

L1500.H15 Lot No. 1210410

Cryopreserved New Zealand Rabbit Hepatocytes
 Female, Pool of 3

Assured Minimum Yield: 5.0×10^6 cells per vial
 Viability: 74.2%

Livers were perfused and subjected to collagenase digestion for the purpose of hepatocyte isolation.

Enzyme Activities		Rate
7-Ethoxycoumarin O-dealkylation	(pmol/million cells/min)	1150 ± 60
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	217 ± 21
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	23.9 ± 2.7

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

To measure metabolic enzyme activities, hepatocytes (1×10^6 /mL) in suspension were incubated in triplicate at $37 \pm 1^\circ\text{C}$ for 30 minutes in Waymouth's medium and 7-ethoxycoumarin (500 μM). Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

Animal Information

Species: Rabbit
 Strain: New Zealand
 Sex: Female
 Age: Sexually mature
 Vendor: Covance Research Products, Denver, PA

Animals were housed in an AAALAC-accredited facility and allowed to acclimate \geq 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: $70^\circ\text{F} \pm 2^\circ\text{F}$
 Humidity: 30-70 %
 Bedding: None used
 Cage: Conventional wire grid cage



Store in liquid nitrogen, vapor phase

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 02 November 2012