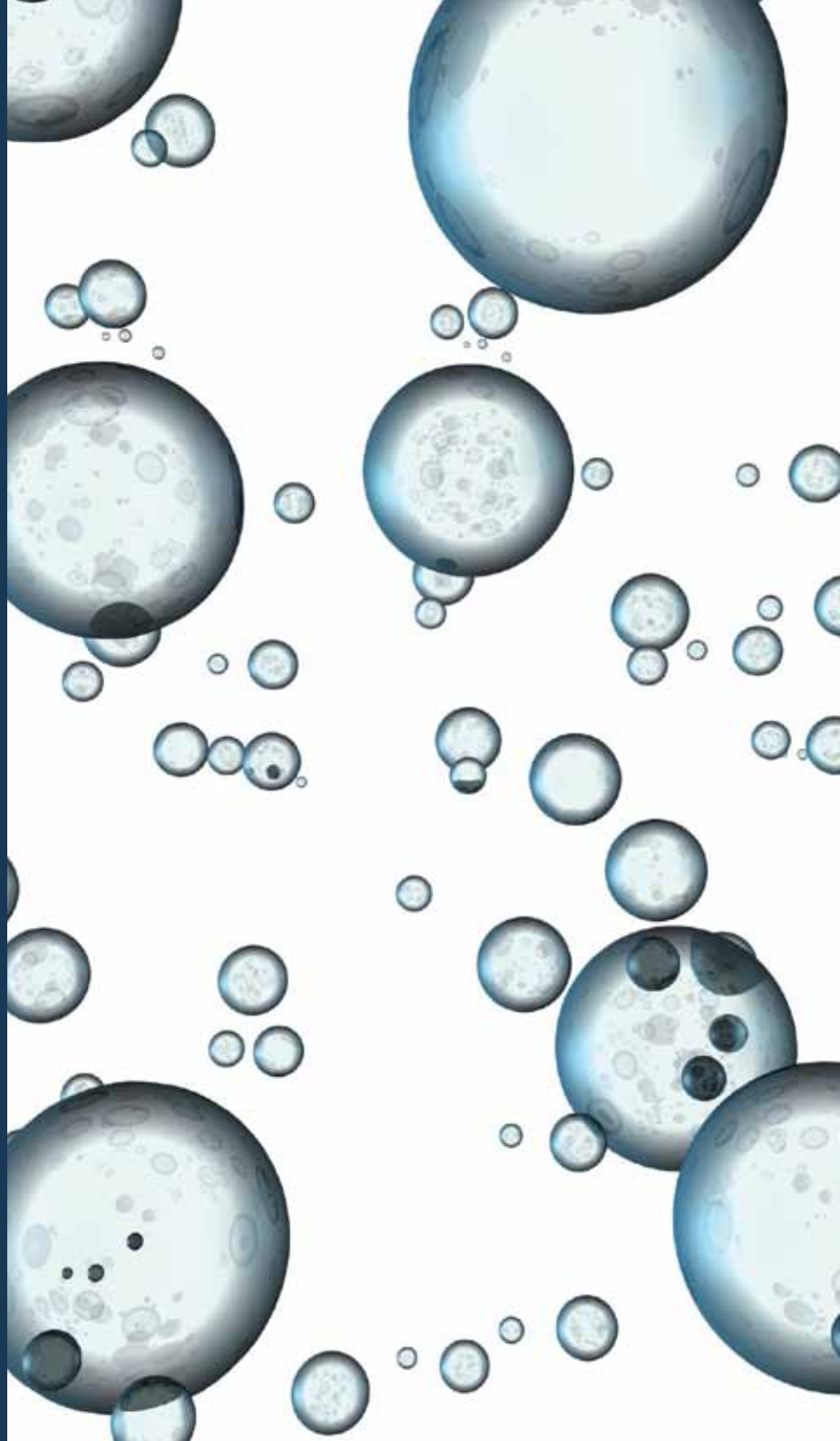


3

3.10 Peptide Synthesis Reagents



Peptide Synthesis Reagents

Eurogentec and AnaSpec, Inc. are leading providers of integrated proteomics solutions for worldwide life science research. With a vision for innovation through synergy, we offer expertise in three primary technologies: peptides, detection reagents, and combinatorial chemistry.

This triple-platform approach allows AnaSpec to provide research solutions that are unavailable to single-platform companies.

We carry a broad product line of biochemicals and reagents for basic research, high-throughput screening and drug discovery.

Proteomic Synthesis Reagents Products that can be discovered on www.eurogentec.com

Resins

- ▣ 2-Chloriotrityl Resins
- ▣ Trityl Resins
- ▣ Fmoc-Rink Amide MBHA Resins
- ▣ HMP Resins
- ▣ MAPS Resins
- ▣ PAM Resins
- ▣ Tentagel Resins
- ▣ Wang Resins
- ▣ AFC, AMC and pNA Resins
- ▣ Seber Resins
- ▣ Miscellaneous Resins

Reagents

- ▣ Coupling Reagents and Additives
- ▣ Protecting Reagents
- ▣ Labeling Reagents
- ▣ Other Peptide Synthesis Reagents
- ▣ Building Blocks

Unusual Amino Acids

- ▣ Analogues of Benzoic Acid
- ▣ Analogues of Phenylalanine and Tyrosine
- ▣ Miscellaneous Aromatic Amino Acids
- ▣ Analogues of Proline
- ▣ Analogues of Tryptophan
- ▣ Miscellaneous Heterocyclic Amino Acids
- ▣ Analogues of Arginine and Lysine
- ▣ Analogues of Alanine, Glycine, Valine and Leucine
- ▣ Analogues of Aspartic and Glutamic Acids
- ▣ Analogues of Cysteine and Methionine
- ▣ Analogues of Serine, Threonine and Statine
- ▣ Miscellaneous Aliphatic Amino Acids
- ▣ β -Amino Acids
- ▣ N- α -Methyl Amino Acids

Amino Alcohols

Amino Acid Cartridges

- ▣ Fmoc-L-Amino Acid Cartridges for ABI Synthesizers (1 mmol)
- ▣ Fmoc-D-Amino Acid Cartridges for ABI Synthesizers (1 mmol)
- ▣ Boc-L-Amino Acid Cartridges for ABI Synthesizers (1 mmol)
- ▣ Boc-D-Amino Acid Cartridges for ABI Synthesizers (1 mmol)
- ▣ Boc-L-Am no Acid Cartridges for ABI Synthesizers (2 mmol)
- ▣ Miscellaneous Amino Acid Cartridges for ABI Synthesizers (1 mmol)
- ▣ Fmoc-L-Amino Acid Cartridges for Protein Technologies' Symphony Peptide Synthesizers

Accessories

- ▣ Filtration Accessories
- ▣ Reaction Vessels and Accessories
- ▣ Miscellaneous Accessories

Reaction Vessel



Our valued customers have chosen AnaSpec/Eurogentec as provider for

Description	Reference	Name	Size
Standard Amino Acids	AS-20141-5	Fmoc - Arg(Pbf) - OH	500 g
Synthesis Reagent	AS-25185-XX	DTT	XX g
Resins	AS-60098	Fmoc - Leu - ethylamide - Sieber resin	1 g
Resins	AS-20084	Fmoc-Rink Amide MBHA resin	10 g
Unusual Amino Acids	AS-64175-1000	(R) - N - Fmoc - 2 - (7' - octenyl) alanine	1 g