



643PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Mammalian GenLysates™

Lyophilized Protein Lysates

(Cat. # See Appendix)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION 3

ITEM(S) SUPPLIED 3

STORAGE & STABILITY 3

PROTOCOL 3

CERTIFICATE OF ANALYSIS 3

PRODUCT WARRANTY: 4

APPENDIX: MAMMALIAN GENLYSATES™ 4

 BRAIN LYSATES 4

 BREAST LYSATES 4

 CERVIX LYSATES 4

 COLON MUSCLE LYSATES 4

 ESOPHAGUS LYSATE 4

 EYE LYSATES 4

 HEART LYSATES 5

 HYPOPHARYNX LYSATES 5

 KIDNEY LYSATES 5

 LIVER LYSATES 5

 LUNG LYSATES 5

 OVARY LYSATES 5

 PANCREAS LYSATES 5

 PROSTATE LYSATES 5

 RECTUM LYSATES 5

 SKELETAL MUSCLE LYSATES 5

 SKIN LYSATES 5

 SPINAL CORD LYSATES 5

 SPLEEN LYSATES 5

 STOMACH LYSATES 6

 TESTIS LYSATES 6

 THYROID LYSATES 6

 UTERUS LYSATES 6

RELATED PRODUCTS 6

INTRODUCTION

GenLysates™ are ready-to-use total protein extracts from a wide selection of tissues and cells. The GenLysates™ are ideal for SDS-PAGE and Western blot analysis. The total proteins extraction was performed according to Laemmli¹. The proteins are extracted in a denaturing buffer containing 62mM Tris.HCl and 2% SDS with a protease inhibitor cocktail to minimize proteolytic damage to proteins.

ITEM(S) SUPPLIED

Description	Size
GenLysate™	150µg protein

See Appendix for catalog number and product description.

STORAGE & STABILITY

GenLysates™ are shipped on blue ice and should be stored at -20°C upon arrival. Stable for 1 year if stored as recommended.

NOTE: After reconstitution, the GenLysate™ must be stored at -70°C.

PROTOCOL

1. Briefly, spin the tube containing GenLysate™ and reconstitute in 30µl deionized water to give 5µg/µl concentration.
2. Add an appropriate volume of sample buffer containing a reducing agent (e.g. 5% β-mercaptoethanol or 5-20mM DTT final concentration) in the reconstituted GenLysate™ tube.
3. The GenLysate™ may also be diluted if needed with an appropriate sample buffer containing a reducing agent.
4. Before loading the sample on the gel, put the tube in a boiling water bath for approx 5 minutes and briefly spin it again. Vortex the tube and the sample is now ready for loading on to the gel.

NOTE: Avoid frequent thawing and freezing of the GenLysate™. It is recommended to make appropriate aliquots of lysate after reconstitution and then freeze them at -70°C.

CERTIFICATE OF ANALYSIS

- The protein concentration of GenLysate™ samples was measured using NI™-Protein Assay (G-Biosciences' Cat. # 786-005) with BSA as a standard.
- The lysate was run on a 4-20% polyacrylamide gel and after electrophoresis the gel was stained with RapidStain™ (G-Biosciences' Cat. # 786-31). The protein bands were sharp and well resolved and there was no smear. Proteins from 5-200kD were visible.
- When appropriate, the protein lysates are also tested by Western blot analysis; protein transfer blots are probed for the presence of actin or tubulin.

PRODUCT WARRANTY:

GenLysates™ are warranted to perform at least once using kits and reagents from G-Biosciences.

APPENDIX: MAMMALIAN GENLYSATES™

Cat. #	Region	Species	Tissue Type
Brain Lysates			
NLH-02	Whole	Human	Normal
SLH-026	Cerebellum	Human	Normal
SLH-025	Frontal Cortex	Human	Normal
SLH-024	Hippocampus	Human	Normal
SLH-023	Medulla	Human	Normal
SLH-0210	Mid Brain	Human	Normal
SLH-029	Olfactory Lobe	Human	Normal
SLH-021	Pons	Human	Normal
SLH-0211	Posterior Cortex	Human	Normal
SLH-028	Striatum	Human	Normal
SLH-0213	Temporal Lobe	Human	Normal
SLH-0212	Thalamus	Human	Normal
NLM-02	Whole	Mouse	Normal
SLM-026	Cerebellum	Mouse	Normal
SLM-022	Entorhinal Cortex	Mouse	Normal
SLM-025	Frontal Cortex	Mouse	Normal
SLM-024	Hippocampus	Mouse	Normal
SLM-023	Medulla	Mouse	Normal
SLM-0210	Mid Brain	Mouse	Normal
SLM-029	Olfactory Lobe	Mouse	Normal
SLM-021	Pons	Mouse	Normal
SLM-0211	Posterior Cortex	Mouse	Normal
SLM-028	Striatum	Mouse	Normal
SLM-0212	Thalamus	Mouse	Normal
PBL-09	Amygdala	Macaque	Normal
PBL-14	Cingulate Gyrus	Macaque	Normal
PBL-08	Cerebellum	Macaque	Normal
PBL-13	Frontal Cortex	Macaque	Normal
PBL-04	Hippocampus	Macaque	Normal
PBL-06	Hypothalamus	Macaque	Normal
PBL-10	Medulla	Macaque	Normal
PBL-11	Occipital Lobe	Macaque	Normal
PBL-12	Parahippocampal Gyrus	Macaque	Normal
PBL-15	Parietal Lobe	Macaque	Normal
PBL-16	Pituitary	Macaque	Normal

Cat. #	Region	Species	Tissue Type
PBL-05	Pons	Macaque	Normal
PBL-03	Striatum	Macaque	Normal
PBL-07	Temporal Lobe	Macaque	Normal
PBL-02	Thalamus	Macaque	Normal
NLR-02	Whole	Rat	Normal
SLR-026	Cerebellum	Rat	Normal
SLR-022	Entorhinal Cortex	Rat	Normal
SLR-025	Frontal Cortex	Rat	Normal
SLR-024	Hippocampus	Rat	Normal
SLR-023	Medulla	Rat	Normal
SLR-0210	Mid Brain	Rat	Normal
SLR-029	Olfactory Lobe	Rat	Normal
SLR-021	Pons	Rat	Normal
SLR-0211	Posterior Cortex	Rat	Normal
SLR-028	Striatum	Rat	Normal
SLR-0212	Thalamus	Rat	Normal
Breast Lysates			
NLH-13	Whole	Human	Normal
TLH-13	Whole	Human	Tumor
Cervix Lysates			
NLH-17	Whole	Human	Normal
TLH-17	Whole	Human	Tumor
Colon Muscle Lysates			
NLH-20	Whole	Human	Normal
TLH-20	Whole	Human	Tumor
Esophagus Lysate			
NLH-22	Whole	Human	Normal
Eye Lysates			
NLB-10	Whole	Rabbit	Normal
SLB-1003	Aqueous Humor	Rabbit	Normal
SLB-1002	Cornea	Rabbit	Normal
SLB-1001	Retina	Rabbit	Normal
NLR-10	Whole	Rat	Normal
SLR-1003	Aqueous Humor	Rat	Normal
SLR-1002	Cornea	Rat	Normal
SLR-1001	Retina	Rat	Normal

Cat. #	Region	Species	Tissue Type
Heart Lysates			
NLH-08	Whole	Human	Normal
SLH-087	Aorta	Human	Normal
SLH-088	Atrium (Left)	Human	Normal
SLH-089	Atrium (Right)	Human	Normal
SLH-090	Vena Cava	Human	Normal
SLH-081	Ventricle (Left)	Human	Normal
SLH-082	Ventricle (Right)	Human	Normal
NLM-08	Whole	Mouse	Normal
ML-10	Whole	Pig	Normal
SLP-087	Aorta	Pig	Normal
SLP-085	Aortic Valve	Pig	Normal
SLP-083	Mitral Valve	Pig	Normal
SLP-084	Pulmonary Valve	Pig	Normal
SLP-086	Tricuspid Valve	Pig	Normal
SLP-081	Ventricle (Left)	Pig	Normal
SLP-082	Ventricle (Right)	Pig	Normal
NLB-08	Whole	Rabbit	Normal
NLR-08	Whole	Rat	Normal
Hypopharynx Lysates			
NLH-15	Whole	Human	Normal
TLH-15	Whole	Human	Tumor
Kidney Lysates			
NLH-04	Whole	Human	Normal
NLM-04	Whole	Mouse	Normal
SLM-042	Cortex	Mouse	Normal
SLM-041	Medulla	Mouse	Normal
NLB-04	Whole	Rabbit	Normal
NLR-04	Whole	Rat	Normal
SLR-042	Cortex	Rat	Normal
SLR-041	Medulla	Rat	Normal
Liver Lysates			
NLH-01	Whole	Human	Normal
TLH-01	Whole	Human	Tumor
ML-05	Whole	Macaque	Normal
NLM-01	Whole	Mouse	Normal
CFM-014	Smooth ER Fraction	Mouse	Normal
CFM-013	Rough ER Fraction	Mouse	Normal
CFM-012	Nuclear Fraction	Mouse	Normal
CFM-011	Mitochondria	Mouse	Normal
CFM-015	Golgi Complex	Mouse	Normal

Cat. #	Region	Species	Tissue Type
ML-07	Whole	Pig	Normal
NLB-01	Whole	Rabbit	Normal
NLR-01	Whole	Rat	Normal
Lung Lysates			
NLH-03	Whole	Human	Normal
TLH-03	Whole	Human	Tumor
NLM-03	Whole	Mouse	Normal
NLB-03	Whole	Rabbit	Normal
NLR-03	Whole	Rat	Normal
Ovary Lysates			
NLH-07	Whole	Human	Normal
TLH-07	Whole	Human	Tumor
NLM-07	Whole	Mouse	Normal
NLR-07	Whole	Rat	Normal
Pancreas Lysates			
NLH-09	Whole	Human	Normal
TLH-09	Whole	Human	Tumor
NLM-09	Whole	Mouse	Normal
NLR-09	Whole	Rat	Normal
Prostate Lysates			
NLH-14	Whole	Human	Normal
TLH-14	Whole	Human	Tumor
Rectum Lysates			
NLH-16	Whole	Human	Normal
TLH-16	Whole	Human	Tumor
Skeletal Muscle Lysates			
NLH-21	Whole	Human	Normal
NLB-21	Whole	Rabbit	Normal
Skin Lysates			
NLH-18	Whole	Human	Normal
TLH-18	Whole	Human	Tumor
Spinal Cord Lysates			
SLH-027	Whole	Human	Normal
SLM-027	Whole	Mouse	Normal
PBL-01	Whole	Macaque	Normal
SLR-027	Whole	Rat	Normal
Spleen Lysates			
NLH-05	Whole	Human	Normal
NLM-05	Whole	Mouse	Normal
NLR-05	Whole	Rat	Normal

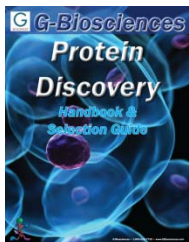
Cat. #	Region	Species	Tissue Type
<i>Stomach Lysates</i>			
NLH-12	Whole	Human	Normal
TLH-12	Whole	Human	Tumor
<i>Testis Lysates</i>			
NLH-06	Whole	Human	Normal
TLH-06	Whole	Human	Tumor
NLM-06	Whole	Mouse	Normal
NLR-06	Whole	Rat	Normal
<i>Thyroid Lysates</i>			
NLH-19	Whole	Human	Normal
TLH-19	Whole	Human	Tumor
<i>Uterus Lysates</i>			
NLH-11	Whole	Human	Normal
TLH-11	Whole	Human	Tumor

REFERENCE

1. Laemmli, U.K. (1970) Nature. 227: 680

RELATED PRODUCTS

Download our Protein Discovery Handbook.



<http://info.gbiosciences.com/complete-protein-discovery-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 8/10/2012 CMH

This page is intentionally left blank



www.GBiosciences.com