

Ordering Information for the Biotechnology Explorer™ Program



Product Name	Catalog Number	AP Biology	Bioinformatics	Biotechnology	Chemistry/Physics	Food and Agricultural Science	Forensic Science	General Biology	Genetics	Health Science and Medicine	Home Science	Immunology	Microbiology	Middle School	Molecular Biology or DNA Analysis	Protein Analysis	STEM
<i>C. elegans</i> Behavior Kit	166-5120EDU	•	•	•				•	•				•		•		•
pGLO™ Bacterial Transformation Kit	166-0003EDU	•		•				•					•		•		•
Green Fluorescent Protein Chromatography Kit	166-0005EDU			•	•			•					•		•	•	•
pGLO Kit SDS-PAGE Extension	166-0013EDU			•	•			•					•			•	•
Secrets of the Rainforest™ Kit	166-0006EDU			•				•		•			•		•	•	•
Microbes and Health Kit	166-5030EDU			•		•		•		•			•	•			•
Biofuel Enzyme Kit	166-5035EDU	•		•	•	•		•								•	•
Size Exclusion Chromatography Kit	166-0008EDU			•	•									•		•	•
Got Protein?™ Kit	166-2900EDU			•	•	•		•								•	•
ELISA Immuno Explorer™ Kit	166-2400EDU			•		•		•		•		•		•		•	•
Comparative Proteomics Kit I: Protein Profiler Module	166-2700EDU	•	•	•												•	•
Comparative Proteomics Kit II: Western Blot Module	166-2800EDU	•		•								•				•	•
Comparative Proteomics Kits I and II: Protein Profiler and Western Blot Modules	166-2850EDU	•	•	•								•				•	•
Rapid Blotting and V3 Western Workflow™ Starter Kit	166-2875EDU	•	•	•								•				•	•
DNA Model Kit	166-7015EDU			•				•	•					•	•		•
Genes in a Bottle™ Kit	166-2300EDU			•				•	•					•	•		•
Forensic DNA Fingerprinting Kit	166-0007EDU	•		•			•	•	•						•		•
Analysis of Precut Lambda DNA Kit	166-0001EDU	•		•				•							•		•
STEM Electrophoresis Classroom Kit (includes IDEA kit reagent refill pack)	166-5090EDU			•		•		•						•			•
Crime Scene Investigator PCR Basics™ Kit	166-2600EDU			•			•	•	•						•		•
PV92 PCR Informatics Kit	166-2100EDU	•	•	•					•						•		•
Fish DNA Barcoding Kit	166-5100EDU		•	•		•			•						•		•
GMO Investigator™ Kit	166-2500EDU			•		•			•						•		•
Crime Scene Investigator PCR Basics Real-Time PCR Starter Kit	166-2660EDU			•			•		•						•		•
GMO Investigator Real-Time PCR Starter Kit	166-2560EDU			•		•			•						•		•
Cloning and Sequencing Explorer Series (includes eight modules)	166-5000EDU	•	•	•		•			•				•		•	•	•
Protein Expression and Purification Series — Centrifugation Purification Process (includes four modules)	166-5040EDU			•	•					•		•	•		•	•	•
Genes in a Bottle — DNA Necklace Kit	Available to order on www.amazon.com										•						
DNA Model Kit	Available to order on www.amazon.com										•						
The Candy Caper — A Crime Scene Detective Kit	Available to order on www.amazon.com										•						