

## L1000 Lot No. 1610101

New Zealand Rabbit Liver Microsomes  
 Untreated, Male, Pool of 8  
 0.5 mL at 20 mg protein / mL  
 Suspension medium: 250 mM sucrose

Specific Content and Enzyme Activities		Content / Rate
Cytochrome P450 content	(nmol/mg protein)	2.214
Cytochrome b <sub>5</sub> content	(nmol/mg protein)	1.730
NADPH-cytochrome c reductase	(nmol/mg protein/min)	194 ± 4
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	3160 ± 140

Characterization is performed when the first lot of a product from a given subcellular fraction (e.g., S9) is prepared. Subsequent lots are subject to a verification test only. 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 microsomes (50 µg/mL) were incubated in triplicate at 37 ± 1°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 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

### Animal Information

Species: Rabbit  
 Strain: New Zealand  
 Sex: Male  
 Age: >9 months  
 Vendor: Covance Research Products, Denver, PA

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

Food: Purina 5326 High Fiber Diet (*ad libitum*)  
 Water: Automatic watering system (*ad libitum*)  
 Light/dark cycle: 16 hours light / 8 hours dark  
 Temperature: 68°F -72°F  
 Humidity: 50-55 %  
 Bedding: None used  
 Cage: Conventional wire grid 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 13 September 2016