



integrated proteomics solutions.

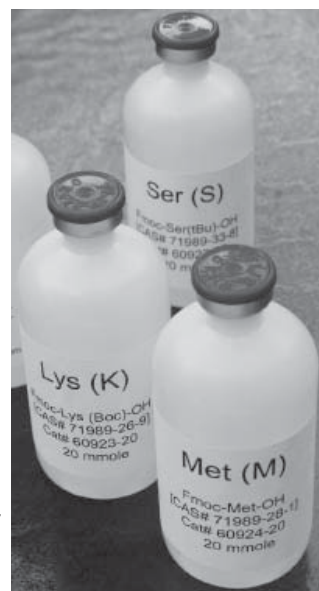
SYNTHESIS ACCESSORIES

AnaSpec, Inc. • 2149 O'Toole Ave. • San Jose, CA 95131 • 800-452-5530 • Fax 408-452-5059 • service@anaspec.com • www.anaspec.com

Amino acid cartridges



For ABI 433A & 431A



For Protein Technology Symphony Synthesizer

Reaction Vessels for Protein Technology Symphony



Cat # 60879 (0.25mmol)
Reaction vessel for use with Rainin Symphony peptide synthesizer.
100 /pk

Cat # 60911 (0.5 mmol) AnaSpec reaction vessel designed for use with Rainin Symphony peptide synthesizer. The top of this reaction vessel easily screws to the bottom to ensure a tight fit.
100 /pk

Filter Paper for ABI 433A & 431A



Product Name	Size	Catalog#
Filter Paper, 0.312" For use in stainless steel in-line filter paper	100 /box	10812
Filter Paper, 1.08", tabbed For use with ABI reaction vessels	25 /box	10814
Filter Paper, 1.1" For use with ABI reaction vessel 0.1 mmol synthesis	25 /box	10813
Filter Paper, 1.25" For use with ABI 40-60 mL reaction vessel 0.25 mmol-1 mmol synthesis.	25 /box	10310