

Choose Your Competent Cells With Confidence

Difficult Cloning Competent Cells*

Large or Litigated DNA		
AGL Cat. #	Product	Size
200314	XL10-Gold Ultracompetent	≥5 x 10 ⁹
200315	XL10-Gold Ultracompetent	≥5 x 10 ⁹
200317	XL10-Gold Kan ^r Ultracompetent	≥5 x 10 ⁹
200159	ElectroTen-Blue Electroporation	≥5 x 10 ¹⁰

Unstable Cloning		
AGL Cat. #	Product	Size
200152	SURE 2 Supercompetent	≥1 x 10 ⁹
200238	SURE	≥5 x 10 ⁸
200227	SURE Electroporation	≥1 x 10 ¹⁰

Toxic Cloning 4-Fold Reduction of Copy Number		
AGL Cat. #	Product	Size
200171	ABLE C	5 x 10 ⁸ - 2 x 10 ⁷

10-Fold Reduction of Copy Number		
AGL Cat. #	Product	Size
200172	ABLE K	5 x 10 ⁸ - 2 x 10 ⁷

Genomic DNA or Methylated cDNA		
AGL Cat. #	Product	Size
200151	XL2-Blue MRF ⁺ Ultracompetent	≥5 x 10 ⁹
200230	XL1-Blue MRF ⁺ Supercompetent	≥1 x 10 ⁹
200248	XL1-Blue MRF ⁺ Kan ^r Supercompetent	≥1 x 10 ⁹
200229	XL1-Blue MR Supercompetent	≥1 x 10 ⁹

Phage Display		
AGL Cat. #	Product	Size
200123	TG1 Electroporation	≥1 x 10 ¹⁰

Difficult Cloning Pack		
AGL Cat. #	Product	
	XL10-Gold Kan ^r Ultracompetent	
230247	SURE 2 Supercompetent	
	ABLE K	

ABLE Kit	
AGL Cat. #	Product
	ABLE C
200170	ABLE K

ABLE Electroporation Kit	
AGL Cat. #	Product
	ABLE C Electroporation
200160	ABLE K Electroporation

General Cloning Competent Cells*

Routine Cloning		
AGL Cat. #	Product	Size
200150	XL2-Blue Ultracompetent	≥5 x 10 ⁹
200236	XL1-Blue Supercompetent	≥1 x 10 ⁹
200228	XL1-Blue Electroporation	≥1 x 10 ¹⁰
200249	XL1-Blue	≥1 x 10 ⁸
200130	XL1-Blue Subcloning Grade	≥1 x 10 ⁶

Conveniently Packaged		
AGL Cat. #	Product	Size
230350	SoloPack Gold Supercompetent	≥1 x 10 ⁹
230325	SoloPack Gold	≥1 x 10 ⁸
200324	96Pack Gold	≥1 x 10 ⁸

Generate Unmethylated DNA		
AGL Cat. #	Product	Size
200247	SCS110	≥5 x 10 ⁶
200239	JM110	≥5 x 10 ⁶

Random Mutagenesis		
AGL Cat. #	Product	Size
200129	XL1-Red	≥1 x 10 ⁶

Random Mutagenesis		
AGL Cat. #	Product	Size
200231	SCS1 Supercompetent	≥1 x 10 ⁹
200232	AG1	≥1 x 10 ⁸
200233	NM522	≥1 x 10 ⁸
200234	JM101	≥1 x 10 ⁸
200235	JM109	≥1 x 10 ⁸

Difficult Cloning Pack		
AGL Cat. #	Product	
	XL2-Blue Ultracompetent	
230248	XL2-Blue MRF ⁺ Ultracompetent	
	SoloPack Gold	

Protein Expression Competent Cells*

General Protein Expression		
AGL Cat. #	Product	Size
230132	BL21-Gold (DE3)	≥1 x 10 ⁸
200131	BL21(DE3)	≥1 x 10 ⁶

Codon Bias Universal Strain		
AGL Cat. #	Product	Size
230280	BL21-CodonPlus (DE3)-RIPL	≥1 x 10 ⁶

Conveniently Packaged		
AGL Cat. #	Product	Size
230350	SoloPack Gold Supercompetent	≥1 x 10 ⁹
230325	SoloPack Gold	≥1 x 10 ⁸
200324	96Pack Gold	≥1 x 10 ⁸

High GC Content		
AGL Cat. #	Product	Size
230255	BL21-CodonPlus (DE3)-RP	≥1 x 10 ⁷
230275	BL21-CodonPlus (DE3)-RP-X	≥1 x 10 ⁷
230250	BL21-CodonPlus-RP	≥1 x 10 ⁷

High AT Content		
AGL Cat. #	Product	Size
230245	BL21-CodonPlus (DE3)-RIL	≥1 x 10 ⁷
230265	BL21-CodonPlus (DE3)-RIL-X	≥1 x 10 ⁷
230240	BL21-CodonPlus-RIL	≥1 x 10 ⁷

Toxic Proteins		
AGL Cat. #	Product	Size
230134	BL21-Gold (DE3) pLysS	≥1 x 10 ⁸
200132	BL21(DE3) pLysS	≥1 x 10 ⁶
230130	BL21-Gold	≥1 x 10 ⁸
200133	BL21	≥1 x 10 ⁶

Solubility & Codon Bias		
AGL Cat. #	Product	Size
230193	ArcticExpress (DE3)-RIL	≥5 x 10 ⁶
230194	ArcticExpress (DE3)-RP	≥5 x 10 ⁶
230195	ArcticExpress RIL	≥5 x 10 ⁶
230196	ArcticExpress RP	≥5 x 10 ⁶

Solubility		
AGL Cat. #	Product	Size
230191	ArcticExpress	≥5 x 10 ⁶
230192	ArcticExpress (DE3)	≥5 x 10 ⁶

Generate Phosphorylated Proteins		
AGL Cat. #	Product	Size
200134	TKB1	≥5 x 10 ⁵

Protein Expression Pack		
AGL Cat. #	Product	Size
	BL21 (DE3)	≥5 x 10 ⁷
230246	BL21-Gold (DE3)	≥5 x 10 ⁵
	BL21-CodonPlus (DE3)-RIPL	
	BL21-Gold (DE3) pLysS	

