

**R3000.H15      Lot No. 1110064**

Cryopreserved Wistar Rat Hepatocytes

**Animal Information**

Tissue was obtained from male Wistar rats (n=4), between 8 and 12 weeks of age. (Charles River Laboratories, Raleigh, NC)

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

**Assured Minimum Yield:**  $7.0 \times 10^6$  per vial**Viability:** 75.6%

(Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and Hepatocyte Isolation Kit.)

**Marker substrate reaction (pmol/million cells/min)**

7-Ethoxycoumarin <i>O</i> -dealkylation	197 ± 17
7-Hydroxycoumarin glucuronidation	237 ± 22
7-Hydroxycoumarin sulfonation	112 ± 4

**CAUTION:** Although strict measures are taken to ensure that livers obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

**Store vials in liquid nitrogen, vapor phase.**

Data sheet prepared 23-Mar-2011