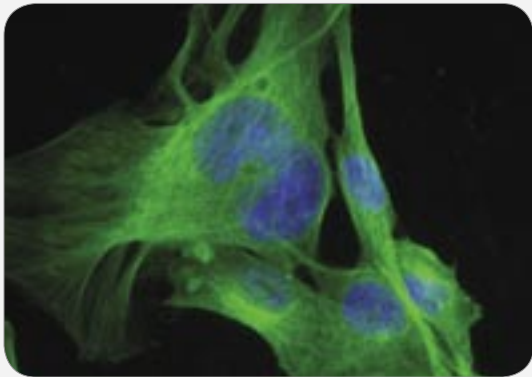
ELISA

Peptide Synthesis - >80% purity (non-phosphopeptide),
>90% purity (phosphopeptide), HPLC & Mass Spec certified

Peptide-Carrier Protein Conjugation - Conjugation to BSA, KLH, others

Affinity Purification

Antibody Conjugation to Fluorophores



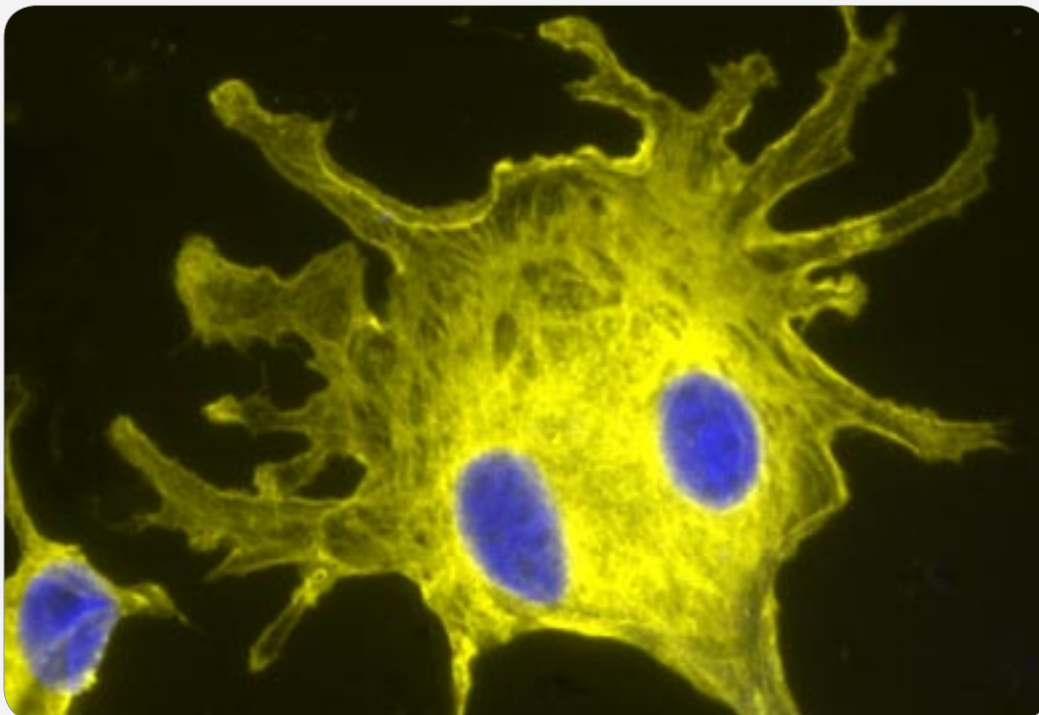
Complete Anti-Peptide Antibody Package includes:

Peptide Synthesis (up to 15 residues)

Antibody Production (first & second bleeds)

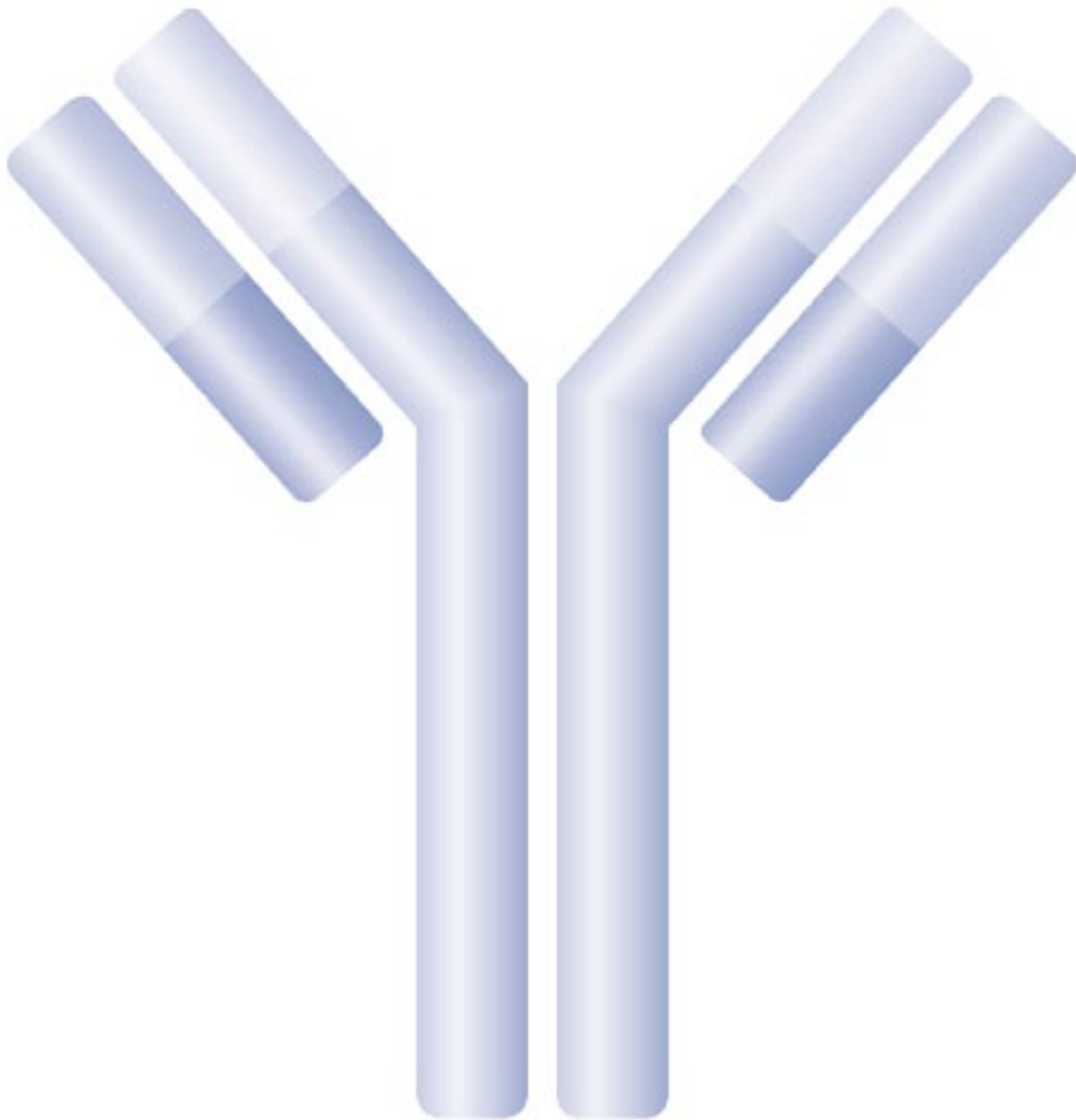
Peptide Conjugation

ELISA Titration



ISO 9001:2000 Company

Added Values to AnaSpec's Custom Antibody Service:



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

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.