

## R1000.IS9 Lot No. 2310231

Sprague Dawley (SD) Rat Intestine S9 Fraction

Untreated, Male, Pool of 200

1.0 mL at 4 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 1 mM EDTA, 20% glycerol, heparin, PMSF, leupeptin, DTT, aprotinin

Enzyme Activities		Rate
NADPH-cytochrome c reductase	(nmol/mg protein/min)	30.3 ± 1.2
Testosterone 6β-hydroxylation	(pmol/mg protein/min)	52.1 ± 4.1
Midazolam 1'-hydroxylation	(pmol/mg protein/min)	32.0 ± 3.0
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	59.7 ± 7.8

Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

Aprotinin, Leupeptin, and Phenylmethylsulfonyl-fluoride were used in the preparation of this S9 fraction. Subcellular fractions were prepared from duodenal and jejunal tissue.

To measure cytochrome P450 (CYP) activity, intestine S9 samples (0.2 mg/mL) were incubated in triplicate at 37 ± 2°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and testosterone (250 μM) or midazolam (30 μM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure UDP-glucuronosyltransferase (UGT) activity, intestine S9 samples (0.2 mg/mL) were incubated in triplicate at 37 ± 2°C for 10 minutes in Tris-HCl (100 mM, pH 7.7 at 37°C), CHAPS (0.5 mM), EDTA (1.0 mM), MgCl<sub>2</sub> (10 mM), D-saccharic acid 1,4-lactone (100 μM), uridine diphosphate-glucuronic acid (10.0 mM) and 4-methylumbelliferone (1 mM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

### 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 ≥ seven days before use.

Food: Purina 5L79 (*ad libitum*)  
 Water: Automatic watering system (*ad libitum*)  
 Light/dark cycle: 5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)  
 Temperature: 70°F ± 2°F  
 Humidity: 30-70 %  
 Bedding: Beta Chip (hardwood), NEPCO, Warrensburg, NY  
 Cage: Polycarbonate Shoebox Cage, conventional cage



## 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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**

Signature and Date: Stephanie Helmstetter 20 September 2023