

# GMP Peptides

15 YEARS OF DISCOVERY  
15  
ANASPEC

With 15 years of peptide synthesis experience, AnaSpec Inc, is committed to meet or exceed the quality standards of GMP, and ISO 9001:2000



■ Cyclic peptides

■ Dye-Labeled peptides

■ FRET Probes

■ Heavy Isotope Labeled

■ Stapled peptides

■ Phosphopeptides

■ Biotinylated peptides

■ Pegylated peptides

■ Unusual Modifications

# Let AnaSpec take you from research to Bulk API production

Solid Phase Synthesis

HPLC Purification

Lyophilization

Cyclic peptides, Dye-labeled peptides, FRET probes, Phospho-peptides, Heavy isotope labeled, Stapled peptides



- 5,000 sq. ft. GMP Facility on-site located in San Jose, CA
- Class 10,000 Cleanroom
- Milligram to Kilogram Scale
- Custom or Catalog sequences
- Customizable product specifications

- Bulk API Manufacturing
- QC Product Release Testing
- High Purity >98%
- Stability Studies
- Full Documentation for Drug Master File
- Excellent Technical Support

AnaSpec holds a California Drug Manufacturing License (License# 47174) and a FDA Registration for GMP Drug Manufacturing (Registration # 2951078), ISO 9001:2000



Delivery schedule dependent upon the specification.  
For inquiries on custom quotations please email [service@anaspec.com](mailto:service@anaspec.com)