

R1000.H15+ Lot No. 1510038

Cryopreserved Sprague Dawley (SD) Rat Hepatocytes
Male, Pool of 12

Assured Minimum Yield: 7.0×10^6 cells per vial
Viability: 82%

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)	249± 15
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	192 ± 4
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	99.5 ± 5.0

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 William's E+ 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: International Genetic Standard (IGS), Sprague Dawley
Sex: Male
Age: ~ 8-12 weeks
Vendor: Charles River, Raleigh, NC

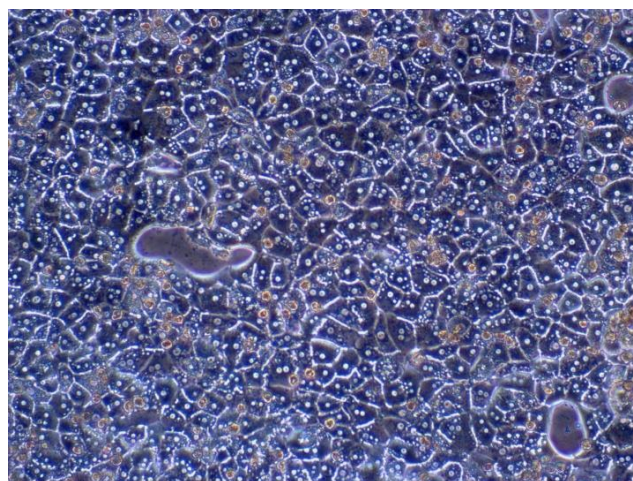
Animals were housed in an AAALAC-accredited facility and allowed to acclimate \geq seven days before use.

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

Recommended seeding density
1.4 million cells/mL



R1000.H15+ 1510038 day 3 of culture



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 05 March 2015