

Recombinant Protein Expression: Extraction and Stabilization

Reference and Selection Guide

Efficiently Extract Protein

Preserve Protein Function

Prevent Protein Degradation

INNOVATION @ WORK



Bacterial

Mammalian

Plant

Yeast

sigma.com/lysis

SIGMA[®]

Protease and Phosphatase Inhibitor Cocktails

sigma.com/protinhib

Sample Type	General Use		Bacterial	Mammalian	Plant	Yeast	Histidine-tagged Proteins	Tissue Culture Media	Phosphatase Inhibitors	
									Tyrosine/Alkaline	Serine/Threonine
Catalog Number	P2714	S8820	P8465	P8340	P9599	P8215	P8849	P1860	P5726	P2850
Specificity of Inhibition	Serine, Cysteine, Aspartic, & Metalloproteases	Serine, Cysteine, Aspartic, & Metalloproteases	Serine, Cysteine, Aspartic, & Metalloproteases, & Aminopeptidases	Serine, Cysteine, Aspartic Proteases & Amidopeptidases	Serine, Cysteine, Aspartic and Metalloproteases, & Aminopeptidases	Serine, Cysteine, Aspartic, & Metalloproteases	Serine, Cysteine, Aspartic and Thermolysin-like Proteases & Aminopeptidases	Serine Cysteine, Aspartic Proteases, & Aminopeptidases	Acid and Alkaline Phosphatases, as well as Tyrosine Protein Phosphatases	Serine, Threonine Phosphatases and various isozymes of Alkaline Phosphatase
Cocktail Format	Lyophilized powder that is water soluble	Tablet that is water soluble	Lyophilized powder with a vial of DMSO	Solution in DMSO	Solution in DMSO	Solution in DMSO	Solution in DMSO	Solution in DMSO	Solution in water	Solution in DMSO
Usage (amount of sample that 1 ml of cocktail will inhibit); typically enough for 100 ml of cell lysate, depending on cell type	1 mg of USP pancreatin; typically enough for 100 ml of cell lysate	1 mg of USP pancreatin (1x solution); one tablet generates 100 ml of protease inhibitor solution	20 ml of cell lysate from 4 g (wet weight) of <i>E. coli</i> cells	100 ml of cell lysate from 20 g (wet weight) of bovine liver or 10 ml of cell lysate from CHO cells at 10 ⁶ cells per ml	100 ml of cell lysate from 30 g (wet weight) of plant tissue	100 ml of cell lysate from 20 g (wet weight) of <i>Saccharomyces cerevisiae</i> cells	100 ml of cell lysate from 20 g (wet weight) of <i>E. coli</i> cells or 10 g (wet weight) of baculovirus infected cells	Dilution of ≥1:200 in tissue culture media to prevent proteolytic degradation of secreted proteins	100,000 x g supernatant from human placenta, bovine liver, rabbit muscle, A431 or Jurkat cell extracts at a protein concentration of approximately 5 mg/ml	
Cocktail Components (and concentrations when available)	AEBSF, 2 mM Aprotinin, 0.3 μM Bestatin, 130 μM EDTA, 1 mM E-64, 14 μM Leupeptin, 1 μM	AEBSF, 2 mM Aprotinin, 0.3 μM Bestatin, 130 μM EDTA, 1 mM E-64, 14 μM Leupeptin, 1 μM	AEBSF, 23 mM Bestatin, 2 mM EDTA, 100 mM E-64, 0.3 mM Pepstatin A, 0.3 mM	AEBSF, 104 mM Aprotinin, 80 μM Bestatin, 4 mM E-64, 1.4 mM Leupeptin, 2 mM Pepstatin A, 1.5 mM	AEBSF Bestatin E-64 Leupeptin Pepstatin A 1,10-Phenanthroline	AEBSF, 100 mM E-64, 1.4 mM Pepstatin A, 2.2 mM 1,10-Phenanthroline, 500 mM	AEBSF Bestatin E-64 Pepstatin A Phosphoramidon	Aprotinin Bestatin E-64 Leupeptin Pepstatin A	Imidazole Sodium Molybdate Sodium Orthovanadate Sodium Tartrate	Bromotetramisole Canthardin Microcystin LR
Additional Special / Specific Applications (Tested on but not limited to)	Optimized and tested for general use	Optimized and tested for general use	Various <i>E. coli</i> strains	Use at 1:100 for undiluted blood samples, liver tissue extracts and CHO cell lysates	Tested on extracts from kidney bean, pea, wheat, tobacco, and <i>Arabidopsis</i>	<i>Saccharomyces cerevisiae</i> cells	<i>E. coli</i> cells; baculovirus infected <i>Spodoptera frugiperda</i> pupal ovary cells	After 48 hours in medium, add fresh inhibitors in fresh medium	Tested on various animal tissues (bovine liver, human placenta, rabbit muscle) and A431 and Jurkat cell lines	Tested on various animal tissues (bovine liver, human placenta, rabbit muscle) and A431 and Jurkat cell lines
Storage and Stability (unopened)	-20 °C ≥ 4 years	2-8 °C ≥ 4 years	-20 °C ≥ 4 years	At least 4 years at -20 °C; 8 months at 2-8 °C; and 2 months at RT	At least 4 years at -20 °C; 8 months at 2-8 °C; and 2 months at RT	At least 4 years at -20 °C; 8 months at 2-8 °C; and 2 months at RT	At least 4 years at -20 °C; 3 months at 2-8 °C; and 2 weeks at RT	At least 4 years at -20 °C; 8 months at 2-8 °C; and 2 months at RT	2-8 °C ≥ 2 years	2-8 °C ≥ 4 years

Protease and Phosphatase Inhibitors Available Individually

Inhibitors	Cat. No.	Application
AESF	A8456	Inhibits serine proteases, such as trypsin and chymotrypsin
Aprotinin, saline solution Aprotinin, affinity purified	A6279 A4529	Inhibits serine proteases such as trypsin, chymotrypsin, plasmin, trypsinogen, urokinase, and kallikrein. Also inhibits human leukocyte elastase, but not pancreatic elastase
Bestatin hydrochloride	B8385	Inhibits aminopeptidases, such as leucine aminopeptidase proteases
Cantharidin	C7632	Inhibits protein phosphatase 2A (PP-2A)
Chymostatin	C7268	Inhibits serine and cysteine proteases
E-64	E3132	Inhibits cysteine proteases such as calpain, papain, and cathepsin B
EDTA disodium salt dihydrate EDTA, 0.5 M solution disodium salt	E5134 E7889	Inhibits metalloproteases; chelates; permeabilizes Gram-negative bacteria
EGTA	E8145	Inhibits metalloproteases; chelates
Imidazole	I0125	Inhibits alkaline phosphatases
Leupeptin hemisulfate Leupeptin TFA salt Leupeptin hydrochloride	L2884 L2023 L9783	Inhibits both serine and cysteine proteases such as calpain, trypsin, papain, and cathepsin B
Microcystin LR	M2912	Inhibits protein phosphatases 1 and 2A (PP-1 and PP-2A)
Pepstatin A Pepstatin A, 90% pure	P4265 P5318	Inhibits acid proteases such as pepsin (human or porcine), renin, cathepsin D, chymosin (bovine rennin), and protease B
1,10-Phenanthroline	P9375	Inhibits metalloproteases
Phosphoramidon	R7385	Inhibits thermolysin and collagenase
PMSF	P7626	Inhibits serine proteases
Sodium molybdate	M1003	Inhibits acid and phosphoprotein phosphatases
Sodium orthovanadate	S6508	Inhibits a number of ATPases, protein tyrosine phosphatases, and other phosphate-transferring enzymes
Sodium tartrate dibasic	S4797	Inhibits alkaline phosphatases

Recombinant Protein Expression: Extraction and Stabilization

Organism

BACTERIA

	WHOLE CELL DETERGENT LYSIS			WHOLE CELL DETERGENT LYSIS
Product	CellLytic™ B	CellLytic B 2x	CellLytic B 10x	CellLytic Express
Catalog number	B7435	B7310	C8740	C1990
Description	A ready-to-use solution containing 40 mM Tris-HCl, pH 8, and a proprietary, non-denaturing detergent formulation of zwitterionic detergents	A more concentrated (2x), ready-to-use solution containing 40 mM Tris-HCl, pH 8, and a proprietary, non-denaturing formulation detergent	A highly concentrated (10x) solution containing a proprietary, non-denaturing detergent formulation and no buffering components	A ready-to-use, all-in-one, non-denaturing proprietary powder formulation of detergent and enzymes optimized for one-step extraction of proteins from bacterial cells (in-culture cell lysis)
Package sizes	50 ml, 500 ml (Sufficient for processing 5 g and 50 g of wet cell paste, respectively)	50 ml, 250 ml (Sufficient for processing up to 10 g and 50 g of wet cell paste, respectively)	10 ml, 50 ml, 100 ml (Sufficient for processing ~10 g, 50 g, and 100 g of wet cell paste, respectively)	25 ml (1.26 g) 10 x 25 ml 6 x 500 ml (25.2 g)
Sample type	<ul style="list-style-type: none"> Bacterial cell pellets (lysis and protein extraction) Isolation of inclusion bodies 	<ul style="list-style-type: none"> Bacterial cell pellets (concentrated at 2x for more difficult to lyse organisms) Isolation of inclusion bodies 	<ul style="list-style-type: none"> Bacterial cell pellets (concentrated at 10x in water; dilute for use into buffer of choice) Isolation of inclusion bodies 	<ul style="list-style-type: none"> Direct lysis of bacterial cell cultures (without centrifugation) Affinity purification of proteins without the need for special equipment
Strain/Cell line compatibility	<ul style="list-style-type: none"> Optimized for <i>E. coli</i> strain BL21 DH5α™ JM109 Other similar cells 	<ul style="list-style-type: none"> Optimized for <i>E. coli</i> strain BL21 DH5α JM109 Other similar cells 	<ul style="list-style-type: none"> Optimized for <i>E. coli</i> strain BL21 DH5α JM109 Other similar cells 	<ul style="list-style-type: none"> Optimized for <i>E. coli</i> strain BL21 DH5α JM109 Other similar cells
Usage	<p><u>Small scale:</u> 0.4 ml of CellLytic B per 1.5 ml culture pellet</p> <p><u>Large scale:</u> 10-20 ml per g of wet cell paste</p>	<p><u>Small scale:</u> 0.2 ml of CellLytic B 2x per 1.5 ml of culture pellet</p> <p><u>Large scale:</u> 5 ml per g of wet cell paste</p>	<p>Test 5, 10, and 20-fold dilutions into the desired buffer; add diluted solution at 10-20 ml per g of wet cell paste</p>	<p><u>Small scale:</u> one 25 ml bottle per 25 ml of culture</p> <p><u>Large Scale:</u> one 500 ml bottle for 500 ml of culture</p> <p>For cultures <25 ml, add 0.05 g of powder per ml of culture</p>
Type of detergent/ Dialyzable?	Zwitterionic/Yes	Zwitterionic/Yes	Zwitterionic/Yes	Zwitterionic/Yes
Protein active?	Yes – Suitable for protein activity assays	Yes – Suitable for protein activity assays	Yes – Suitable for protein activity assays	N/A; typically resin-bound protein-activity dependent on isolation/elution methods
Compatible purification and detection products	<ul style="list-style-type: none"> Antibody/EZview™ Affinity Gels FLAG® HIS-Select®/IMAC GST Protease Inhibitor Cocktails Lysozyme solution (L3790) 	<ul style="list-style-type: none"> Antibody/EZview Affinity Gels FLAG HIS-Select/IMAC GST Protease Inhibitor Cocktails Lysozyme solution (L3790) 	<ul style="list-style-type: none"> Antibody/EZview Affinity Gels FLAG HIS-Select/IMAC GST Protease Inhibitor Cocktails Lysozyme solution (L3790) 	<ul style="list-style-type: none"> Antibody/EZview Affinity Gels FLAG HIS-Select/IMAC GST Protease Inhibitor Cocktails Lysozyme solution (L3790)
Quantitation and staining capabilities	<ul style="list-style-type: none"> Protein assay Protein staining 	<ul style="list-style-type: none"> Protein assay Protein staining 	<ul style="list-style-type: none"> Protein assay Protein staining 	<ul style="list-style-type: none"> Protein assay Protein staining
Applications	<ul style="list-style-type: none"> Affinity purification SDS-PAGE Western blotting Dot blotting Immunoprecipitation 	<ul style="list-style-type: none"> Affinity purification SDS-PAGE Western blotting Dot blotting Immunoprecipitation 	<ul style="list-style-type: none"> Affinity purification SDS-PAGE Western blotting Dot blotting Immunoprecipitation 	<ul style="list-style-type: none"> Affinity purification SDS-PAGE Western blotting Dot blotting Immunoprecipitation
Related products containing this reagent	See CB0050 & CB0500 kits	See B7435 & CB0500 kits	See B7435 & CB0500 kits	CellLytic Express Tablets (C5491)
Physical form	Ready-to-use solution	Concentrated solution (2x)	Concentrated solution (10x)	Easy-to-dissolve powder
Storage and stability	Room temperature 2 years	Room temperature 2 years	Room temperature 2 years	-20 °C 2 years

CELL DETERGENT & ENZYMATIC LYSIS		INCLUSION BODY SOLUBILIZATION	INTEGRATED LYSIS & AFFINITY PURIFICATION – WHOLE CELLS		WHOLE CELLS
Cellytic Express Tablets	Cellytic B Plus Kit	Cellytic IB	HIS-Select® iLAP® HC Nickel Coated Plate	HIS-Select iLAP 5 ml Column	Cellytic M
C5491	CB0050 & CB0500	C5236	H9412	H9913	C2978
A ready-to-use, all-in-one, non-denaturing proprietary tablet formulation of detergent and enzymes optimized for one-step extraction of proteins from bacterial cells (in-culture cell lysis)	Kits containing Cellytic B Cell Lysis Reagent (B7435), lysozyme, Benzonase®, and a Protease Inhibitor Cocktail to extract soluble proteins and remove cell debris from inclusion bodies	A ready-to-use solution designed to solubilize inclusion bodies formed in bacteria over-expressing recombinant proteins	A single-use integrated lysis and affinity purification plate that combines cell lysis and purification steps into one. Plate contains proprietary coatings for both lysis and affinity purification of histidine-tagged proteins.	A single-use integrated lysis and affinity purification column designed for one-step purification of histidine-tagged proteins directly from a 5 ml bacterial culture	A ready-to-use solution containing a mild detergent in a bicine buffer designed for efficient whole-cell protein extraction from cultured mammalian cells
25 ea 100 ea	1 kit: 50 ml (CB0050) or 500 ml (CB0500) Cellytic B for processing 5 g or 50 g of cell paste, respectively	25 ml 100 ml	1 each 5 x 1 each	1 each 25 each	50 ml 250 ml
<ul style="list-style-type: none"> • Direct lysis of bacterial cell cultures (without centrifugation) • Affinity purification of proteins without the need for special equipment 	Both Gram-negative and difficult to lyse Gram-positive bacteria	Bacterial inclusion bodies	Bacterial culture (direct lysis, no centrifugation necessary)	Bacterial culture (direct lysis, no centrifugation necessary)	Both suspension and adherent cells (no cell scraping necessary; efficient and rapid cell lysis and protein solubilization)
<ul style="list-style-type: none"> • Optimized for <i>E. coli</i> strain BL21 • DH5α • JM109 • Other similar cells 	<ul style="list-style-type: none"> • Optimized for <i>E. coli</i> strain BL21 • Other similar bacteria 	<ul style="list-style-type: none"> • Various <i>E. coli</i> strains 	<ul style="list-style-type: none"> • Optimized for <i>E. coli</i> strain BL21 • Other similar bacteria 	<ul style="list-style-type: none"> • Optimized for <i>E. coli</i> strain BL21 • Other similar bacteria grown in Terrific Broth 	<ul style="list-style-type: none"> • HeLa • HL-60 • PC-12 • CHO • Bovine Aorta Endothelial Cells (BAEC) • Jurkat • COS • A431
Use 1 tablet for 1 ml of culture; for high OD cultures, additional tablets can be used	10 ml per g of wet cell paste (~250 ml culture); dilute 1:10 for inclusion body isolation	8 ml per g of purified inclusion body pellet	200 µl of overnight bacterial culture per well	5 ml of an overnight bacterial culture per column	125 µl for 10 ⁶ -10 ⁷ cells; 500-1000 µl for a 100 mm plate; 200-400 µl for a 35 mm plate
Zwitterionic/Yes	Zwitterionic/Yes	N/A (strong denaturant in mild neutral buffer)	Zwitterionic/Yes	Zwitterionic/Yes	Zwitterionic/Yes
N/A; resin-bound protein – activity dependent on isolation/elution methods	Yes – Suitable for protein activity assays	N/A; need to renature, therefore, dependent on the nature of the specific protein	Activity dependent on method used for isolation and elution of plate-bound protein	Activity dependent on method used for isolation and elution of resin-bound protein	Yes – Suitable for protein activity assays
<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease Inhibitor Cocktails • Lysozyme solution (L3790) 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease Inhibitor Cocktails • Lysozyme solution (L3790) 	Dependent on the nature of the recombinant protein after dialysis and refolding	<ul style="list-style-type: none"> • HIS-Select/IMAC 	<ul style="list-style-type: none"> • HIS-Select/IMAC 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails • Lysozyme solution (L3790)
<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay – only after renaturation/dialysis • Protein staining – after dialysis 	<ul style="list-style-type: none"> • Protein assay – depending on elution method • Protein staining 	<ul style="list-style-type: none"> • Protein assay – depending on elution method • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining
<ul style="list-style-type: none"> • Affinity purification • SDS-PAGE • Western blotting • Dot blotting • Immunoprecipitation 	<ul style="list-style-type: none"> • Affinity purification • SDS-PAGE • Western blotting • Dot blotting • Immunoprecipitation 	<ul style="list-style-type: none"> • SDS-PAGE (following dialysis of the denaturant) • Protein refolding studies 	<ul style="list-style-type: none"> • Western blotting • Dot blotting 	<ul style="list-style-type: none"> • Western blotting • Dot blotting 	<ul style="list-style-type: none"> • Reporter gene assays • Kinase assays • Phosphatase assays • Immunoassays • DNA-protein interaction assays (gel-shift assays)
Cellytic Express powder (C1990), HIS-Select iLAP 5 ml column (H9913)	N/A	N/A	N/A	N/A; see Cellytic Express Tablet, C5491	CELLMM2, IP0020, IP0010, EHM1, HMP1
Easy-to-dissolve tablets	Solutions ready to mix	Ready-to-use solution	96-well plate coated with proprietary lysis and affinity capture reagents	5 ml chromatography column containing 5 Cellytic Express tablets and 1 HIS-Select resin pellet	Ready-to-use solution
-20 °C 2 years	Cellytic B at room temperature; the rest of the kit at -20 °C 1 year	Room temperature ≥1 year	2-8 °C 4 years	2-8 °C 2 years	Room temperature 2 years

TISSUES	MEMBRANE	NUCLEAR	WHOLE CELLS	
CellLytic MT	CellLytic MEM	CellLytic NUCLEAR™ Extraction Kit	RIPA buffer	Mammalian Cell Lysis Kit
C3228	CE0050	NXTRACT	R0278	MCL1
A ready-to-use solution consisting of a dialyzable mild detergent, bicine, and 150 mM NaCl for efficient extraction of tissue proteins for analysis	A kit containing Lysis and Separation Buffer, Wash Buffer, 4 M NaCl Solution, and Protease Inhibitor Cocktail for isolation of hydrophobic and raft-associated proteins	A complete system for preparing nuclear and cytoplasmic protein extracts from mammalian cultured cells or tissues, allowing extraction of functional nuclear proteins	A ready-to-use solution for general cell lysis, which contains 50 mM Tris-HCl, pH 8, 150 mM NaCl, 1% IGEPAL® CA-630, 0.5% sodium deoxycholate, and 0.1% SDS	A kit containing all reagents required for the solubilization of mammalian cells and tissue for subsequent protein analysis. Contains: 5x buffer (Tris-EDTA), 5x NaCl, 5x SDS, 5x deoxycholic acid, 5x IGEPAL CA-630, and protease inhibitor cocktail
50 ml 500 ml	1 kit: 50 ml of lysis and separation buffer, sufficient for 80 samples	1 kit: 100 extractions of 100 µl of packed cells (10 ⁶ -10 ⁷ cells), or 10 extractions of 1 ml of packed cells (10 ⁸ -10 ¹⁰ cells)	50 ml 500 ml	1 kit: extraction of 250 ml of total cell suspension, or extraction of cells from 250 plates (100 mm diameter)
<ul style="list-style-type: none"> • Mammalian tissues • Mammalian cell lines (for efficient extraction of proteins that require higher salt in the lysis buffer) 	Cultured mammalian cells (isolation based on phase separation, not high-speed centrifugation)	Mammalian cells and fresh or frozen tissues (Reagents allow for flexible, application-specific methods of extracting functional nuclear proteins)	Both adherent and suspension cultured mammalian cells (Rapid, efficient cell lysis and solubilization of proteins)	Mammalian cell lines (same application as RIPA buffer (R0278))
<ul style="list-style-type: none"> • Rat tissue (brain, liver, muscle, kidney, heart, and spleen) • Mouse tissue (brain, kidney, and muscle) 	<ul style="list-style-type: none"> • HeLa • HEK-293 • NIH3T3 • COS • CHO 	<ul style="list-style-type: none"> • HeLa • Jurkat • Rat liver • Bovine Aorta Endothelial Cells (BAEC) • Mouse brain and lung • Rabbit muscle • CHO • PC-12 • COS 	<ul style="list-style-type: none"> • Mammalian cell lines. 	<ul style="list-style-type: none"> • HeLa • COS • PC-12 • Bovine Aorta Endothelial Cells (BAEC) • CHO • Jurkat
20 ml per g of tissue (1:20 ratio)	600 µl of lysis and separation buffer for every 10 ⁶ -10 ⁷ cells	500 µl of 1x lysis buffer for every 100 µl of packed cell volume (5x volume)	1 ml of RIPA Buffer is sufficient to lyse cells from one 100 mm culture dish (0.5 to 5 x 10 ⁷ cells)	1 ml per sample For adherent or suspension cells, use 10 ⁶ -10 ⁷ cells/ml; for tissue, use 5-20 mg of tissue/ml
Zwitterionic/Yes	Non-ionic/No	Non-ionic/No – only partially	Non-ionic/No – only partially Anionic/Yes	Non-ionic/No – only partially Anionic/Yes
Yes – Suitable for protein activity assays	Yes – Suitable for protein activity assays	Yes – Suitable for protein activity assays	N/A	N/A
<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails • Lysozyme solution (L3790) 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease Inhibitor Cocktails
<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay – after sample is diluted • Protein staining 	<ul style="list-style-type: none"> • Protein assay – if sample is diluted • Protein staining
<ul style="list-style-type: none"> • Reporter gene assays • Kinase assays • Phosphatase assays • Immunoassays • DNA-protein interaction assays (gel-shift assays) 	<ul style="list-style-type: none"> • SDS-PAGE • Western blotting • Dot blotting • Immunoprecipitation 	<ul style="list-style-type: none"> • Electrophoresis Mobility Shift Assay (EMSA) • DNase I footprinting analysis • Transcription assays • Further purification 	<ul style="list-style-type: none"> • Affinity purification 	<ul style="list-style-type: none"> • Western blotting • Dot blotting • Immunoprecipitation • Affinity purification
N/A	N/A	N/A	N/A	N/A
Ready-to-use solution	Solutions ready to mix	Solutions ready to mix	Ready-to-use solution	Concentrated solutions ready to mix
Room temperature 2 years	-20 °C; after opening the kit, store the Protease Inhibitor Cocktail at -20 °C and the rest at 2-8 °C 2 years	-20 °C; see bulletin for storage of individual components when opened 3 years	2-8 °C ≥ 2 years	-20 °C 2 years

PLANT

YEAST

Organism

WHOLE CELLS & TISSUES	NUCLEAR	DETERGENT	DETERGENT & ENZYMATIC LYSIS	Organism
CellLytic P	CellLytic PN Kit	CellLytic Y	CellLytic Y Plus Kit	Product
C2360	CELLYTPN1	C4482	CYP1	Catalog number
A ready-to-use non-ionic detergent formulation for convenient and efficient plant cell lysis and protein solubilization	A kit containing complete reagents for rapid isolation of nuclei and extraction of functional nuclear proteins from plant leaves	A phosphate-free and non-denaturing reagent for effective yeast cell lysis and protein solubilization	A kit for highly efficient protein extraction from yeast cells, which contains Reaction Buffer, DTT, Lyticase, Extraction Buffer, Triton® X-100, and Protease Inhibitor cocktail	Description
50 ml 250 ml	1 kit: 30 extractions from 20 g of fresh or frozen plant leaves	50 ml 500 ml	1 kit: spheroplast formation and protein extraction from 10-40 g of yeast cell paste	Package sizes
Fresh or frozen plant leaves (Extraction and solubilization of protein from less than one gram to hundreds of grams)	Plant leaves (Isolation of nuclei and extraction of nuclear proteins from a few grams to hundreds of grams)	Fast and efficient lysis and solubilization of proteins from yeast cells	This kit provides a convenient and highly efficient method of spheroplast formation and protein extraction from yeast cells	Sample type
<ul style="list-style-type: none"> • Tobacco • Tomato • Spinach • <i>Arabidopsis</i> 	<ul style="list-style-type: none"> • Tobacco • Tomato • Spinach • <i>Arabidopsis</i> 	<ul style="list-style-type: none"> • Various strains of <i>Saccharomyces cerevisia</i> [Y187, Y190, W303(a), S288C, SP1, Σ1278b, BJ2168, cdc25] • <i>Schizosaccharomyces pombe</i> 	<ul style="list-style-type: none"> • <i>Saccharomyces cerevisiae</i> • <i>Pichia pastoris</i> • <i>Schizosaccharomyces pombe</i> 	Strain/Cell line compatibility
1 ml of reagent is sufficient for extraction of 0.5-1 g of plant leaves	3 ml of 1x nuclei isolation buffer per gram of leaf tissue (typically use 60 ml per isolation for 20 g leaf tissue)	2.5-5 ml per g of yeast cells	300 µl of reaction buffer and 500 µl of extraction buffer per 2 x 10 ⁹ cells	Usage
Non-ionic/No – only partially	Non-ionic/No – only partially	Non-ionic/No – only partially	Non-ionic/No – only partially	Type of detergent/ Dialyzable?
Yes – Suitable for protein activity assays	N/A; presence of DTT in protein extraction buffer may interfere with protein activity	Yes – Suitable for protein activity assays	Yes – Suitable for protein activity assays	Protein active?
<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails for plant tissue 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels – must be done in absence of DTT • FLAG (no DTT) • HIS-Select/IMAC • GST • Protease and Phosphatase Inhibitor Cocktails for plant tissue 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels – must be done in absence of DTT • FLAG (no DTT) • HIS-Select/IMAC • GST • Protease Inhibitor Cocktails 	<ul style="list-style-type: none"> • Antibody/EZview Affinity Gels • FLAG • HIS-Select/IMAC • GST • Protease Inhibitor Cocktails 	Compatible purification and detection products
<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining 	<ul style="list-style-type: none"> • Protein assay • Protein staining 	Quantitation and staining capabilities
<ul style="list-style-type: none"> • Western blotting • DNA-protein interaction assays (Gel-shift) • Enzymatic assays (e.g., kinase assays and phosphatase assays) 	<ul style="list-style-type: none"> • Western blotting • DNase I footprinting • Gel-shift assays • Source of chromatin, genomic DNA and RNA 	<ul style="list-style-type: none"> • Reporter gene assays • Western blotting • Immunoprecipitation • Affinity purification • Gel-shift assays 	<ul style="list-style-type: none"> • Reporter gene assays • Western blotting • Immunoprecipitation • Affinity purification • Gel-shift assays 	Applications
N/A	N/A	FLAG Y purification kit (CELLYM2), HIS-Select Y Purification Kit (HYP1)	N/A	Related products containing this reagent
Ready-to-use solution	Solutions ready to mix	Ready-to-use solution	Solutions ready to mix	Physical form
Room temperature 2 years	2-8 °C; Filter Mesh 100 is stored at room temperature 2 years	Room temperature 2 years	-20 °C; see bulletin for storage of individual components once opened 2 years	Storage and stability

Argentina

SIGMA-ALDRICH DE ARGENTINA S.A.
Free Tel: 0810 888 7446
Tel: (+54) 11 4556 1472
Fax: (+54) 11 4552 1698

Australia

SIGMA-ALDRICH PTY LTD.
Free Tel: 1800 800 097
Free Fax: 1800 800 096
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Austria

SIGMA-ALDRICH HANDELS GmbH
Tel: (+43) 1 605 81 10
Fax: (+43) 1 605 81 20

Belgium

SIGMA-ALDRICH NV/SA.
Free Tel: 0800 14747
Free Fax: 0800 14745
Tel: (+32) 3 899 13 01
Fax: (+32) 3 899 13 11

Brazil

SIGMA-ALDRICH BRASIL LTDA.
Free Tel: 0800 701 7425
Tel: (+55) 11 3732 3100
Fax: (+55) 11 5522 9895

Canada

SIGMA-ALDRICH CANADA LTD.
Free Tel: 1800 565 1400
Free Fax: 1800 265 3858
Tel: (+1) 905 829 9500
Fax: (+1) 905 829 9292

China

SIGMA-ALDRICH (SHANGHAI)
TRADING CO. LTD.
Free Tel: 800 819 3336
Tel: (+86) 21 6141 5566
Fax: (+86) 21 6141 5567

Czech Republic

SIGMA-ALDRICH S.R.O.
Tel: (+420) 246 003 200
Fax: (+420) 246 003 291

Denmark

SIGMA-ALDRICH DENMARK A/S
Tel: (+45) 43 56 59 10
Fax: (+45) 43 56 59 05

Finland

SIGMA-ALDRICH FINLAND OY
Tel: (+358) 9 350 9250
Fax: (+358) 9 350 9255

France

SIGMA-ALDRICH CHIMIE S.à.r.l.
Free Tel: 0800 211 408
Free Fax: 0800 031 052
Tel: (+33) 474 82 28 00
Fax: (+33) 474 95 68 08

Germany

SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800 51 55 000
Free Fax: 0800 64 90 000
Tel: (+49) 89 6513 0
Fax: (+49) 89 6513 1160

Greece

SIGMA-ALDRICH (O.M.) LTD.
Tel: (+30) 210 994 8010
Fax: (+30) 210 994 3831

Hungary

SIGMA-ALDRICH Kft
Ingyenes zöld telefon: 06 80 355 355
Ingyenes zöld fax: 06 80 344 344
Tel: (+36) 1 235 9055
Fax: (+36) 1 235 9050

India

SIGMA-ALDRICH CHEMICALS
PRIVATE LIMITED
Telephone
Bangalore: (+91) 80 6621 9600
New Delhi: (+91) 11 4165 4255
Mumbai: (+91) 22 2570 2364
Hyderabad: (+91) 40 6684 5488
Fax
Bangalore: (+91) 80 6621 9650
New Delhi: (+91) 11 4165 4266
Mumbai: (+91) 22 2579 7589
Hyderabad: (+91) 40 6684 5466

Ireland

SIGMA-ALDRICH IRELAND LTD.
Free Tel: 1800 200 888
Free Fax: 1800 600 222
Tel: (+353) 1 404 1900
Fax: (+353) 1 404 1910

Israel

SIGMA-ALDRICH ISRAEL LTD.
Free Tel: 1 800 70 2222
Tel: (+972) 8 948 4100
Fax: (+972) 8 948 4200

Italy

SIGMA-ALDRICH S.r.l.
Numero Verde: 800 827018
Tel: (+39) 02 3341 7310
Fax: (+39) 02 3801 0737

Japan

SIGMA-ALDRICH JAPAN K.K.
Tokyo Tel: (+81) 3 5796 7300
Tokyo Fax: (+81) 3 5796 7315

Korea

SIGMA-ALDRICH KOREA
Free Tel: (+82) 80 023 7111
Free Fax: (+82) 80 023 8111
Tel: (+82) 31 329 9000
Fax: (+82) 31 329 9090

Malaysia

SIGMA-ALDRICH (M) SDN. BHD
Tel: (+60) 3 5635 3321
Fax: (+60) 3 5635 4116

Mexico

SIGMA-ALDRICH QUÍMICA, S.A. de C.V.
Free Tel: 01 800 007 5300
Free Fax: 01 800 712 9920
Tel: 52 722 276 1600
Fax: 52 722 276 1601

The Netherlands

SIGMA-ALDRICH CHEMIE BV
Free Tel: 0800 022 9088
Free Fax: 0800 022 9089
Tel: (+31) 78 620 5411
Fax: (+31) 78 620 5421

New Zealand

SIGMA-ALDRICH NEW ZEALAND LTD.
Free Tel: 0800 936 666
Free Fax: 0800 937 777
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Norway

SIGMA-ALDRICH NORWAY AS
Tel: (+47) 23 17 60 60
Fax: (+47) 23 17 60 50

Poland

SIGMA-ALDRICH Sp. z o.o.
Tel: (+48) 61 829 01 00
Fax: (+48) 61 829 01 20

Portugal

SIGMA-ALDRICH QUÍMICA, S.A.
Free Tel: 800 202 180
Free Fax: 800 202 178
Tel: (+351) 21 924 2555
Fax: (+351) 21 924 2610

Russia

SIGMA-ALDRICH RUS, LLC
Tel: +7 (495) 621 6037
Fax: +7 (495) 621 5923

Singapore

SIGMA-ALDRICH PTE. LTD.
Tel: (+65) 6779 1200
Fax: (+65) 6779 1822

South Africa

SIGMA-ALDRICH
SOUTH AFRICA (PTY) LTD.
Free Tel: 0800 1100 75
Free Fax: 0800 1100 79
Tel: (+27) 11 979 1188
Fax: (+27) 11 979 1119

Spain

SIGMA-ALDRICH QUÍMICA, S.A.
Free Tel: 900 101 376
Free Fax: 900 102 028
Tel: (+34) 91 661 99 77
Fax: (+34) 91 661 96 42

Sweden

SIGMA-ALDRICH SWEDEN AB
Tel: (+46) 8 742 4200
Fax: (+46) 8 742 4243

Switzerland

SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800 80 00 80
Free Fax: 0800 80 00 81
Tel: (+41) 81 755 2828
Fax: (+41) 81 755 2815

United Kingdom

SIGMA-ALDRICH COMPANY LTD.
Free Tel: 0800 717 181
Free Fax: 0800 378 785
Tel: (+44) 1747 833 000
Fax: (+44) 1747 833 313
SAFC (UK) Free Tel: 0800 71 71 17

United States

SIGMA-ALDRICH
P.O. Box 14508
St. Louis, Missouri 63178
Toll-Free: 800 325 3010
Toll-Free Fax: 800 325 5052
Call Collect: (+1) 314 771 5750
Tel: (+1) 314 771 5765
Fax: (+1) 314 771 5757

Internet

sigma-aldrich.com

World Headquarters

3050 Spruce St., St. Louis, MO 63103
(314) 771-5765
sigma-aldrich.com

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052

Technical Service (800) 325-5832 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries SAFC™ (800) 244-1173

The SIGMA-ALDRICH Group

SIGMA-ALDRICH

Accelerating Customers' Success
through Leadership in Life Science,
High Technology and Service

©2007 Sigma-Aldrich Co. All rights reserved.

SIGMA, SAFC, SAFC, SIGMA-ALDRICH, ISOTECH, ALDRICH, FLUKA, and SUPELCO are trademarks belonging to Sigma-Aldrich Co. and its affiliate Sigma-Aldrich Biotechnology, L.P. Riedel-de Haën® is a trademark under license from Riedel-de Haën GmbH. Sigma brand products are sold through Sigma-Aldrich, Inc. CellLytic™, EZView™, HIS-Select®, ILAP®, FLAG® and NuCLEAR™ are registered trademarks of Sigma-Aldrich Biotechnology. DH5α™ is a registered trademark of Invitrogen. IGEPAL® is a registered trademark of Rhone-Poulenc AG Co. Triton® is a registered trademark of Union Carbide Corp. Benzozase® is a registered trademark of Merck KGaA, Darmstadt, Germany. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.

JFQ
01938-504230
0017