

Instruction Manual

Catalog Number	PK-AB718-1477
Description	Adipose tissue lysate was prepared by homogenization in lysis buffer (10 mM HEPES pH 7.9, 1.5 mM MgCl ₂ , 10 mM KCl, 1 mM ethylenediaminetetraacetic acid, 10% glycerol, 1% NP-40, and a cocktail of protease inhibitors). Tissue and cell debris was removed by centrifugation. Protein concentration was determined with Bio-Rad protein assay. The product was boiled for 5 min in 1 x SDS sample buffer (50 mM Tris-HCl pH 6.8, 12.5% glycerol, 1% sodium dodecylsulfate, 0.01% bromophenol blue) containing 5% β-mercaptoethanol.
Quantity	100 µg
Cell Line / Source	Rat adipose tissue from outbred Sprague Dawley Albino rat
Species	Rat
Buffer	Product is supplied in 1x SDS sample buffer containing 5% β-mercaptoethanol.
Growth Media	NA
Formulation	Product is supplied at a concentration of 2 mg/ml. During shipment, small volumes will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Store at 4°C for continuous use. Store at -20°C for three months or at -70°C for up to one year. Avoid repeated freezing and thawing (store working aliquots at -70°C). Under proper storage conditions the shelf life is 6-12 months from the date of receipt.
Tested Application	Western Blot

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.