

R6000.H15+ Lot No. 1610230

Cryopreserved Wistar Han Rat Hepatocytes
 Male, Pool of 8

Assured Minimum Yield: 7.0×10^6 cells per vial
 Viability: 78%

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

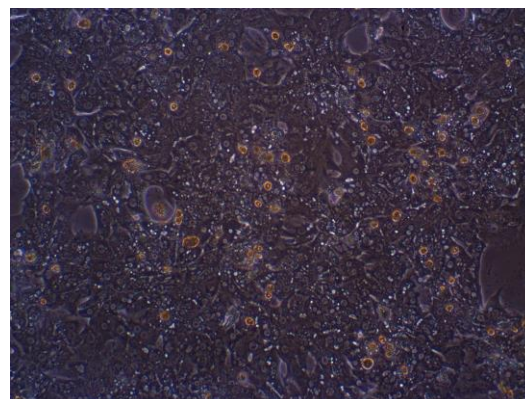
Enzyme Activities		Rate
7-Ethoxycoumarin O-dealkylation	(pmol/million cells/min)	257 ± 31
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	212 ± 15
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	114 ± 5

Values for enzyme activities were determined at a single substrate concentration and are mean \pm 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 Opti^{INCUBATE} 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: Rat
 Strain: Wistar Han
 Sex: Male
 Age: ~ 8-12 weeks
 Vendor: Charles River, Raleigh, NC
 Animals were housed in an AAALAC-accredited facility and allowed to acclimate \geq 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 %
 Bedding: Beta Chip (hardwood), NEPCO, Warrensburg, NY
 Cage: Polycarbonate Shoebox Cage, conventional cage



R6000.H15+ 1610230 day 3 of culture

Plate format	Recommended Seeding Density (million cells/mL)	Recommended Seeding/Feeding Volume Per Well
6 well format	1.2	1.7 mL
12 well format	1.2	650 μL
24 well format	1.2	330 μL
48 well format	1.2	150 μL
96 well format	Not Tested	



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 25 July 2016