

Catalog Number	Description
200123	TG1 Electro-Comp Cells 5 x 0.1 ml
200124	TKX1 Competent Cells
200129	XL1-Red Competent Cells 5 x 0.2 ml
200130	XL1-Blue Subclon Grade Comp Cell 8x0.5ml
200131	BL21(DE3) Competent Cells
200132	BL21(DE3) pLysS Comp Cells
200133	BL21 Competent Cells
200134	TKB1 Competent Cells
200138	XL-1 Blue MRF ⁺ Kan Libr Pack Comp Cells
200150	XL2-Blue UltraComp Cells 10 x 0.1 ml
200151	XL2-Blue MRF UltraComp Cells 10 x 0.1ml
200152	SURE 2 SuperComp Cells 10 x 0.1 ml
200154	BJ5183 Electroporation-Competent Cells
200157	BJ5183-AD-1 Electroporation Comp Cells
200158	XL1-Blue MRF Electro-Comp Cells 5x0.1ml
200159	ElectroTen-Blue Electro-Comp cells
200160	Able Electro-Comp Cells 5 x 0.1 ml
200161	Able C Electro-Comp Cells 5 x 0.1 ml
200162	Able K Electro-Comp Cells 5 x 0.1 ml
200170	Able Comp Cells(Able C and K) 5 x 0.2ml
200171	Able C Competent Cells 5 x 0.2 ml
200172	Able K Competent Cells 5 x 0.2 ml
200192	BacterioMatch II Vldtn Rprtr Comp Cell
200201	Gigapack III Gold Packaging Extract,4rxn
200202	Gigapack III Gold Packaging Extract,7rxn
200203	Gigapack III Gold Packaging Extract11rxn
200204	Gigapack III Plus Packaging Extract,4rxn
200205	Gigapack III Plus Packaging Extract,7rxn
200206	Gigapack III Plus Packaging Extract11rxn
200207	Gigapack III XL Packaging Extract, 4 rxn
200208	Gigapack III XL Packaging Extract, 7 rxn
200209	Gigapack III XL Packaging Extract,11 rxn
200227	SURE Electro-Comp Cells 5 x 0.1 ml
200228	XL1-Blue Electro-Comp Cells 5 x 0.1 ml
200229	XL1-Blue MR SuperComp Cells 5 x 0.2 ml
200230	XL1-Blue MRF SuperComp Cells 5 x 0.2 ml
200231	SCSI SuperCompetent Cells 5 x 0.2 ml
200232	AG1 Competent Cells 5 x 0.2 ml
200233	NM522 Competent Cells 5 x 0.2 ml
200234	JM101 Competent Cells 5 x 0.2 ml
200235	JM109 Competent Cells 5 x 0.2 ml
200236	XL1-Blue SuperCompetent Cells 5 x 0.2 ml
200238	SURE Competent Cells 5 x 0.2 ml
200239	JM110 Competent Cells 5 x 0.2 ml
200247	SCS110 Competent Cells 5 x 0.2 ml
200248	XL1-Blue MRF Kan SuperComp Cell 5x0.2ml

200249	XL1-Blue Competent Cells 5 x 0.2 ml
200285	Mammalian Transfection Kit
200314	XL10-Gold Competent Cells, 5 x 100 ul
200315	XL10-Gold Competent Cells, 10 x 100 ul
200317	XL10-Gold Kanr Comp Cells 10 x 0.1 ml
200324	96Pack Gold Competent Cells
200385	Mammalian Transfection Kit-Calcium Phosphate
200388	Transfection MBS Mammalian Kit
200488	ViraPack Transfection Kit
200513	QuikChange Multi Site-Directed Mutagenesis (Commercial), 30 rxn
200514	QuikChange Multi Site-Directed Mutagenesis (Academic), 30 rxn
200515	QuikChange Multi Site-Directed Mutagenesis (Academic), 10 rxn
200516	QuikChange XL Site-Directed 30 rxn
200517	QuikChange XL Site-Directed 10 rxn
200518	QuikChange Site-Directed 30 rxn
200519	QuikChange Site-Directed 10 rxn
200521	QuikChange II XL site-directed 10rx
200522	QuikChange II XL site-directed 30rx
200523	QuikChange II site-directed 10 rxn
200524	QuikChange II site-directed 30 rxn
200531	QuikChange Multi Site-Directed Mutagenesis (Commercial), 10 rxn
200550	GeneMorph II Random Mutagenesis
200552	GeneMorph II EZClone Domain Mutagenesis
200555	QuikChange II-E Site-Directed Mutagenesis, 10 rxns
204130	GeneJammer Transfection Reagent
204131	GeneJammer Transfection Reagent
204132	GeneJammer Transfection Reagent
210513	QC Lightning MultiSite-Directed Mutagenesis (Academic) 30 rxn
210514	QC Lightning Multi Site-Directed Mutagenesis (Commercial) 30 rxn
210515	QC Lightning Multi Site-Directed Mutagenesis (Academic) 10 rxn
210516	QC Lightning Multi Site-Directed Mutagenesis (Commercial) 10 rxn
210518	QuikChange Lightning 10 rxn, -20C
210519	QuikChange Lightning 30 rxn
211170	pCMV-Tag 1 Vector
211172	pCMV-Tag2 Vector
211173	pCMV-Tag3 Vector
211174	pCMV-Tag4 Vector
211175	pCMV-Tag5 Vector
211521	pET-3 a-d Vectors
211523	pET-11a-d Vectors
211621	pET-3 a-d Expression system
211623	pET-11 a-d Expression System
212205	pBluescript II SK(+) Phagemid Kit
212206	pBluescript II SK(-) Phagemid Kit
212207	pBluescript II KS(+) Phagemid Kit
212208	pBluescript II KS(-) Phagemid Kit
212209	pBK-CMV Vector

212215	pBC SK(+) Phagemid
212216	pBC SK(-) Phagemid
212217	pBC KS(+) Phagemid
212218	pBC KS(-) Phagemid
212220	pCMV-Script Vector Kit
212224	pCMV-Script XR Predigested Vector
212240	pBluescript II XR Predigested Vector
212250	pBluescript II RI Predigested Vector
213201	pMC1Neo and pMC1Neo Poly A Vectors
214301	pCAL-c Vector
214302	pCAL-n Vector
214303	Calmodulin Affinity Resin
214310	pCAL-n-EK Vector
214311	pCAL-n-FLAG Vector
214501	pDual Expression Vector
214502	pDual Expression Vector
214503	pDual GC Vector 1 ug/ul
216201	pSG5 Expression Vector
217450	LacSwitch II Mammalian Expression System
217451	pESC-HIS Vector
217452	pESC-LEU Vector
217453	pESC-TRP Vector
217454	pESC-URA Vector
217455	pESC Vectors, all four
217460	pERV3 Vector
217461	pEGSH Vector
217467	Ponasterone A
217468	Complete Control Vector Kit
217561	pFB-Neo Vector
217563	pFB Vector
217564	pFB-ERV Vector
217566	pVPack-GP Vector
217567	pVPack-VSV-G Vector
217568	pVPack-Ampho Vector
217569	pVPack-Eco Vector
217570	pVPack-10A1 Vector
219000	PathDetect c-JUN Trans-Reporting System
219001	PathDetect pFR-CAT Trans-Reporter Plsmd
219002	PathDetect pFR-B-Gal Trans-RprtrPlsmd
219004	PathDetect pFR-SEAP Trans-Reportr Plasmid
219005	PathDetect Elk1 Trans-Reporting System
219010	PathDetect CREB Trans-Reporting System
219015	PathDetect CHOP Trans-Reporting System
219020	Luciferase Assay Kit
219026	PathDetect pFA-ATF2 Plasmid
219031	PathDetect pFA-cFos Plasmid
219036	PathDetect pFA-CMV Plasmid

219040	PathDetect pMUT-ELK1 Plasmid
219041	PathDetect pMUT-cJun Plasmid
219042	PathDetect pMUT-CREB Plasmid
219050	PathDetect pFR-Luc Plasmid
219053	PathDetect pFA-Jun Plasmid
219054	PathDetect pFA-CHOP Plasmid
219056	PathDetect pFC-dbd Plasmid
219059	PathDetect pFC-MEKK Plasmid
219062	PathDetect pFA-Elk1 Plasmid
219065	PathDetect pFC-MEK1 Plasmid
219068	PathDetect pFA-CREB Plasmid
219071	PathDetect pFC-PKA Plasmid
219073	PathDetect AP-1 Cis-Reporting System
219074	PathDetect pAP1-Luc Plasmid
219075	PathDetect CRE Cis-Reporting System
219076	PathDetect pCRE-Luc Plasmid
219077	PathDetect NF-kB Cis-Reporting System
219078	PathDetect pNFkB-Luc Plasmid
219079	PathDetect SRE Cis-Reporting System
219080	PathDetect pSRE-Luc Plasmid
219081	PathDetect SRF Cis-Reporting System
219082	PathDetect pSRF-Luc Plasmid
219083	PathDetect p53 Cis-Reporting System
219084	PathDetect pFC-p53 Plasmid
219085	PathDetect p53-Luc Plasmid
219086	PathDetect pFC-MEK3 Plasmid
219087	PathDetect pLuc-MCS Plasmid
219088	PathDetect pNFAT-Luc Plasmid
219089	PathDetect pISRE-Luc Plasmid
219090	PathDetect pCIS-CK Plasmid
219091	PathDetect pGAS-Luc Plasmid
219092	PathDetect ISRE Cis-Reporting System
219093	PathDetect GAS Cis-Reporting System
219094	PathDetect NFAT Cis-Reporting System
219095	PathDetect TARE Cis-Reporting System
230130	BL21-Gold Competent Cells
230132	BL21-Gold (DE3) Competent Cells
230134	BL21-Gold(DE3)pLysS CompCells
230191	ArcticExpress Competent Cells
230192	ArcticExpress (DE3) Competent Cells
230193	ArcticExpress (DE3) RIL Competent Cells
230194	ArcticExpress (DE3) RP Competent Cells
230195	ArcticExpress RIL Competent Cells
230196	ArcticExpress RP Competent Cells
230240	BL21-CodonPlus -RIL Competent Cells
230245	BL21-CodonPlus (DE3)-RIL Comp Cells
230246	Protein Expression Competent Cell Pack

230247	Difficult Cloning Competent Cell Pack
230248	Routine Cloning Competent Cell Pack
230250	BL21-CodonPlus -RP Competent Cells
230255	BL21-CodonPlus (DE3)-RP Comp Cells
230265	BL21-CodonPlus (DE3)-RIL-X Comp Cells
230275	BL21-CodonPlus (DE3)-RP-X Comp cells
230280	BL21-CodonPlus (DE3)-RILP Comp Cells
230325	SoloPack Gold Competent Cells
230350	SoloPack Gold SuperCompetent Cells
235200	Lambda CE6 Induction Kit
240005	pAdEasy-1 Vector
240006	pShuttle Vector
240007	pShuttle-CMV Vector
240008	pShuttle-CMV lacZ Vector
240009	AdEasy Adenoviral Vector System
240010	AdEasy XL Adenoviral System
240027	Vitality pFB-hrGFP Retroviral Vector
240028	pCFB-EGSH Retroviral Vector
240029	pFB-Neo-LacZ Retroviral Vector
240030	pFB-Luc Retroviral Vector
240031	Vitality pIRES-hrGFP-1 Vector
240032	Vitality pIRES-hrGFP-2 Vector
240035	Vitality phrGFP-C Mammalian ExprsnVctr
240036	Vitality phrGFP-N Mammalian ExprsnVctr
240039	PathDetect pTARE-Luc Plasmid
240042	Vitality phr-GFP-Mito Subcellular Vctr
240043	Vitality phrGFP-Nuc Subcellular Vector
240044	Vitality phrGFP-Golgi SubcellularVectr
240049	PathDetect pAP-1-hrGFP Cis-ReprtrPlsmd
240050	PathDetect pCRE-hrGFP Cis-ReprtrPlasmid
240051	PathDetect pNF-kB-hrGFP Cis-RprtrPlsmd
240053	PathDetect pNFAT-hrGFP Cis-ReprtrPlsmd
240059	Vitality phrGFP-1 Vector
240062	Vitality phrGFP Promoterless Vector
240063	Vitality phrGFP-Perox Subcellular Vctr
240071	AAV Helper-Free System
240073	AAV-293 Cells
240074	pAAV-hrGFP Vector
240075	pAAV-IRES-hrGFP Vector
240081	pShuttle-IRES-hrGFP-1 vector
240082	pShuttle-IRES-hrGFP-2 vector
240085	AD-293 Cells
240099	InterPlay TAP Purification Buffer Kit
240101	InterPlay N-terminal Mammalian TAP Vectr
240102	InterPlay C-Terminal Mammalian TAP Vectr
240103	Kit,InterPlay N-Terminal MammalianTAPSy
240104	Kit,InterPlay C-TerminalMammalianTAP Sy

240107	InterPlay Mammalian TAP Purification Kit
240109	AAV-HT1080 Cells
240111	PathDetect c/EBP Cis-Reporting System
240112	PathDetect c/EBP Cis-Reporting Plasmid
240113	PathDetect DR1/RXRE Cis-Reporting System
240114	PathDetect DR1/RXRE Cis-Reporting Plasmid
240115	PathDetect DR3/VDR Cis-Reporting System
240116	PathDetect DR3/VDR Cis-Reporting Plasmid
240119	PathDetect DR5/RARE Cis-Reporting System
240120	PathDetect DR5/RARE Cis-Reporting Plasmid
240129	PathDetect Egr-1 Cis-Reporting System
240130	PathDetect Egr-1 Cis-Reporting Plasmid
240131	PathDetect LILRE Cis-Reporting System
240132	PathDetect LILRE Cis-Reporting Plasmid
240133	PathDetect GRE Cis-Reporting System
240134	PathDetect GRE Cis-Reporting Plasmid
240135	PathDetect DR4/TR Cis-Reporting System
240136	PathDetect DR4/TR Cis-Reporting Plasmid
240141	Vitality full length hrGFP Antibody
240142	Vitality full length hrGFP Antibody
240143	Vitality hrGFP II-1MammalianExprssVctr
240144	Vitality hrGFP II-CMammalianExprssVctr
240145	Vitality hrGFP II-NMammalianExprssVctr
240157	Vitality IRES-hrGFP IIMammlnExprssVctr
240162	VeriFlex N-terminal SBP Vector Set
240163	VariFlex N-terminal SBP System
240164	VariFlex N-terminal SBP-SET Vector Set
240165	VariFlex N-terminal SBP-SET System
240172	VariFlex N-Terminal SET Vector Set
240174	VariFlex C-Terminal SBP Vector
240175	VariFlex C-Terminal SBP System
240176	VariFlex C-Terminal SBP-SET Vector Set
240177	VariFlex C-terminal SBP-SET Expression System
240184	VariFlex C-terminal SET Vector Set
240195	pCMV-3Tag-1 Epitope TaggingMammlnExpVctr
240196	pCMV-3Tag-2 Epitope TaggingMammlnExpVctr
240197	pCMV-3Tag-3 Epitope TaggingMammlnExpVctr
240198	pCMV-3Tag-4 Epitope TaggingMammlnExpVctr
240200	pCMV-3Tag-6 vectors
240202	pCMV-3Tag-7 vectors
240203	pCMV-3Tag-8 vectors
240204	pCMV-3Tag-9 vectors
240205	StrataClone PCR Cloning Kit
240207	StrataClone Blunt PCR Cloning Kit
240213	InterPlay Adenoviral N-Terminal TAP Sys
240214	InterPlay Adenoviral N-terminal TAP Vctr
240215	InterPlay Adenoviral C-Terminal TAP Sys

240216	InterPlay Adenoviral C-Terminal Vectors
240218	StrataClone Ultra Blunt PCR Cloning Kit
240228	StrataClone Mammalian Untagged
240229	StrataClone Mammalian N-Flag
240230	StrataClone Mammalian C-Flag
240231	StrataClone Mammalian N-cMyc
240232	StrataClone Mammalian C-cMyc
240241	Vitality hrGFP monoclonal antibody
240242	Vitality hrGFP monoclonal antibody 100ul
240243	AdEasy Virus Purification Kit
241612	Kt,LambdaEMBL3/Gigapack III Gold Clnin
241711	Kt,Lambda EMBL3/Gigapack II XL Clonin
247201	Lambda DASH II Undigested Vector Kit
247211	Lambda DASH II/BamH I Vector Kit
247212	Lambda DASH II/EcoR I Vector Kit
247613	Kt,Lambda DASH II/BamH I/Gigapack III
247614	Kit,Lambda DASH II/EcoR I/Gigapack III
247713	Kit,LambdaDASH II/BamHI/Gigapack III X
247714	Kit,LambdaDASH II/EcoRI/Gigapack III XL
248211	Lambda FIX II/Xho I Partial Fill-in Vect
248612	Kit,Lambda FIX II/Gigapack III Cloning
248712	Kit, LambdaFIX II/Gigapack III XL Clonin
251301	SuperCos I Vector Kit
5185-5984	Mult Aff Rem Column 4.6 x 50mm Hu-6
5185-5985	Mult Aff Rem Column 4.6 x 100mm Hu-6
5185-5986	Mult Aff Rem Reagent Kit
5185-5987	Buffer A, Equil/Load/Wash, 1L
5185-5988	Buffer B, Elution 1L
5185-5989	HSA Standard
5185-5990	Spin Filters
5185-5991	Spin Concentrators, 5K MWCO, 4ul, 25/pk
5188-2750	Peptide Cleanup C18 Spin Tubes, pk/50
5188-5202	Concentrators, Spin 50K MWCO
5188-5217	Mult Aff Rem Column 4.6 x 50mm Ms-3
5188-5218	Mult Aff Rem Column 4.6 x100mm MS-3
5188-5230	Mult Aff Rem Spin Cart. 0.45mL Hu-6
5188-5231	mRP 4.6 x 50 mm Column
5188-5239	Cleanup C18 Pipette Tips, pk/96
5188-5249	Luer-Lok Adapters
5188-5250	Syringes, Plastic, 5ml Luer-Lok
5188-5251	Microtube, Screwtop, 1.5ml, pk/100
5188-5252	Spin Cartridge Caps and Plugs
5188-5253	Needles, PTFE Luer-Lok
5188-5254	Mult Aff Rem Spin Cartridge Reagent Kit
5188-5289	Mult Aff Rem Spin cart. 0.45mL Ms-3
5188-5332	Mult Aff Rem Column 4.6 x 50mm Hu- HC
5188-5333	Mult Aff Rem Column 4.6 x 100mm Hu-6HC

5188-5334	Mult Aff Rem spin Cart, Hu-Albumin
5188-5336	10 x 100 High Capacity Hu6 MARS
5188-5341	Mult Aff Rem Spin Cart 0.45mL Hu-6HC
5188-6408	Multiple Affinity Spin Cartridge Hu-PL7
5188-6409	Multiple Affinity 4.6 x 50 Hu-PL7
5188-6410	Multiple Affinity 4.6 x 100 Hu-PL7
5188-6411	Multiple Affinity 10 x 100mm, Hu-7
5188-6510	mRP 0.5 x 100mm Column
5188-6557	Mult Aff Rem Column,4.6 x 50mm,Hu-14
5188-6558	Mult Aff Rem Column,4.6 x 100mm,Hu-14
5188-6559	Mult Aff Rem Column,10 x 100mm,Hu-14
5188-6560	Mult Aff Rem Spin Cart Hu-14,0.45mL
5188-6562	Mult Aff Rem Column, 4.6 x 50mm, Hu-Albumin
5188-8825	Mult. Aff. Rem. Spin Cartr. HSA/IgG, 0.4
5188-8826	Mult. Aff. Rem. Col. HSA/IgG, 4.6 x 50 m
5190-7995	MARS Col 2um Replacement Frit
G5900A	QuikChange HT 150nt, 10 sites
G5900B	QuikChange HT 150nt, 20 sites
G5901A	QuikChange HT 200nt, 10 sites
G5901B	QuikChange HT 200nt, 20 sites
G5902A	QuikChange HT 150nt-academic, 10sites
G5902B	QuikChange HT 150nt-academic, 20sites
G5903A	QuikChange HT 200nt-academic, 10sites
G5903B	QuikChange HT 200nt-academic, 20sites