

# **PromegaExpress™ Product List**

## Table of Contents

<u>Amplification</u> .....	2
<u>Biochemicals and Florescent Dyes</u> .....	2
<u>Cell Signaling and Drug Screening Assays</u> .....	3
<u>DNA and RNA Purification</u> .....	4
<u>Genetic Identity</u> .....	5
<u>Genetic Reporters and Transfection Systems</u> .....	6
<u>Imaging and Immunological Detection</u> .....	7
<u>Industrial and Environmental Monitoring</u> .....	8
<u>in vitro Transcription /Translation Regulation</u> .....	8
<u>Molecular Weight Markers</u> .....	8
<u>Protein Expression, Purification and Analysis</u> .....	9
<u>Restriction/Modifying Enzymes and RNase Inhibitors</u> .....	10
<u>Sequencing Mutagenesis and Labeling</u> .....	11
<u>Vectors and Cloning Tools</u> .....	11

<b>Amplification</b>	
Access RT-PCR System	Random Primers
AccessQuick RT-PCR System	PCR Master Mix
dATP	PCR Nucleotide Mix
dCTP	pGEM-T Easy Vector System I
dGTP	pGEM-T Vector System I
dTTP	Beta-Actin Primer Pair
Set of dATP, dCTP, dGTP, dTTP	CNTF Primer Pair
dUTP	NT-3 Primer Pair
dNTP Mix	p75 Primer Pair
GoTaq Colorless Master Mix	TrkB Primer Pair
GoTaq DNA Polymerase	Promega Barrier Tips
GoTaq Flexi DNA Polymerase	Plexor One-Step qRT-PCR System
GoTaq Green Master Mix	Plexor qPCR System
GoTaq PCR Core System I	Plexor Two-Step qRT-PCR System
GoTaq PCR Core System II	Reverse Transcription 10X Buffer
5X Colorless GoTaq Flexi Reaction Buffer	Reverse Transcription System
5X Colorless GoTaq Reaction Buffer	TaqBead Hot Start Polymerase, 1.25u/bead
5X Green GoTaq Flexi Reaction Buffer	Tfl DNA Polymerase
5X Green GoTaq Reaction Buffer	Tli DNA Polymerase
Magnesium Chloride Solution	Tth DNA Polymerase
ImProm-II Reverse Transcriptase	Oligo(dT)15 Primer

<b>Biochemicals and Fluorescent Dyes</b>	
5M Sodium Chloride	DABCYL Acid
Acrylamide	DABCYL, SE
Acrylamide/N,N'-Methylenebisacrylamide Solution	DABSYL chloride
5(6)-CFDA, SE	Rhodamine 110
5(6)-FAM, SE	SDS Solution (10% w/v)
5-FAM, SE	Sodium Chloride
5-FITC	Sodium Dodecyl Sulphate
5-ROX, SE	SSC Buffer, 20X
5-TAMRA, SE	Dil
6-FAM, SE	JC-1
6-FITC	TAE Buffer
6-ROX, SE	TBE Buffer, 10X
6-TAMRA, SE	TCEP
Ammonium Persulfate	TE Buffer, 1X
Ammonium Sulfate	5-TMRIA (Tetramethylrhodamine-5-iodoacetamide)
NBD chloride	Fluorescein-5-maleimide
NBD fluoride	SBF-Cl
Biotin, SE (Biotin, succinimidyl ester)	Tris Base
Bisacrylamide	Tris-HCl
Boric Acid	Triton X-100

<b>Biochemicals and Fluorescent Dyes</b>	
Bovine Serum Albumin, Acetylated	Tween 20
Calcein, Acetoxymethyl Ester	Urea, 99% Pure
CTC---MaxxPure Grade	1,5-IAEDANS
Dextran Sulphate	7-Hydroxycoumarin-3-carboxylic acid
DTT	7-Hydroxycoumarin-3-carboxylic acid, succinimidyl ester
EDTA, 0.5M (pH 8.0)	7-Methoxycoumarin-3-carboxylic acid
EDTA, Disodium Salt	7-Methoxycoumarin-3-carboxylic acid, succinimidyl ester
Ethidium Bromide Solution	EDANS
Formamide	Fluorescamine
Glycerol	5(6)-ROX (5(6)-Carboxy-X-rhodamine)
Glycine	5(6)-TAMRA (5(6)-Carboxytetramethylrhodamine)
Guanidine Thiocyanate	5-Amino TMR (5-Aminotetramethylrhodamine)
Guanidine-HCl	5-FAM (5-Carboxyfluorescein)
HEPES (free acid)	5-ROX, TEA Salt
2'-Fluoro-dATP, Li Salt	5-TAMRA
2'-Fluoro-dCTP, Li Salt	6-FAM
2'-Fluoro-dGTP, Li Salt	6-ROX
2'-Fluoro-dUTP, Li Salt	6-TAMRA
2'-O-Methyl-rATP, Li Salt	Fluorescein Cadaverine
2'-O-Methyl-rCTP, Li Salt	5-aminoallyl-dUTP, Li Salt
2'-O-Methyl-rGTP, Li Salt	MOPS
2'-O-Methyl-rUTP, Li Salt	Nuclease-Free Water
PEG 8000 Powder	

<b>Cell Signaling and Drug Screening Assays</b>	
4alpha-PMA	ProFluor Src-Family Kinase Assay
Apo-ONE Homogeneous Caspase-3/7 Assay	ProFluor Tyrosine Phosphatase Assay
BacTiter-Glo Microbial Cell Viability Assay	Proteasome-Glo 3-Substrate System
Calpain-Glo Protease Assay	Proteasome-Glo Caspase-Like Assay
cAMP-Dependent Protein Kinase Peptide Inhibitor	Proteasome-Glo Cell-Based Assay
cAMP-Glo Assay	Proteasome-Glo Chymotrypsin-Like Assay
Casein Kinase I	DPPIV-Glo Protease Assay
Casein Kinase I Peptide Substrate	rhEGF
Casein Kinase II Peptide Substrate	Griess Reagent System
CaspACE Assay System	mNGF, 2.5S
CaspACE FITC-VAD-FMK In Situ Marker	rhBDNF
Caspase Inhibitor Ac-DEVD-CHO	rhFGF, Basic
Caspase Inhibitor Z-VAD-FMK, 20mM	rhGDNF
Caspase-Glo 3/7 Assay	rhIGF-I
Caspase-Glo 8 Assay	rhIL-4
Caspase-Glo 9 Assay	rhNT-3
cdc2 Protein Kinase Peptide Substrate	rhTNF-alpha
CellTiter 96 AQueous MTS Reagent Powder	InCELLect AKAP St-Ht31 Inhibitor Peptide

<b>Cell Signaling and Drug Screening Assays</b>	
CellTiter 96 AQueous Cell Proliferation Assay	InCELLect St-Ht31P Control Peptide
CellTiter 96 AQueous One Solution Cell Proliferation Assay	Kemptide (PKA) Peptide Substrate
CellTiter 96 Non-Radioactive Cell Proliferation Assay	Kinase-Glo Luminescent Kinase Assay
CellTiter-Blue Cell Viability Assay	Kinase-Glo Plus Luminescent Kinase Assay
CellTiter-Glo Luminescent Cell Viability Assay	LY 294002
cAMP, 1mM	MAO-Glo Assay
cGMP, 1mM	MAO-Glo Assay
cGMP-Dependent Protein Kinase Peptide Substrate	MEK Inhibitor U0126
NADPH Regeneration System	MultiTox-Fluor Multiplex Cytotoxicity Assay
P450-Glo CYP1A1 Assay	Myristoylated Protein Kinase C Peptide Inhibitor
P450-Glo CYP1A2 Assay	Neurogranin (28-43) (PKC) Peptide Substrate
P450-Glo CYP1B1 Assay	Serine/Threonine Phosphatase Assay System
P450-Glo CYP2C19 Assay	Tyrosine Phosphatase Assay System
P450-Glo CYP2C8 Assay	Olomoucine (cdc2 Protein Kinase Inhibitor)
P450-Glo CYP2C9 Assay	PD 98059
P450-Glo CYP2D6 Assay	PDE-Glo Phosphodiesterase Assay
P450-Glo CYP3A4 Assay	PepTag cAMP-Dependent Protein Kinase Assay
P450-Glo CYP3A7 Assay	Proteasome-Glo Trypsin-Like Assay
CytoTox 96 Non-Radioactive Cytotoxicity Assay	psiCHECK-1 Vector
CytoTox-Fluor Cytotoxicity Assay	psiCHECK-2 Vector
CytoTox-ONE Homogeneous Membrane Integrity Assay	SAM2 96 Biotin Capture Plate
DeadEnd Colorimetric TUNEL System	SAM2 Biotin Capture Membrane
DNA-Dependent Protein Kinase Peptide Substrate	Kinase-Glo Luminescent Kinase Assay
PepTag Non-Radioactive PKC Assay	SB 203580
Pgp-Glo Assay System	SignaTECT Protein Kinase Assay System
PMA	SignaTECT Protein Tyrosine Kinase Assay System
PPase-2A	rhLung Beta Tryptase
ProFluor PKA Assay	rhSkin Beta Tryptase
ProFluor Ser/Thr PPase Assay	

<b>DNA and RNA Purification</b>	
Maxwell 16 DNA Purification Kits	Wizard DNA Clean-Up System
Maxwell 16 Cell RNA Purification Kits	DNA Rehydration Solution
Maxwell 16 LEV Total RNA Purification Kit	Nuclei Lysis Solution
Maxwell 16 Polyhistidine Protein Purification Kit	Protein Precipitation Solution
CelluACE XG System	Proteinase K
Direct Purification Buffer	RNase A Solution
DNA Quantitation System	Wizard Genomic DNA Purification Kit
MagneSil KF Genomic System	Cell Lysis Solution

<b>DNA and RNA Purification</b>	
384-Well Plate, Conical	Cell Resuspension Solution
384-Well Plate, Flat	Elution Buffer
Alcohol Wash, Blood	Neutralization Solution
Anti-Foam Reagent	Wizard MagneSil Plasmid Purification System
Collection Plates (4-pack)	Wizard MagneSil Sequencing Reaction Clean-Up System
MagneSil Blood Genomic, Max Yield System	4/40 Wash Solution
Salt Wash, Blood	Endotoxin Removal Resin
eLysis Buffer, Large Volume System	Wizard MagneSil Tfx System
Lysis Buffer	Wash Buffer, Plant
MagneSil ONE, Fixed Yield Blood Genomic System	Wizard Magnetic 96 DNA Plant System
MagneSil BLUE	Precipitation Solution, Food
MagneSil GREEN	Wizard Magnetic DNA Purification System for Food
MagneSil RED	Column Wash Solution (CWB)
MagneSil Total RNA mini-Isolation System	Alkaline Protease Solution
Biotinylated Oligo(dT) Probe (50pmol/ul)	Wizard Plus SV Minipreps DNA Purification System
PolyATtract mRNA Isolation System I	Wizard SV 96 Binding Plates
PolyATtract mRNA Isolation System IV	Wizard SV 96 Cell Lysis Solution
PolyATtract GTC Extraction Buffer	Wizard SV 96 Cell Resuspension Solution
PolyATtract System 1000	Wizard SV 96 Lysate Clearing Plates
PureYield Plasmid Midiprep System	Wizard SV 96 Neutralization Solution
PureYield RNA Midiprep System	Wizard SV 96 Plasmid DNA Purification System
Red Blood Cell Lysis Solution (CLB)	Wizard SV 96 Wash Solution
RNA Lysis Buffer (RLA)	Wizard SV 9600 Plasmid DNA Purification System
RNA Wash Solution (RWA)	Wizard SV Wash Solution
ReadyAmp Genomic DNA Purification System	Wizard SV 96 Genomic DNA Purification System
RNAgents Denaturing Solution	Wizard SV Lysis Buffer
RNAgents Total RNA Isolation System	Membrane Binding Solution
Streptavidin	Wizard SV 96 PCR Clean-Up System
Streptavidin MagneSphere Paramagnetic Particles	Wizard SV Gel and PCR Clean-Up System
SV 96 Total RNA Isolation System	Wizard SV Genomic DNA Purification System
SV Total RNA Isolation System	PureYield Plasmid Maxiprep System

<b>Genetic Identity</b>	
AluQuant Human DNA Quantitation System	1.1ml, Square-Well, V-Bottom Deep Well Plate
Disposable Polypropylene Test Tubes (8 x 50mm)	1.2ml, Round-Bottom Deep Well Plate
ENLITEN rLuciferase/Luciferin Reagent	10ml, 24-Well Deep Well Plate
ART Pipet Tips	2.2ml, Square-Well Deep Well Plate
Agarose	Deep Well Heat Transfer Block
Blue Dextran Loading Solution	Pyramid-Bottom Reservoir
Bromophenol Blue Loading Solution	Shaker Integration Plate
Gel Tracking Dye	Internal Lane Standard 600
Gold ST*R 10X Buffer	K562 DNA High Molecular Weight

<b>Genetic Identity</b>	
Mineral Oil	Plexor HY System
STR 10X Buffer	Water, Amplification Grade
STR 2X Loading Solution	PowerPlex 1.1 and 2.1 Systems
Differex System	D16S539 Add-In for PowerPlex 1.1
DNA IQ Reference Sample Kit for Maxwell 16	PowerPlex 1.2 System
DNA IQ Spin Baskets	PowerPlex 16 Monoplex Systems
DNA IQ System	9947A DNA
Tissue and Hair Extraction Kit (for use with DNA IQ)	PowerPlex 16 BIO System
Fluorescent Ladder (CXR), 60-400 Bases	PowerPlex 16 System
FFFL Allelic Ladder Mix (Fluorescein)	PowerPlex Matrix Standards
GammaSTR Allelic Ladder Mix (Fluorescein)	PowerPlex ES System
GammaSTR Multiplex (Fluorescein)	9948 Male DNA
Amelogenin	PowerPlex Y System
CSF1PO	HPRTB
CTT Allelic Ladder Mix	LPL
FESFPS	TH01
FFv Allelic Ladder Mix	TPOX
GenePrint SilverSTR III System	vWA

<b>Genetic Reporters and Transfection Systems</b>	
Antibiotic G-418 Sulfate Solution	pGL4.15[luc2P/Hygro] Vector
Beetle Luciferin, Potassium Salt	pGL4.16[luc2CP/Hygro] Vector
Beta-Galactosidase Enzyme Assay System	pGL4.17[luc2/Neo] Vector
Reporter Lysis 5X Buffer	pGL4.18[luc2P/Neo] Vector
Beta-Glo Assay System	pGL4.19[luc2CP/Neo] Vector
Bright-Glo Luciferase Assay System	pGL4.20[luc2/Puro] Vector
CAT Enzyme Assay System	pGL4.21[luc2P/Puro] Vector
Chloramphenicol Acetyltransferase	pGL4.22[luc2CP/Puro] Vector
Chroma-Glo Luciferase Assay System	pGL4.23[luc2/minP] Vector
pCBG68-Basic Vector	pGL4.24[luc2P/minP] Vector
pCBG68-Control Vector	pGL4.25[luc2CP/minP] Vector
pCBG99-Basic Vector	pGL4.26[luc2/minP/Hygro] Vector
pCBG99-Control Vector	pGL4.27[luc2P/minP/Hygro] Vector
pCBR-Basic Vector	pGL4.28[luc2CP/minP/Hygro] Vector
pCBR-Control Vector	pGL4.29[luc2P/CRE/Hygro] Vector
CodeBreaker siRNA Transfection Reagent	pGL4.30[luc2P/NFAT-RE/Hygro] Vector
2-Methylcoelenterazine	pGL4.31[luc2P/Gal4UAS/Hygro] Vector
Black Lite Coelenterazine	pGL4.70[hRluc] Vector
Coelenteramine	pGL4.71[hRlucP] Vector
Dual-Glo Luciferase Assay System	pGL4.72[hRlucCP] Vector
Dual-Luciferase Reporter 1000 Assay System	pGL4.73[hRluc/SV40] Vector
Passive Lysis 5X Buffer	pGL4.74[hRluc/TK] Vector
EnduRen Live Cell Substrate	pGL4.75[hRluc/CMV] Vector
Glo Lysis Buffer, 1X	pGL4.76[hRluc/Hygro] Vector
GloResponse CRE-luc2P HEK293 Cell Line	pGL4.77[hRlucP/Hygro] Vector
GloResponse NFAT-RE-luc2P HEK293 Cell Line	pGL4.78[hRlucCP/Hygro] Vector
Luciferase 1000 Assay System	pGL4.79[hRluc/Neo] Vector

<b>Genetic Reporters and Transfection Systems</b>	
Luciferase Assay System	pGL4.80[hRlucP/Neo] Vector
Luciferase Cell Culture Lysis 5X Reagent	pGL4.81[hRlucCP/Neo] Vector
Luciferin-EF	pGL4.82[hRluc/Puro] Vector
n-Butyryl CoA	pGL4.83[hRlucP/Puro] Vector
ONE-Glo Luciferase Assay System	pGL4.84[hRlucCP/Puro] Vector
pGL4.10[luc2] Vector	ProFection Mammalian Transfection System
pGL4.11[luc2P] Vector	pSP-luc+NF Fusion Vector
pGL4.12[luc2CP] Vector	pSV-beta-Galactosidase Control Vector
pGL4.13[luc2/SV40] Vector	Renilla Luciferase Assay System
pGL4.14[luc2/Hygro] Vector	GLprimer1 (clockwise)
phRL-TK Vector	GLprimer2 (counterclockwise)
phRL-TK(Int-) Vector	RVprimer3 (clockwise)
Tfx Reagents Transfection Trio	RVprimer4 (counterclockwise)
Tfx-20 Reagent	Steady-Glo Luciferase Assay System
Tfx-50 Reagent	phRG-B Vector
TransFast Transfection Reagent	phRG-TK Vector
Transfectam Reagent for Eukaryotic Cells	Disposable Cuvettes, 1,000/pkg
	Range Extender

<b>Imaging and Immunological Detection</b>	
Anti-Human IgG (H+L)	Anti-pS473 Akt pAb
Anti-Mouse IgG (H+L)	Anti-pT183 MAPK pAb, Rabbit
Anti-Rabbit IgG	Anti-Rat CNTF pAb
Anti-Rat IgG (H+L), AP Conjugate	Anti-TGFbeta1 pAb
BCIP/NBT Color Development Substrate	Anti-TrkB In pAb
Donkey Anti-Goat IgG	Anti-VACHT pAb
Anti-ACTIVE CaM KII pAb, Rabbit, (pT286)	AttoPhos AP Fluorescent Substrate System
Anti-ACTIVE Caspase-3 pAb	AttoPhos Buffer
Anti-ACTIVE JNK pAb, Rabbit, (pTPpY)	AttoPhos Substrate
Anti-ACTIVE MAPK pAb, Rabbit, (pTEpY)	BDNF Emax ImmunoAssay System
Anti-ACTIVE p38 pAb, Rabbit, (pTGpY)	Block & Sample 5X Buffer
Anti-beta-Galactosidase, Purified Monoclonal Antibody	Blot-Qualified BSA
Anti-BetaIII Tubulin mAb	E. coli Extract for Background Reduction
Anti-Chicken IgY	Tween
Anti-Cytochrome c mAb	Chicken IgY, Control Immunoglobulin
Anti-ERK 1/2 pAb, Rabbit	Donkey Anti-Rabbit IgG
Anti-GFAP pAb	EGGstract IgY Purification System
Anti-Human BDNF pAb	GDNF Emax ImmunoAssay System
Anti-Human GDNF pAb	Monster Green Fluorescent Protein pHMGFP Vector
Anti-Human NT-3 pAb	NGF Emax ImmunoAssay System
Anti-Human p75 pAb	NT-3 Emax ImmunoAssay System
Anti-Human Tryptase mAb Biotin	ProtoBlot II AP System with Stabilized Substrate
Anti-Luciferase pAb	Streptavidin Alkaline Phosphatase
Anti-NGF mAb	TGF beta 1 Emax ImmunoAssay System
Anti-Pan Trk pAb	TGF beta 2 Emax ImmunoAssay System

<b>Imaging and Immunological Detection</b>	
Anti-PARP p85 Fragment pAb	TGF beta Sample 10X Buffer
HaloTag Oregon Green Ligand	TMB One Solution
HaloTag PEG-Biotin Ligand	HaloTag Alexa Fluor 488 Ligand
HaloTag pHT2 Vector	HaloTag Amine (O2) Ligand
HaloTag Succinimidyl Ester (O2) Ligand	HaloTag Amine (O4) Ligand
HaloTag Succinimidyl Ester (O4) Ligand	HaloTag Biotin Ligand
HaloTag Thiol (O4) Ligand	HaloTag Coumarin Ligand
HaloTag TMR Ligand	HaloTag diAcFAM Ligand
Anti-HaloTag pAb	HaloTag Iodoacetamide (O2) Ligand
Western Blue Stabilized Substrate for Alkaline Phosphatase	HaloTag Iodoacetamide (O4) Ligand
	TMB Stabilized Substrate for Horseradish Peroxidase

<b>Industrial and Environmental Monitoring</b>	
ENLITEN ATP Assay System	ISOQUANT Isoaspartate Detection Kit
ENLITEN Total ATP Rapid Biocontamination Detection Kit	

<b><i>in vitro</i> Transcription/Transcription Regulation</b>	
rhAP1 (c-Jun)	RiboMAX Large Scale RNA Production System
ChipShot Direct Labeling and Clean-Up System	Riboprobe Combination System
ChipShot Indirect Labeling and Clean-Up System	rATP, rCTP, rGTP, rUTP, each
Gel Shift Binding 5X Buffer	Riboprobe System Buffers
rATP	RQ1 RNase-Free DNase
rCTP	Transcription Optimized 5X Buffer
rGTP	Riboprobe System
rUTP	T7 RiboMAX Express Large Scale RNA Production System
pGEM Express Positive Control Template	AP1 Consensus Oligonucleotide
Primer Extension System---AMV Reverse Transcriptase	AP2 Consensus Oligonucleotide
Ribo m7G Cap Analog	CREB Consensus Oligonucleotide
SP1 Consensus Oligonucleotide	NF-kappaB Consensus Oligonucleotide
TFIID Consensus Oligonucleotide	OCT1 Consensus Oligonucleotide
GeneClip U1 Hairpin Cloning Systems	siSTRIKE U6 Hairpin Cloning System
	T7 RiboMAX Express RNAi System

<b>Molecular Weight Markers</b>	
100bp DNA Ladder	Ethidium Bromide Solution
100bp DNA Step Ladder	Lambda DNA/EcoRI + HindIII Markers
10bp DNA Step Ladder	Lambda DNA/EcoRI Markers
1kb DNA Ladder	Lambda DNA/HindIII Markers
1kb DNA Step Ladder	pGEM DNA Markers
200bp DNA Step Ladder	phiX174 DNA/HaeIII Markers
25bp DNA Step Ladder	phiX174 DNA/HinfI Markers
50bp DNA Step Ladder	PCR Markers
Analytical Marker DNA Wide Range	200bp DNA Step Ladder

<b>Molecular Weight Markers</b>	
BenchTop 100bp DNA Ladder	25bp DNA Step Ladder
BenchTop 1kb DNA Ladder	50bp DNA Step Ladder
BenchTop PCR Markers	Human Genomic DNA
BenchTop pGEM DNA Markers	Mouse Genomic DNA
BenchTop PhiX174 DNA/HaeIII Markers	phiX174 DNA/Hinfl Dephosphorylated Markers
Blue/Orange Loading Dye, 6X	Supercoiled DNA Ladder, 2-10kb
Broad Range Protein Molecular Weight Markers	

<b>Protein Expression, Purification and Analysis</b>	
CheckMate Mammalian Two-Hybrid System	HaloCHIP System
CheckMate Negative Control Vectors	HisLink 96 Protein Purification System
CheckMate Positive Control Vectors	HisLink Protein Purification Resin
CheckMate/Flexi Vector Mammalian Two-Hybrid System	HisLink Spin Protein Purification System
pFN10A (ACT) Flexi Vector	MagneGST Glutathione Particles
pFN11A (BIND) Flexi Vector	MagneGST Protein Purification System
PinPoint Vector Sequencing Primer	MagZ Protein Purification System
PinPoint Xa Protein Purification System	pAdVantage Vector
Factor Xa Protease	pCI Mammalian Expression Vector
FastBreak Cell Lysis Reagent, 10X	pCI-neo Mammalian Expression Vector
10X Flexi Enzyme Blend (Sgfl and Pmel)	pCMVTNT Vector
Carboxy Flexi System, Transfer	pGEM Beta-Gal Control DNA
Flexi System, Entry/Transfer	PinPoint Xa-1 Vector
HaloLink Magnetic Beads	ProTEV Protease
HaloLink Resin	pSI Mammalian Expression Vector
pF1A T7 Flexi Vector	pTNT Vector
pF1K T7 Flexi Vector	Coumermycin A1
pF3A WG (BYDV) Flexi Vector	Novobiocin Sodium Salt
pF3K WG (BYDV) Flexi Vector	pFC8A (HaloTag) CMV Flexi Vector
pF4A CMV Flexi Vector	pFC8K (HaloTag) CMV Flexi Vector
pF4K CMV Flexi Vector	pF12A RM Flexi Vector
pF5A CMV-neo Flexi Vector	pF12K RM Flexi Vector
pF5K CMV-neo Flexi Vector	pReg neo Vector
pF9A CMV hRluc-neo Flexi Vector	Regulated Mammalian Expression System
pFC7A (HQ) Flexi Vector	Sequencing Grade Modified Trypsin
pFC7K (HQ) Flexi Vector	L-Rhamnose Monohydrate
pFN2A (GST) Flexi Vector	SoftLink Soft Release Avidin Resin
pFN2K (GST) Flexi Vector	TetraLink Tetrameric Avidin Resin
pFN6A (HQ) Flexi Vector	Magnesium Acetate
pFN6K (HQ) Flexi Vector	Potassium Chloride
Gel Drying Film	Trypsin Gold, Mass Spectrometry Grade
Gel Drying Kit	MagneHis Ni-Particles
	MagneHis Protein Purification System

Restriction/Modifying Enzymes and RNase Inhibitors		
4-CORE Buffer Pack	EcoICRI	RNasin Ribonuclease Inhibitor
AatII	EcoRI	RsaI
Acc65I	EcoRV	S1 Nuclease
AccB7I	Exonuclease III	SacI
AccI	FokI	SacII
AccIII	HaeII	SalI
AgarACE Enzyme	HaeIII	Sau3AI
AgeI	HhaI	Scal
Alkaline Phosphatase,	HincII	SfiI
CIAP Buffer Pack	HindIII	SgfI
AluI	HinfI	Shrimp Alkaline Phosphatase
Alw26I	HpaI	ssDNA Binding Protein
Alw44I	HpaII	SinI
AMV Reverse Transcriptase	Hsp92I	Smal
ApaI	Hsp92II	SnaBI
AvaI	I-PpoI	SP6 RNA Polymerase
AvaII	KpnI	SpeI
BalI	Rapid Ligation Buffer	SphI
BamHI	LigaFast	SspI
BanI	MboI	StuI
BanII	MboII	StyI
BbuI	MluI	T3 RNA Polymerase
BclI	M-MLV Reverse Transcriptase	T4 DNA Ligase
BglI	MspA1I	T4 DNA Ligase Buffer Pack
BglII	MspI	T4 DNA Polymerase
BsaMI	MULTI-CORE Buffer Pack	T4 PNK Buffer Pack
Bsp1286I	Mung Bean Nuclease	T4 Polynucleotide Kinase
BsrSI	NaeI	T4 RNA Ligase
BssHII	NarI	T7 RNA Polymerase
Bst98I	NciI	TaqI
BstEII	NcoI	rTerminal Deoxynucleotidyl Transferase
BstOI	NdeI	Terminal Transferase Buffer Pack
BstXI	NdeII	Thioredoxin
BstZI	NgoMIV	Topoisomerase I
Bsu36I	NheI	Tru9I
CfoI	NotI	TSAP
Clal	NruI	Tth1111
Csp45I	NsiI	Turbo NaeI
CspI	PstI	Turbo NarI
Ddel	PvuI	VspI
DNA Polymerase I	PvuII	XbaI
Klenow Fragment	Ribonuclease H	XhoI
Dpnl	RNase ONE Ribonuclease	XhoII
Dral	RNasin Plus RNase Inhibitor	XmaI
EclHKL	rRNasin Ribonuclease Inhibitor	XmnI

<b>Sequencing, Mutagenesis and Labeling</b>	
Ampicillin Repair Oligonucleotide	Prime-a-Gene Labeling System
pALTER-MAX Vector	pTarget Sequencing Primer
Automatic Processor Compatible (APC) Film	pUC/M13 Primer, Forward
DNA 5' End-Labeling System	pUC/M13 Primer, Reverse
Erase-a-Base System (minus vectors & bacterial strain)	SP6 Promoter Primer
fmol DNA Cycle Sequencing System	T3 Promoter Primer
GeneEditor Antibiotic Selection Mix	T7 EEV Promoter Primer
Bottom Strand Selection Oligonucleotide	T7 Promoter Primer
Labeling 5X Buffer	SILVER SEQUENCE DNA Sequencing System
	Top Strand Selection Oligonucleotide

<b>Vectors and Cloning Tools</b>	
Herring Sperm DNA	pGEM-7Zf(-) Vector
IPTG, Dioxane-Free	pGEM-7Zf(+) Vector
Lambda DNA	pGEM-9Zf(-) Vector
pBR322 Vector	phiX174, RF DNA
pGEM-11Zf(-) Vector	pSP64 Poly(A) Vector
pGEM-11Zf(+) Vector	pSP72 Vector
pGEM-13Zf(+) Vector	pSP73 Vector
pGEM-3Z Vector	R408 Helper Phage
pGEM-3Zf(-) Vector	R408 Helper Phage DNA
pGEM-3Zf(+) Vector	Sephacryl S-400
pGEM-4Z Vector	Spin Columns
pGEM-5Zf(-) Vector	Universal RiboClone cDNA Synthesis System
pGEM-5Zf(+) Vector	Unmethylated Lambda DNA
EcoRI Adaptors	X-Gal