



CUSTOM Peptides FOR HTS

AnaSpec, Inc., with years of peptide synthesis experience has now developed a new technology to synthesize peptides on microtiter plate. This new technology significantly reduces the cost of peptide synthesis at small scale level. AnaSpec now offers this custom peptide service:

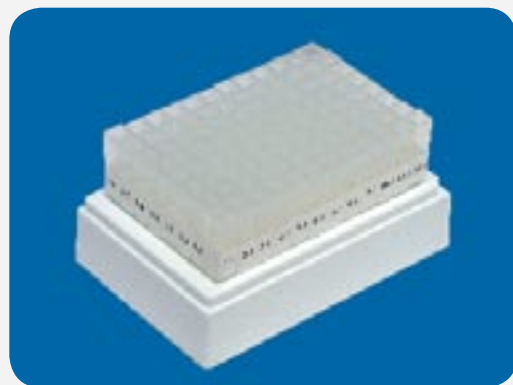
1. Up to 15 mer
2. 5-10 mg unpurified peptide
3. Amide at C-terminus
4. 4 HPLC analyses per plate
5. 48 peptides minimum order
6. 3-4 weeks delivery

The following options are also available:

1. Peptide length > 15 mer
2. 10-15 mg unpurified peptide
3. Purified peptide
4. Other C-terminus modifications
5. HPLC analysis on all peptides
6. Modified peptides - dye labeled, biotinylated, etc.



ISO 9001:2008 Company



Peptides on a 96-well plate.