

## M1000.H15+ Lot No. 1710136

Cryopreserved CD1 Mouse Hepatocytes  
 Male, Pool of 15

Assured Minimum Yield:  $2.0 \times 10^6$  cells per vial  
 Viability: 96%

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) | 162 ± 12 |
| 7-Hydroxycoumarin glucuronidation | (pmol/million cells/min) | 348 ± 13 |
| 7-Hydroxycoumarin sulfonation     | (pmol/million cells/min) | 119 ± 4  |

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 Opti<sup>INCUBATE</sup> medium and 7-ethoxycoumarin (500 µM). Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

### Animal Information

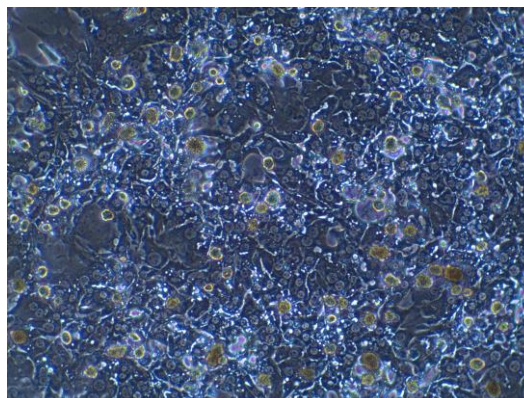
Species: Mouse  
 Strain: CD1  
 Sex: Male  
 Age: ~ 8-12 weeks  
 Vendor: Charles River, Raleigh, NC

Animals were housed in an AAALAC-accredited facility and allowed to acclimate ≥ 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



M1000.H15+ 1710136 day 3 of culture

|                | Recommended Seeding Density (million cells/mL) | Recommended Seeding/Feeding Volume Per Well |
|----------------|--|---|
| 6 well format  | 0.6  | 1.7 mL                                      |
| 12 well format | 0.6  | 650 µL                                      |
| 24 well format | 0.6  | 330 µL                                      |
| 48 well format | 0.6  | 150 µL                                      |
| 96 well format | 0.6  | 50 µL                                       |



## 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 18 August 2017