

## R1000.S9 Lot No. 2210317

Sprague Dawley (SD) Rat Liver S9 Fraction Untreated, Male, Pool of 1061
1.0 mL at 20 mg protein / mL

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

Specific Content and Enzyme Activities		Content / Rate
Cytochrome P450 content	(nmol/mg protein)	0.231
Cytochrome b₅ content	(nmol/mg protein)	0.175
7-Ethoxycoumarin <i>O</i> -dealkylation	(pmol/mg protein/min)	409 ± 25
Glucuronidation of 4-methylumbelliferone	(nmol/mg protein/min)	90.9 ± 7.7
CDNB <sup>a</sup>	(nmol/mg protein/min)	579 ± 14

<sup>&</sup>lt;sup>a</sup> 1-Chloro-2,4-dinitrobenzene-glutathione conjugation by glutathione S-transferase.

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

To measure cytochrome P450 (CYP) activity, liver S9 samples (0.2 mg/mL) were incubated in triplicate at  $37 \pm 2^{\circ}$ 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 7-ethoxycoumarin (500  $\mu$ M), 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, liver S9 samples (0.1 mg/mL) were incubated in triplicate at  $37 \pm 2^{\circ}$ C for 10 minutes in Tris-HCl (100 mM, pH 7.7 at  $37^{\circ}$ C), CHAPS (0.5 mM), EDTA (1.0 mM), MgCl<sub>2</sub> (10 mM), D-saccharic acid 1,4-lactone (100  $\mu$ M), uridine diphosphate-glucuronic acid (8.0 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.

To measure glutathione S-transferase activity (GST), liver S9 samples (5 to  $50 \mu g/mL$ ) were incubated in triplicate at  $37 \pm 2^{\circ}C$  for 10 minutes in potassium phosphate buffer (100 mM, pH 6.5), glutathione (1 mM), and CDNB (1 mM), at the final concentrations indicated. Reaction rates are determined by photometric kinetic measurements at 340 nm.

## **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:  $70^{\circ}\text{F} \pm 2^{\circ}\text{F}$ Humidity:  $30-70^{\circ}$ 

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.