

Added Values to AnaSpec's Custom Synthesis Service:

1. Pre-synthesis sequence examination
2. Free consultation on design of peptides
3. Expert experience in synthesis of unusual peptides
4. Quality Assurance
 - a) ISO 9001-2000 Certified
 - b) In-process and final QC. Peptides are checked three times
by mass spec and HPLC
- 5) On-line status checking of synthesis progress
- 6) Dedicated customer support staff to maintain close communication with customers
- 7) Prompt delivery
- 8) Computerized record keeping

