

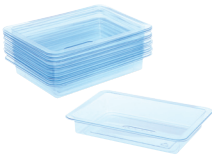






Reduce. Reuse. Refresh.

Refresh kit components reduce packaging waste, reuse components, and refresh your kits, and you'll save storage space by purchasing individual items. Bio-Rad now has many individual components for Biotechnology Explorer kits available for purchase.

	pGLO Bacterial Transformation Kit	Microbes and Health Kit/Microbial Culturing Module	Genes in a Bottle Kit	Analysis of Precut Lambda DNA Kit	Restriction Digestion and Analysis of Lambda DNA Kit	Forensic DNA Fingerprinting Kit	Crime Scene Investigator PCR Basics Kit	PV92 PCR Informatics Kit	GMO Investigator Kit	Description	Catalog No.
 <p>Petri Dishes</p>	•									pGLO Bacterial Transformation Kit Refill Package	166-0555EDU
	•									pGLO Plasmid, 20 µg, lyophilized	166-0405EDU
	•									Arabinose, 600 mg, lyophilized	166-0406EDU
	•	•								Ampicillin, 30 mg, lyophilized	166-0407EDU
	•	•								<i>E. coli</i> Strain HB101 K-12, lyophilized, 1	166-0408EDU
	•									Transformation Solution, 15 ml	166-0409EDU
 <p>Inoculation Loops</p>	•									LB Broth, 10 ml	166-0421EDU
	•	•								LB Nutrient Agar Powder, 20 g, makes forty 60 mm agar plates	166-0600EDU
	•	•								LB Nutrient Agar Powder, 500 g, makes one thousand 60 mm agar plates	166-0472EDU
	•	•								Petri Dishes, 60 mm, sterile, 500	166-0470EDU
 <p>Gel Staining Trays</p>	•	•								Inoculation Loops, 10 µl, sterile, 80	166-0471EDU
	•			•	•	•	•	•	•	Jellyfish Foam Floating Racks, 8 racks, 12 microcentrifuge tube wells	166-0479EDU
	•									Long-Wave UV Lamp, 1	166-0500EDU
	•									Long-Wave UV Pen Light, 1	166-0530EDU
 <p>Disposable Plastic Transfer Pipets</p>				•	•	•	•	•	•	TAE, 50x, 1 L	161-0743EDU
				•	•	•	•	•	•	TAE, 50x, 5 L cube	161-0773EDU
				•	•	•	•	•	•	Sample Loading Dye, 5x, 10 ml	161-0767EDU
				•	•	•	•	•	•	Certified Molecular Biology Agarose, 25 g	161-3100EDU
				•	•	•	•	•	•	Certified Molecular Biology Agarose, 125 g	161-3101EDU
				•	•	•	•	•	•	Certified Molecular Biology Agarose, 500 g	161-3102EDU
				•	•	•	•	•	•	Fast Blast DNA Stain, 500x, 100 ml	166-0420EDU
				•	•	•	•	•	•	Small Fast Blast DNA Stain Electrophoresis Reagent Pack	166-0450EDU
				•	•	•	•	•	•	Medium Fast Blast DNA Stain Electrophoresis Reagent Pack	166-0455EDU
				•	•	•	•	•	•	Large Fast Blast DNA Stain Electrophoresis Reagent Pack	166-0460EDU
 <p>1.5 ml Microcentrifuge Tubes</p>			•							Genes in a Bottle DNA Extraction Refill Package	166-2001EDU
			•							Lysis Buffer, 150 ml	166-2002EDU
	•	•					•	•	•	MasterMix for PCR, 2x, 90 units, 1.2 ml	166-5009EDU
	•								•	Disposable Plastic Transfer Pipets, sterile, 500	166-0474EDU
 <p>Conical Centrifuge Tubes</p>			•							Disposable Plastic Transfer Pipets, nonsterile, 500	166-0480EDU
	•									Colored 1.5 ml Microcentrifuge Tubes, 6 colors, 600	166-0473EDU
	•			•	•	•	•	•	•	EZ Micro Test Tubes, 1.5 ml, natural, 500	223-9480EDU
	•									EZ Micro Test Tubes, 2 ml, natural, 500	223-9430EDU
									•	Conical Tubes, 1.5 ml, with installed O-ring screwcaps, sterile, 500	224-0110EDU
 <p>Cell Culture Tubes</p>			•							Conical Centrifuge Tubes, 15 ml, 50	166-0475EDU
		•								Cell Culture Tubes, 17 x 100, 14 ml, sterile, 25	166-0476EDU
									•	0.2 ml Tubes With Domed Caps, natural, 1000	TWI-0201EDU
									•	PCR Tube Capless Adapters, 500	223-9500EDU
									•	96-Place PCR Tube Rack and Cover, 5	TRC-0501EDU