

## Amino acid cartridges for ABI 433A & 431A



## Reaction Vessels for Protein Technology's Symphony



Cat # 60879 (0.25mmol)  
Reaction vessel for use with Rainin Symphony peptide synthesizer.  
100 /pk retail \$540

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 retail \$750

## Filter Paper for ABI 433A & 431A



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