

## Z6000.H15 Lot No. 1410242

Cryopreserved Gottingen Minipig Hepatocytes  
 Male, Pool of 3

Assured Minimum Yield:  $4.0 \times 10^6$  cells per vial  
 Viability: 78%

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

<b>Enzyme Activities</b>		<b>Rate</b>
7-Ethoxycoumarin O-dealkylation	(pmol/million cells/min)	127 ± 13
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	27.6 ± 2.0
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	No peak detected

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 ( $1 \times 10^6$  /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: Minipig  
 Strain: Gottingen  
 Sex: Male  
 Age: Sexually mature  
 Vendor: Marshall Bioresources, New York

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

Food: Harlan Swine Feed  
 Water: Automatic watering system (*ad libitum*)  
 Light/dark cycle: 12 hour light cycle  
 Temperature: 61-81°F  
 Humidity: 30-70 %



## 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 10 December 2014