

## R1000.ES9 Lot No. 2110123

Sprague Dawley Rat Skin S9 Fraction

Untreated, Male, Pool of 10

1.0 mL at 5 mg protein / mL

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

<b>Enzyme Activities</b>		<b>Rate</b>
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	4.03 ± 0.16
Testosterone 6β-hydroxylation	(pmol/mg protein/min)	7.03 ± 0.40
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	4.19 ± 0.49
Clopidogrel hydrolysis	(pmol/mg protein/min)	141 ± 10
6α-Methylprednisolone 21-hemisuccinate hydrolysis	(pmol/mg protein/min)	658 ± 52

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, skin 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), 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, skin 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 (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 carboxylesterase activity, skin S9 samples (0.15 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), and clopidogrel hydrogen sulfate (150 μM), at the final concentrations indicated. Metabolite formation was determined by LC-MS/MS methods with deuterated metabolites as internal standards.

To measure carboxylesterase activity, skin S9 samples (0.15 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), and 6α-methylprednisolone 21-hemisuccinate (750 μM), at the final concentrations indicated. Metabolite formation was determined by 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 ( <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 ± 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.

Datasheet prepared 21 May 2021