

Map of AP-Aligned Bio-Rad Kits with Investigative Laboratories



Cover more than one AP Biology Big Idea with these AP-aligned Bio-Rad kits.

Big Idea 1								Big Idea 2					Big Idea 3						Big Idea 4									
	pGLO™ Bacterial Transformation Kit	pGLO Transformation and Inquiry Kit for AP Biology	Biofuel Enzyme Kit	PV92 PCR Informatics Kit	Comparative Proteomics Kit I: Protein Profiler Module	Fish DNA Barcoding Kit	Cloning and Sequencing Explorer Series		pGLO Bacterial Transformation Kit	pGLO Transformation and Inquiry Kit for AP Biology	Biofuel Enzyme Kit	Cloning and Sequencing Explorer Series	<i>C. elegans</i> Behavior Kit		pGLO Bacterial Transformation Kit	pGLO Transformation and Inquiry Kit for AP Biology	PV92 PCR Informatics Kit	Comparative Proteomics Kit I: Protein Profiler Module	Cloning and Sequencing Explorer Series	Forensic DNA Fingerprinting Kit/ Lambda DNA Kits		pGLO Bacterial Transformation Kit	pGLO Transformation and Inquiry Kit for AP Biology	Biofuel Enzyme Kit	Comparative Proteomics Kit I: Protein Profiler Module	<i>C. elegans</i> Behavior Kit	Forensic DNA Fingerprinting Kit/ Lambda DNA Kits	
LAB 1								LAB 4						LAB 7								LAB 10						
LAB 2								LAB 5						LAB 8								LAB 11						
LAB 3								LAB 6						LAB 9								LAB 12						
																						LAB 13						
RELATED LABS								RELATED LABS						RELATED LABS							RELATED LABS							
														Genes in a Bottle™ Kit (166-2300EDU)						Protein Expression and Purification Series (166-5040EDU, 166-5045EDU, 166-5050EDU, and 166-5070EDU)								
														Crime Scene Investigator PCR Basics™ Kits (166-2600EDU and 166-2650EDU)														
														pGLO Kit SDS-PAGE Extension (166-0013EDU)														
														IDEA Kits (166-5075EDU and 166-5077EDU)														

- pGLO Bacterial Transformation Kit Catalog #166-0003EDU
- Biofuel Enzyme Kit Catalog #166-5035EDU
- PV92 PCR Informatics Kit Catalog #166-2100EDU
- Comparative Proteomics Kit I: Protein Profiler Module Catalog #166-2700EDU
- Fish DNA Barcoding Kit Catalog #166-5100EDU
- Cloning and Sequencing Explorer Series Catalog #166-5000EDU
- *C. elegans* Behavior Kit Catalog #166-5120EDU
- Forensic DNA Fingerprinting Kit Catalog #166-0007EDU
- pGLO Transformation and Inquiry Kit for AP Biology Catalog #166-0335EDU
- Analysis of Precut Lambda DNA Kit Catalog #166-0001EDU
- Restriction Digestion and Analysis of Lambda DNA Kit Catalog #166-0002EDU