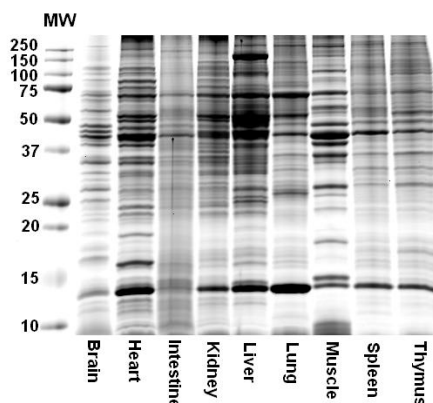




Product Data Sheet

Product Name:	Mouse Muscle Tissue Lysate (<i>Mus musculus</i>)
Catalog Number:	29508
Lot Number:	See label on the vial
Product Description:	Muscle from C57/BL6 mice was collected, homogenized in modified RIPA buffer, and cleared by centrifugation. Total protein concentration of the lysate was determined using BCA protein assay (BSA was used as a standard). Lysate was mixed with 2X Laemmli sample buffer containing reducing reagent and heated for 5 minutes at 90 °C. Lysate is supplied as 200 µg per vial at 2.0 mg/ml total protein concentration in 1 x Laemmli sample buffer.
RIPA Buffer:	150 mM NaCl, 50 mM Tris-Cl, 1% Triton X-100, 0.5% sodium deoxycholate, 0.1% sodium dodecyl sulfate, 1 mM EDTA, 1 mM PMSF, 5 µg/ml Pepstatin A, 2 µg/ml Leupeptin, 1 mM sodium orthovanadate, 1 mM sodium fluoride.
Application:	Mouse muscle lysate is suitable for SDS-PAGE electrophoresis and/or consecutive Western immunoblotting procedure. 5-20 µg of the total protein per lane is recommended for mini gels.



SDS-PAGE image of mouse tissue lysates (10 µg protein/lane). Proteins were separated on 8-16% Tris-HCl gradient gel.

Storage: Store at -80 °C for up to 6 months; avoid repeating freeze-thaw cycles.

Related Products:

Mouse Brain Tissue Lysate Cat. #29507
Mouse Heart Tissue Lysate Cat. #29511
Mouse Intestine Tissue Lysate Cat. #29513
Mouse Kidney Tissue Lysate Cat. #55891
Mouse Liver Tissue Lysate Cat. #29509
Mouse Lung Tissue Lysate Cat. #29510
Mouse Spleen Tissue Lysate Cat. #29505
Mouse Thymus Tissue Lysate Cat. #55890

This product is for in vitro research use only.