

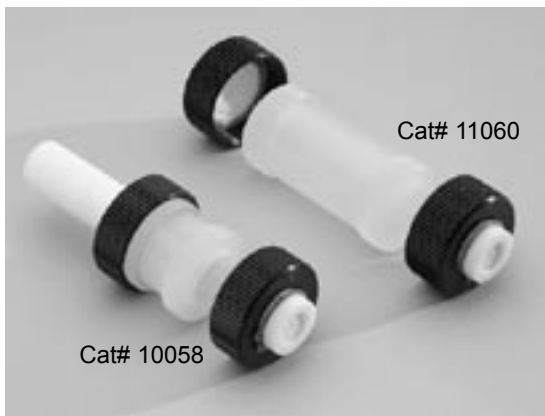


integrated proteomics solutions.

REACTION VESSELS

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

AnaSpec's Reaction Vessels for Use in ABI 433A & 431A Synthesizers



0.10 mmol Reaction Vessel, cat# 10058

0.25 mmol Reaction Vessel, cat# 11059

0.50 mmol Reaction Vessel, cat# 11060

Reaction Vessels for Protein Technology's Symphony



Cat # 60879 (0.25mmol) Cat # 60911 (0.5 mmol)



Cat # 62724 (0.25mmol) Cat # 62725 (0.5 mmol)

Reaction Vessel + Cleavage Unit - **NEW**

Reaction Vessels for Manual Synthesis



15 mL Reaction Vessel, Manual Synthesis, Cat# 23565

25 mL Reaction Vessel, Manual Synthesis, Cat# 23564

80 mL Reaction Vessel, Manual Synthesis, Cat# 20371

200 mL Reaction Vessel, Manual Synthesis, Cat# 20372

500 mL Reaction Vessel, Manual Synthesis, Cat# 25184

1000 mL Reaction Vessel, Manual Synthesis, Cat# 20373