

R1000.RC Lot No. 1310186

Sprague Dawley (SD) Rat Kidney Cytosol

Untreated, Male, Pool of 150

1.0 mL at 5 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 2 mM EDTA

Animal Information

Species:	Rat
Strain:	International Genetic Standard (IGS), Sprague Dawley
Sex:	Male
Age:	~8 weeks
Vendor:	Charles River, Raleigh, NC

Animals were housed in an AAALAC-accredited facility and allowed to acclimate \geq seven days before use.

Food:	Purina 5L79 (<i>ad libitum</i>)
Water:	Automatic watering system (<i>ad libitum</i>)
Light/dark cycle:	5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)
Temperature:	70°F \pm 2°F
Humidity:	30-70 %
Cage:	Beta Chip (hardwood), NEPCO, Warrensburg, NY

Subcellular fractions were prepared from whole kidney.



Store at -80°C

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for *in vitro* use only.

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 19 August 2013