

프로메가 2018 여름이벤트

SUMMER BIG SALE



행사기간 | 2018년 8월 06일 ~ 9월 14일



*본 행사지에 표기된 금액은 부가세 별도 금액입니다.
*단가 계약 및 입찰은 행사에서 제외됩니다.

타Br을 아직 사용하신다면?
일회용 Gel extractor를 사용해 보세요!
내 몸은 소중한니까~

15ul minimum elution이
가능합니다!!

프로메가 Steady seller "T-vector"!!

A260/A230 ratio가 1.8-2.0으로
안정적이고, ethanol을 사용하지 않아,
NGS 목적에 적합합니다.

프로메가 FFPE kit는 xylene을
사용하지 않고, mineral oil을 사용하여
파라핀을 제거합니다.

Cat. No.	제품명	소비자가	이벤트가	할인율
Clean-up				
A2121	x-tracta™ Gel Extractor, 25/pack	47,000	24,000	
A2122	x-tracta™ Gel Extractor, 100/pack	146,000	73,000	
A9281	Wizard® SV Gel and PCR Clean-up System, 50 preps	116,000	58,000	50%
A9282	Wizard® SV Gel and PCR Clean-up System, 250 preps	476,000	238,000	
A9285	Wizard® SV Gel and PCR Clean-up System, 1,000 preps	1,521,000	761,000	그레이트!
Cloning				
A1360	pGEM®-T Easy Vector System I	252,000	189,000	
A1380	pGEM®-T Easy Vector System II	422,000	317,000	
A3600	pGEM®-T Vector System I	207,000	156,000	
A3610	pGEM®-T Vector System II	380,000	285,000	25%
M1801	T4 DNA Ligase, 100u	73,000	55,000	
M1804	T4 DNA Ligase, 500u	291,000	219,000	
M8221	LigaFast™ Rapid DNA Ligation System, 30 rxn	126,000	95,000	
M8225	LigaFast™ Rapid DNA Ligation System, 150 rxn	412,000	309,000	
V3941	X-Gal, 100mg	133,000	67,000	
V3951	IPTG, Dioxane-Free, 5g	324,000	162,000	50%
V3955	IPTG, Dioxane-Free, 1g	87,000	44,000	그레이트!
gDNA				
A2051	ReliaPrep™ gDNA Tissue Miniprep System, 100 Preps	388,000	291,000	
A2052	ReliaPrep™ gDNA Tissue Miniprep System, 250 Preps	886,000	665,000	
A5081	ReliaPrep™ Blood gDNA Miniprep System, 100 Preps	317,000	238,000	25%
A5082	ReliaPrep™ Blood gDNA Miniprep System, 250 Preps	725,000	544,000	
A2352	ReliaPrep™ FFPE gDNA Miniprep System, 100 Preps	465,000	349,000	
A1120	Wizard® Genomic DNA Purification Kit, 100 Isolations	240,000	192,000	
A1125	Wizard® Genomic DNA Purification Kit, 500 Isolations	600,000	480,000	
A1620	Wizard® Genomic DNA Purification Kit, 100 isolations ×10ml	1,481,000	1,185,000	20%
A2360	Wizard® SV Genomic DNA Purification System, 50 preps	176,000	141,000	
A2361	Wizard® SV Genomic DNA Purification System, 250 preps	728,000	583,000	
Plasmid purification				
A1330	Wizard® Plus SV Minipreps DNA Purification System, 50 preps	103,000	52,000	50%
A1460	Wizard® Plus SV Minipreps DNA Purification System, 250 preps	417,000	209,000	
A1222	PureYield™ Plasmid Miniprep System, 250 preps	631,000	505,000	
A1223	PureYield™ Plasmid Miniprep System, 100 preps	282,000	226,000	
A2492	PureYield™ Plasmid Midiprep System, 25 preps	294,000	236,000	
A2495	PureYield™ Plasmid Midiprep System, 100 Preps	1,079,000	864,000	
A2392	PureYield™ Plasmid Maxiprep System, 10 preps	253,000	203,000	20%
A2393	PureYield™ Plasmid Maxiprep System, 25 preps	598,000	479,000	
A6742	PureYield™ Plasmid Midiprep Start-Up Kit, 50 prep	1,506,000	1,205,000	
A6732	PureYield™ Plasmid Maxiprep Start-Up Kit, 100 preps	1,621,000	1,297,000	
Marker				
G3191	RNA Markers, 50ul	197,000	99,000	
G4471	10bp DNA Step Ladder, 32.5ug (50 lanes)	192,000	96,000	
G4511	25bp DNA Step Ladder, 100ug (55 lanes)	192,000	96,000	
G4521	50bp DNA Step Ladder, 340ug/ml 90ug (52 lanes)	225,000	113,000	
G5711	DNA Ladder, 1kb, 500ul (100 lanes)	203,000	102,000	
G6951	100bp DNA Step Ladder, 100ug (100 lanes)	280,000	140,000	50%
G7521	BenchTop pGEM® DNA Markers, 250ul	238,000	119,000	
G7541	BenchTop 1kb DNA Ladder, 600ul	224,000	112,000	
G8291	BenchTop 100bp DNA Ladder, 300ul (50 lanes)	198,000	99,000	
H5041	Ethidium Bromide Solution, 10ml	70,000	35,000	그레이트!
dNTP				
C1141	PCR Nucleotide Mix, 10mM, 200ul	111,000	78,000	
C1145	PCR Nucleotide Mix, 10mM, 1,000ul	428,000	300,000	
U1240	4 dNTP's, 40uMoles each	473,000	332,000	30%
U1511	dNTP Mix, 10mM, 200ul	38,000	27,000	
U1515	dNTP Mix, 10mM, 1,000ul	90,000	63,000	

Mini
Midi
Maxi

Fibrous tissue에 사용 가능한 lysis
buffer가 별도로 들어 있습니다.

Solution type

Vacuum방법에 필요한 manifold,
eluator 그리고 pump까지 모두
포함되어 있습니다.



타Br 구매하세요?
위의 X-tracta를 같이
구매 드립니다!!

어머! 이런건
이때 사다 들으세요~

최고 성능의 Miniprep kit!
50% 할인가로
경험해 보세요.

50% 할인할 때
프로메가 마커 한번
사용해 보세요~



Dye-based ←

Probe-based ←

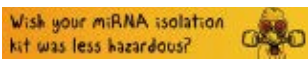
RT mix에는 RNasin Plus
RNase inhibitor가 포함되어 있어.
따로 inhibitor 처리를 하지 않아도 되요~

기존에 M-MLV, AMV, ImProm-II를
사용했다면, 성능과 편리성이 향상된
GoScript RT mix를 사용해 보시는 건 어떨까요?

이 제품은 RT를 15분만에 할 수 있는
Great faster 제품입니다!!

요 제품은~
One-step RT-PCR 제품입니다~

Inhibitor를 사용하시다니!!
철저히 RNA 보로를 하시는군요? 그뤼잇!
그런데 RT mix에는 inhibitor가 포함되어 있어
별도로 사용할 필요가 없다는 사실!!를 아시나요?



유기용매를 사용하지 않고,
miRNA를 다양한 샘플에서
분리정제할 수 있는 제품입니다.
또한 miRNA 이외의 small RNA 분리
정제용으로도 사용할 수 있어요.

Cat. No.	제품명	소비자가	이벤트가	할인율	
Real-time PCR					
A6001	GoTaq® qPCR Master Mix, 5ml (5 x 1ml)	294,000	206,000	30%	
A6002	GoTaq® qPCR Master Mix, 25ml (25 x 1ml)	1,375,000	963,000		
A6010	GoTaq® 2-Step RT-qPCR System 200 rxn	624,000	437,000		
A6020	GoTaq® 1-Step RT-qPCR System, 200 rxn	684,000	479,000		
A6100	GoTaq® Probe qPCR Master Mix, 100 rxn	65,000	46,000		
A6101	GoTaq® Probe qPCR Master Mix, 200 rxn	168,000	118,000		
A6102	GoTaq® Probe qPCR Master Mix, 1,000 rxn	731,000	512,000		
A6110	GoTaq® Probe 2-Step RT-qPCR System, 200 rxn	400,000	280,000		
A6120	GoTaq® Probe 1-Step RT-qPCR System, 200 rxn	253,000	178,000		
Reverse Transcription enzyme					
A2790	GoScript™ Reverse Transcription Mix, Oligo(dT), 50 rxn	326,000	245,000		25%
A2791	GoScript™ Reverse Transcription Mix, Oligo(dT), 100 rxn	436,000	327,000		
A2800	GoScript™ Reverse Transcription Mix, Random Primers, 50 rxn	326,000	245,000		
A2801	GoScript™ Reverse Transcription Mix, Random Primers, 100 rxn	436,000	247,000		
A5000	GoScript™ Reverse Transcription System, 50 rxn	326,000	261,000	20%	
A5001	GoScript™ Reverse Transcription System, 100 rxn	628,000	503,000		
A5003	GoScript™ Reverse Transcriptase, 100 rxn	187,000	150,000		
A5004	GoScript™ Reverse Transcriptase, 500 rxn	720,000	576,000		
A3800	ImProm-II™ Reverse Transcription System, 100 rxn	718,000	575,000		
A3802	ImProm-II™ Reverse Transcriptase, 100 rxn	198,000	159,000		
A3803	ImProm-II™ Reverse Transcriptase, 500 rxn	780,000	624,000		
M1701	M-MLV Reverse Transcriptase, 10,000u	99,000	80,000		
M1705	M-MLV Reverse Transcriptase, 50,000u	374,000	300,000		
M5101	AMV Reverse Transcriptase, 300u	148,000	119,000		
M5108	AMV Reverse Transcriptase, 1,000u	398,000	319,000		
A3500	Reverse Transcription System	719,000	576,000		
One-step RT-PCR					
A1701	AccessQuick™ RT-PCR System, 20 rxn	184,000	129,000	30%	
A1702	AccessQuick™ RT-PCR System, 100 rxn	707,000	495,000		
RNase inhibitor					
N2111	RNasin® RNase Inhibitor, 2,500u	187,000	131,000	30%	
N2115	RNasin® RNase Inhibitor 10,000u	491,000	344,000		
N2511	Recombinant RNasin®RNase Inhibitor, 2,500u	162,000	114,000		
N2515	Recombinant RNasin®RNase Inhibitor, 10,000u	415,000	291,000		
N2611	RNasin® Plus RNase Inhibitor, 2500u	164,000	115,000		
N2615	RNasin® Plus RNase Inhibitor, 10,000u	494,000	346,000		
RT Primer					
C1101	Oligo(dT) 15 Primer, 500ug/ml 20ug	83,000	67,000	20%	
C1181	Random Primers, 20ug	53,000	43,000		
RNA purification_DNase					
M6101	RQ1 RNase-Free DNase, 1,000u	76,000	61,000	20%	
RNA purification					
Z6011	ReliaPrep™ RNA Cell Miniprep System, 50 preps	308,000	231,000	25%	
Z6012	ReliaPrep™ RNA Cell Miniprep System, 250 preps	1,375,000	1,032,000		
Z6111	ReliaPrep™ RNA Tissue Miniprep System, 50 preps	308,000	231,000		
Z6112	ReliaPrep™ RNA Tissue Miniprep System, 250 preps	1,375,000	1,032,000		
Z1002	ReliaPrep™ FFPE Total RNA Miniprep System, 100 preps	881,000	661,000		
Z6211	ReliaPrep™ miRNA Cell and Tissue Miniprep System, 50 preps	460,000	345,000		
Z3100	SV Total RNA Isolation System, 50 preps	294,000	236,000	20%	
Z3105	SV Total RNA Isolation System, 250 preps	1,307,000	1,046,000		

그뤼잇!



ReliaPrep RNA Kit의 column은 이렇게 생겼어요!
Membrane의 직경이 좀더 10-15μm 농축을 할 수 있습니다~



모든 G2 Polymerase는
30% 할인입니다~

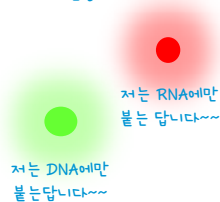
아시나요?

모든 GoTaq 제품에는 gel loading dye가
포함된 Green buffer와 loading dye가 없는
Colorless buffer 두 종류가 들어 있습니다.
실험 목적에 맞게 요렇게 저렇게 사용하세요~

아~ 그래서 master mix가 Green과
Colorless 두 종류였군요? ㅎㅎ



Purity 확인은 흡광방법으로!
정량은 형광방법인 Quantus로
진행하세요~



Cat. No.	제품명	소비자가	이벤트가	할인율
Taq DNA polymerase				
M7841	GoTaq® G2 DNA Polymerase, 100u	49,000	35,000	
M7845	GoTaq® G2 DNA Polymerase, 500u	151,000	106,000	
M7848	GoTaq® G2 DNA Polymerase, 2,500u	661,000	463,000	
M7801	GoTaq® G2 Flexi DNA Polymerase, 100u	49,000	35,000	
M7805	GoTaq® G2 Flexi DNA Polymerase, 500u	151,000	106,000	
M7806	GoTaq® G2 Flexi DNA Polymerase, 2,500u	661,000	463,000	
M7808	GoTaq® G2 Flexi DNA Polymerase, 10,000u	2,240,000	1,568,000	
M7822	GoTaq® G2 Green Master Mix, 100 rxn	54,000	38,000	
M7823	GoTaq® G2 Green Master Mix, 1,000 rxn	448,000	314,000	
M7832	GoTaq® G2 Colorless Master Mix, 100 rxn	54,000	38,000	30%
M7833	GoTaq® G2 Colorless Master Mix, 1,000 rxn	448,000	314,000	
M7401	GoTaq® G2 Hot Start Polymerase, 100u	81,000	57,000	
M7405	GoTaq® G2 Hot Start Polymerase, 500u	252,000	177,000	
M7406	GoTaq® G2 Hot Start Polymerase, 2,500u	1,127,000	789,000	
M7408	GoTaq® G2 Hot Start Polymerase, 10,000u	4,200,000	2,940,000	
M7422	GoTaq® G2 Hot Start Green Master Mix, 100 rxn	90,000	63,000	
M7423	GoTaq® G2 Hot Start Green Master Mix, 1,000 rxn	709,000	497,000	
M7432	GoTaq® G2 Hot Start Colorless Master Mix, 100 rxn	90,000	63,000	
M7433	GoTaq® G2 Hot Start Colorless Master Mix, 1,000 rxn	709,000	497,000	
M3001	GoTaq® DNA Polymerase, 100u	67,000	54,000	
M3005	GoTaq® DNA Polymerase, 500u	258,000	207,000	
M3008	GoTaq® DNA Polymerase, 2500u	1,220,000	976,000	
M8291	GoTaq® Flexi DNA Polymerase, 100u	51,000	41,000	
M8295	GoTaq® Flexi DNA Polymerase, 500u	212,000	170,000	
M8296	GoTaq® Flexi DNA Polymerase, 2500u	940,000	752,000	
M7122	GoTaq® Green Master Mix, 100 rxn	65,000	52,000	
M7123	GoTaq® Green Master Mix, 1,000 rxn	543,000	435,000	
M7132	GoTaq® Colorless Master Mix, 100 rxn	65,000	52,000	20%
M7133	GoTaq® Colorless Master Mix, 1,000 rxn	543,000	435,000	
M5001	GoTaq® Hot Start Polymerase, 100u	107,000	86,000	
M5005	GoTaq® Hot Start Polymerase, 500u	464,000	372,000	
M5008	GoTaq® Hot Start Polymerase, 10,000u	6,790,000	5,432,000	
M5122	GoTaq® Hot Start Green Master Mix, 100 rxn	145,000	116,000	
M5123	GoTaq® Hot Start Green Master Mix, 1,000 rxn	1,146,000	917,000	
M5132	GoTaq® Hot Start Colorless Master Mix, 100 rxn	145,000	116,000	
M5133	GoTaq® Hot Start Colorless Master Mix, 1,000 rxn	1,146,000	917,000	
Special Taq				
M4021	GoTaq® Long PCR Master Mix	241,000	169,000	30%
M7741	Pfu DNA Polymerase, 100u	159,000	128,000	20%
M7745	Pfu DNA Polymerase, 500u	586,000	469,000	
Quantitation				
E2670	QuantiFluor® dsDNA System	377,000	283,000	
E3190	QuantiFluor® ssDNA System	425,000	319,000	
E3310	QuantiFluor® RNA System	425,000	319,000	
E4870	QuantiFluor® ONE dsDNA System, 500rxn	438,000	329,000	25%
E4871	QuantiFluor® ONE dsDNA System, 100rxn	149,000	112,000	
E4942	0.5mL PCR Tubes, 200 pack	104,000	78,000	
E6150	Quantus™ Fluorometer	2,839,000	2,130,000	
E5150	Quantus™ NGS Starter Package	2,860,000	2,145,000	

Restriction enzyme 전품목 50% 할인





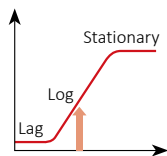
몸에 해로운 EtBr 대신 사용할 수 있는
fluorescent dye예요~~기존 사용하셨던
UV Transilluminator로 확인할 수 있습니다!

SDS-polyacrylamide gel
만들 때 필요한 chemical 입니다~~
미리미리 구비해 두세요!!



Western blot.
washing할 때 필요하시죠? <--

앗! Induction 해야 할 시간이에요!! <--



프로메가에도 G-418이 있습니다~~
모르셨죠? <--

Cat. No.	제품명	소비자가	이벤트가	할인율
Biochemical				
V3121	Agarose, LE, Analytical Grade, 100g	255,000	128,000	
V3125	Agarose, LE, Analytical Grade, 500g	689,000	345,000	
V2111	Agarose, LMP, Analytical Grade, 25g	237,000	119,000	
V2831	Agarose, LMP, Preparative Grade for Large Fragments, 25g	323,000	162,000	
V3841	Agarose, LMP, Preparative Grade for Small Fragments, 25g	209,000	105,000	
H1181	Diamond™ Nucleic Acid Dye, 500ul	105,000	53,000	
H5001	Boric Acid, 500g	94,000	47,000	
H5003	Boric Acid 1000g	148,000	74,000	
H5031	EDTA Disodium Salt 100g	73,000	37,000	
H5032	EDTA Disodium Salt 500g	192,000	96,000	
H5051	Formamide, 100ml	102,000	51,000	
H5052	Formamide, 500ml	218,000	109,000	
H5071	Glycine 500g	126,000	63,000	
H5073	Glycine, 1kg	167,000	84,000	
H5113	Sodium Dodecyl Sulfate(SDS), 100g	111,000	56,000	50%
H5114	Sodium Dodecyl Sulfate(SDS), 500g	253,000	127,000	
H5115	Sodium Dodecyl Sulfate(SDS), 1kg	549,000	275,000	
V6551	Sodium Dodecyl Sulfate(SDS) Solution, 100ml	64,000	32,000	
V6553	Sodium Dodecyl Sulfate(SDS) Solution, 500ml	294,000	147,000	
H5121	Tris-HCl (Tris-Hydrochloride), 100g	85,000	43,000	
H5123	Tris-HCl (Tris-Hydrochloride), 500g	196,000	98,000	
H5125	Tris-HCl (Tris-Hydrochloride), 2500g	867,000	434,000	
V3111	Acrylamide Molecular Grade, 100g	83,000	42,000	
V3115	Acrylamide Molecular Grade, 500g	250,000	125,000	
V3131	Ammonium Persulfate, 25g (APS)	34,000	17,000	
V3141	Bisacrylamide, 25g (N,N'-Methylenebisacrylamide)	52,000	26,000	
V3143	Bisacrylamide, 125g (N,N'-Methylenebisacrylamide)	162,000	81,000	
H5151	Tween® 20, 500ml	143,000	72,000	
H5152	Tween® 20, 100ml	97,000	49,000	
H5131	Tris Base, 500g	146,000	73,000	
H5133	Tris Base, 100g	46,000	23,000	
H5135	Tris Base, 2,500g	558,000	279,000	
H5141	Triton® X-100 500ml	161,000	81,000	
H5142	Triton® X-100 100ml	79,000	40,000	
H5252	Ammonium Sulfate 5kg	228,000	114,000	
H5271	Sodium Chloride 500g	60,000	30,000	
H5273	Sodium Chloride 1kg	105,000	53,000	
H5302	HEPES (free acid), 100g	116,000	58,000	
H5303	HEPES (free acid), 500g	471,000	236,000	
H5433	Glycerol, 1000ml	89,000	45,000	
S2011	Coelenterazine h, 250ug	193,000	97,000	
V3011	PEG 8000 Powder (Polyethylene Glycol), 500g	122,000	61,000	
V3151	DTT, 5g (Dithiothreitol)	142,000	71,000	
V3155	DTT, 25g (Dithiothreitol)	508,000	254,000	
V3171	Urea, 1kg	105,000	53,000	50%
V3175	Urea, 5kg	344,000	172,000	
V3951	IPTG, Dioxane-Free, 5g	324,000	162,000	
V3955	IPTG, Dioxane-Free, 1g	87,000	44,000	
V4221	5M Sodium Chloride, 1,000ml	71,000	36,000	
V4231	0.5M EDTA, pH 8.0, 100ml	22,000	11,000	
V4233	0.5M EDTA, pH 8.0, 400ml	88,000	44,000	
V4251	TBE Buffer, 10X Molecular Biology Grade, 1,000ml	97,000	49,000	
V4261	SSC Buffer, 20X, 1L	108,000	54,000	
V4271	TAE Buffer, 10X, 1,000ml	91,000	46,000	
V4281	TAE, 40X (Tris-acetate-EDTA), 1,000ml	149,000	75,000	
V6231	TE Buffer, 1X, 100ml	49,000	25,000	
V6232	TE Buffer, 1X, 500ml	194,000	97,000	
V7983	Antibiotic G-418 Sulfate, 5g	739,000	370,000	
V8091	Antibiotic G-418 Sulfate Solution, 20ml	232,000	116,000	





BEST

MTS assay

MTT assay

BEST

세포 환원력을 이용하여 최대 72시간까지 real-time으로 세포의 Viability를 발광 기법으로 측정하는 제품입니다.

LDH assay

DNA 염색하는 형광 dye 기법으로 최대 72시간까지 real-time으로 세포 독성 변화 모니터링이 가능합니다.

PS-Annexin V 결합을

microplate에서 최대 48시간까지 real-time으로 모니터링할 수 있습니다.

TIP!! "Glo" 제품은 lysis 능력이 시약 내에 이미 포함되어 있는 별도의 lysis 단계가 필요 없는 homogenous 방법으로 발광 signal이 긴 제품입니다.

Cat. No.	제품명	소비자가	이벤트가	할인율
Cell Viability Assay				
G3580	CellTiter 96® AQueous One Solution Cell Proliferation Assay, 1,000 assays	382,000	249,000	
G3581	CellTiter 96® AQueous One Solution Cell Proliferation Assay, 5,000 assays	1,260,000	819,000	
G5421	CellTiter 96® AQueous Non-Radioactive Cell Proliferation Assay, 1,000 assays	345,000	225,000	35%
G5430	CellTiter 96® AQueous Non-Radioactive Cell Proliferation Assay, 5,000 assays	1,140,000	741,000	
G4000	CellTiter 96® Non-Radioactive Cell Proliferation Assay, 1,000 assays	374,000	244,000	
G4100	CellTiter 96® Non-Radioactive Cell Proliferation Assay, 5,000 assays	1,110,000	722,000	
G7570	CellTiter-Glo® Luminescent Cell Viability, 10ml	160,000	128,000	
G7571	CellTiter-Glo® Luminescent Cell Viability, 10 x 10ml	728,000	583,000	
G7572	CellTiter-Glo® Luminescent Cell Viability, 100ml	576,000	461,000	
G7573	CellTiter-Glo® Luminescent Cell Viability, 10 x 100ml	4,712,000	3,770,000	
G9241	CellTiter-Glo® 2.0 Assay, 10ml	169,000	136,000	20%
G9242	CellTiter-Glo® 2.0 Assay, 100ml	620,000	496,000	
G9243	CellTiter-Glo® 2.0 Assay, 500ml	2,630,000	2,104,000	
G9681	CellTiter-Glo® 3D Cell Viability Assay, 10ml	173,000	139,000	
G9682	CellTiter-Glo® 3D Cell Viability Assay, 10 x 10ml	803,000	643,000	
G9683	CellTiter-Glo® 3D Cell Viability Assay, 100ml	624,000	500,000	
G9711	RealTime-Glo™ MT Cell Viability Assay, 100 reactions	213,000	160,000	
G9712	RealTime-Glo™ MT Cell Viability Assay, 10 x 100 reactions	1,040,000	780,000	25%
G9713	RealTime-Glo™ MT Cell Viability Assay, 1,000 reactions	941,000	706,000	
Cytotoxicity Assay				
G1780	CytoTox 96® Non-Radioactive Cytotoxicity Assay, 1,000 assays	405,000	284,000	30%
G8741	CellTox™ Green Cytotoxicity Assay, 10ml	159,000	120,000	
G8742	CellTox™ Green Cytotoxicity Assay, 50ml	580,000	435,000	25%
G8743	CellTox™ Green Cytotoxicity Assay, 100ml	923,000	693,000	
G8731	CellTox™ Green Express Cytotoxicity Assay, 200ul	867,000	651,000	
Apoptosis Assay				
G8090	Caspase-Glo® 3/7 Assay, 2.5ml	181,000	145,000	
G8091	Caspase-Glo® 3/7 Assay, 10ml	566,000	453,000	
G8092	Caspase-Glo® 3/7 Assay, 100ml	2,960,000	2,368,000	
G8200	Caspase-Glo® 8 Assay, 2.5ml	181,000	145,000	20%
G8201	Caspase-Glo® 8 Assay, 10ml	566,000	453,000	
G8210	Caspase-Glo® 9 Assay, 2.5ml	181,000	145,000	
G8211	Caspase-Glo® 9 Assay, 10ml	566,000	453,000	
JA1011	RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay, 100 assays	681,000	511,000	
JA1012	RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay, 1,000 assays	3,405,000	2,554,000	25%
JA1000	RealTime-Glo™ Annexin V Apoptosis Assay, 100 assays	585,000	439,000	
JA1001	RealTime-Glo™ Annexin V Apoptosis Assay, 1,000 assays	2,925,000	2,194,000	
G3250	DeadEnd™ Fluorometric TUNEL System, 60 reactions	780,000	624,000	
G7130	DeadEnd™ Colorimetric TUNEL System, 40 reactions	768,000	615,000	20%
G7360	DeadEnd™ Colorimetric TUNEL System, 20 reactions	557,000	446,000	

RealTime-Glo™ Annexin V Apoptosis & Necrosis Assay

Healthy Cell
PS confined to inner leaflet
Cell membrane intact
Luminescence (RLU) negative
Fluorescence (RFU) negative

Early Apoptosis
PS translocation to outer leaflet
Cell membrane intact
Luminescence (RLU) POSITIVE
Fluorescence (RFU) negative

Secondary Necrosis
PS on inner and outer leaflet
Cell membrane compromised
Luminescence (RLU) POSITIVE
Fluorescence (RFU) POSITIVE

RealTime-Glo™ Annexin V Assay 원리

Detection 시약에는 NanoLuc® luciferase의 complementary domain을 포함한 두 개의 Annexin V fusion protein (Annexin V-LgBIT & Annexin V-SmBIT)이 포함되어 있습니다.

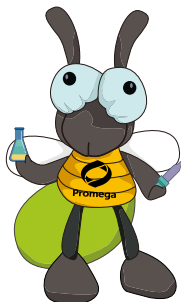
또한 세포막을 투과하지 못하는 pro-fluorescent DNA dye도 포함되어 있습니다.

발광 (luminescence)은 PS와 Annexin V fusion 단백질이 결합 (functional luciferase 형성)할 때까지 낮게 유지됩니다.

형광 (fluorescent signal)은 late apoptosis 단계에서 세포막 손실에 의해 생성됩니다.

* Assay 제품명 앞의 바 색깔에 따라 █ 흡광, █ 형광, █ 발광 assay 기법을 확인하실 수 있습니다.

Cat. No.	제품명	소비자가	이벤트가	할인율	
Luciferase Assay					
BEST Firefly luciferase assay	E1500	Luciferase Assay System, 100 assays	142,000	114,000	
	E1501	Luciferase Assay System, 10-Pack, 1,000 assays	919,000	736,000	
	E4030	Luciferase Assay System with Reporter Lysis Buffer, 100 assays	144,000	116,000	
	E6110	ONE-Glo™ Luciferase Assay System, 10ml	202,000	162,000	
	E6120	ONE-Glo™ Luciferase Assay System, 100ml	1,300,000	1,040,000	
	E8110	ONE-Glo™ EX Luciferase Assay System, 10ml	202,000	162,000	
	E8120	ONE-Glo™ EX Luciferase Assay System, 100ml	1,170,000	936,000	
	E8130	ONE-Glo™ EX Luciferase Assay System, 10 x 10ml	1,400,000	1,120,000	
	E2510	Steady-Glo® Luciferase Assay System, 10ml	149,000	120,000	
	E2520	Steady-Glo® Luciferase Assay System, 100ml	1,100,000	880,000	
E2610	Bright-Glo™ Luciferase Assay System, 10ml	189,000	152,000		
E2620	Bright-Glo™ Luciferase Assay System, 100ml	1,130,000	904,000	20%	
NanoLuc® luciferase assay	N1110	Nano-Glo® Luciferase Assay, 10ml	207,000	166,000	
	N1120	Nano-Glo® Luciferase Assay, 100ml	1,175,000	940,000	
	N1130	Nano-Glo® Luciferase Assay, 10 x 10ml	1,410,000	1,128,000	
Firefly & Renilla (control) Dual luciferase assay	E1910	Dual-Luciferase® Reporter Assay System, 100 assays	302,000	242,000	
	E1960	Dual-Luciferase® Reporter Assay System 10-Pack, 1,000 assays	1,929,000	1,544,000	
	E1980	Dual-Luciferase® Reporter 1000 Assay System, 1,000 assays	1,778,000	1,423,000	
	E2920	Dual-Glo® Luciferase Assay System, 10ml	334,000	268,000	
Firefly & NanoLuc Dual luciferase assay	E2940	Dual-Glo® Luciferase Assay System, 100ml	1,990,000	1,592,000	
	N1610	Nano-Glo® Dual-Luciferase® Reporter Assay System, 10ml	315,000	252,000	
	N1620	Nano-Glo® Dual-Luciferase® Reporter Assay System, 100ml	1,816,000	1,453,000	
N1630	Nano-Glo® Dual-Luciferase® Reporter Assay System, 10 x 10ml	1,956,000	1,565,000		
Transfection					
E2311	FuGENE® HD Transfection Reagent, 1ml	612,000	459,000		
E2312	FuGENE® HD Transfection Reagent, 5 x 1ml	2,775,000	2,082,000		
E2691	FuGENE® 6 Transfection Reagent, 1ml	562,000	422,000		
E2692	FuGENE® 6 Transfection Reagent, 5 x 1ml	2,370,000	1,778,000	25%	
E2693	FuGENE® 6 Transfection Reagent, 0.5ml	338,000	254,000		
E4981	ViaFect™ Transfection Reagent, 0.75ml	408,000	306,000		
E4982	ViaFect™ Transfection Reagent, 2 x 0.75ml	679,000	510,000		
in vivo Imaging					
E1602	Beetle Luciferin, Potassium Salt, 50mg	271,000	190,000	30%	
E1603	Beetle Luciferin, Potassium Salt, 250mg	554,000	388,000		
E1605	Beetle Luciferin, Potassium Salt, 1g	845,000	423,000	50%	
P1041	VivoGlo™ Luciferin, In Vivo Grade, 50mg	279,000	196,000	30%	
P1042	VivoGlo™ Luciferin, In Vivo Grade, 250mg	567,000	397,000		
P1043	VivoGlo™ Luciferin, In Vivo Grade, 1g	966,000	483,000	50%	
Protease					
V5111	Sequencing Grade Modified Trypsin, 5 x 20ug	133,000	107,000		
V5113	Sequencing Grade Modified Trypsin, Frozen, 5 x 20ug	66,000	53,000		
V5117	Sequencing Grade Modified Trypsin, 1 x 100ug	119,000	96,000		
V5280	Trypsin Gold, Mass Spectrometry Grade, 100ug	179,000	144,000		
V5072	Trypsin/Lys-C Mix, Mass Spec Grade, 1 x 100ug	251,000	201,000		
V5073	Trypsin/Lys-C Mix, Mass Spec Grade, 5 x 20ug	271,000	217,000		
V1671	rLys-C, Mass Spec Grade, 15ug	217,000	174,000		
V1061	Chymotrypsin, Sequencing Grade, 25ug	109,000	88,000	20%	
V1062	Chymotrypsin, Sequencing Grade, 100ug	336,000	269,000		
V1621	Asp-N, Sequencing Grade, 2ug	245,000	196,000		
V1651	Glu-C, Sequencing Grade, 50ug	245,000	196,000		
V1881	Arg-C, Sequencing Grade, 10ug	345,000	276,000		
V4831	PNGase F, 500 units	208,000	167,000		
V3021	Proteinase K, 100mg	187,000	150,000		
2nd Antibody					
W4011	Anti-Rabbit IgG (H+L), HRP Conjugate, 300ul	163,000	82,000	50%	
W4021	Anti-Mouse IgG (H+L), HRP Conjugate, 300ul	165,000	83,000		
ECL Solution					
W1001	ECL Western Blotting Substrate, 250ml	248,000	186,000	25%	
W1015	ECL Western Blotting Substrate, 500ml	344,000	224,000	35%	



Mass Spec Grade 제품인 Trypsin Gold와 rLys-C mixture 제품으로 proteolysis 기능이 증가되어 재현성 높은 실험 결과를 제공합니다.



빠르고 편리하며 정확하게
plate-based caspase 1
활성 측정이 가능합니다.

Cat. No.	제품명	소비자가	이벤트가	할인을
Inflammation Assay				
G9951	Caspase-Glo® 1 Inflammasome Assay, 10ml	630,000	504,000	20%
G9952	Caspase-Glo® 1 Inflammasome Assay, 5 x 10ml	2,675,000	2,140,000	
Oxidative Stress Assay				
V6611	GSH/GSSG-Glo™ Assay, 10ml	716,000	573,000	20%
V6612	GSH/GSSG-Glo™ Assay, 50ml	2,870,000	2,296,000	
V6911	GSH-Glo™ Glutathione Assay, 10ml	589,000	472,000	
V6912	GSH-Glo™ Glutathione Assay, 50ml	2,300,000	1,840,000	
G8820	ROS-Glo™ H202 Assay, 10ml	540,000	432,000	
G8821	ROS-Glo™ H202 Assay, 50ml	2,260,000	1,808,000	
G2930	Griess Reagent System	237,000	190,000	
Energy Metabolism Assay				
G9071	NAD/NADH-Glo™ Assay, 10ml	651,000	521,000	20%
G9072	NAD/NADH-Glo™ Assay, 50ml	2,647,000	2,118,000	
G9081	NADP/NADPH-Glo™ Assay, 10ml	651,000	521,000	
G9082	NADP/NADPH-Glo™ Assay, 50ml	2,647,000	2,118,000	
J6021	Glucose-Glo™ Assay, 5ml	540,000	432,000	
J6022	Glucose-Glo™ Assay, 50ml	2,290,000	1,832,000	
J5021	Lactate-Glo™ Assay, 5ml	598,000	479,000	
J5022	Lactate-Glo™ Assay, 50ml	2,484,000	1,988,000	
J7021	Glutamate-Glo™ Assay, 5ml	598,000	479,000	
J7022	Glutamate-Glo™ Assay, 50ml	2,484,000	1,988,000	
J8021	Glutamine/Glutamate-Glo™ Assay, 5ml	641,000	513,000	
J8022	Glutamine/Glutamate-Glo™ Assay, 50ml	2,844,000	2,276,000	
J1341	Glucose Uptake-Glo™ Assay, 5ml	504,000	404,000	
J1342	Glucose Uptake-Glo™ Assay, 10ml	871,000	697,000	
J1343	Glucose Uptake-Glo™ Assay, 50ml	3,485,000	2,788,000	

* Assay 제품명 앞의 바 색깔에 따라 ■ 흡광, ■ 형광, ■ 발광 assay 기법을 확인하실 수 있습니다.

논문인증 이벤트



프로메가 제품을 사용한 국내 연구자 분들의 연구 논문을 찾고 있습니다.
Material & Method에 프로메가 제품 사용을 표기한 SCI 논문(First Author, 2016년 이후의 논문)을 네이버 카페 논문인증 게시판에 올려주세요.

1. 제목: 논문 제목
2. 내용: 사용하신 제품을 작성
3. 첨부: 해당 논문을 첨부 파일로 올려주세요

보내주신 연구자분들께 성함을 정성껏 각인한 LAMY 펜을 감사 선물로 보내드립니다.

- 행사기간** 2018년 9월 30일까지, 선착순 30명
- 참여요건** Material & Method에 프로메가 제품 사용을 표기한 SCI 논문
First Author, 2016년 이후의 논문
- 선물** 참여하신 분의 성함을 각인한 LAMY 펜



프로메가 코리아 • www.promega.com • Telephone 1588-3718 FAX 02-2638-5418

제품 기술 문의 TechServiceKR@promega.com • 제품 주문 문의 CustServiceKR@promega.com