

**P2500.H15A**      **Lot 0610395**

Cryopreserved Cynomolgus Monkey Hepatocytes

**Animal Information**

Tissue was obtained from female cynomolgus monkeys (n=2), approximately 3 and 4 years of age. (Charles Rivers Laboratories, Sierra Division, Sparks, NV)

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

**Assured Minimum:** 4.0 x 10<sup>6</sup> per vial**Viability:** 84.3%

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

<b>Enzyme</b>	<b>Marker substrate reaction (pmol/cells per inc/min)</b>	
Phase I	7-Ethoxycoumarin <i>O</i> -Dealkylation	139
Phase II	7-Hydroxycoumarin Glucuronidation	96.3
Phase II	7-Hydroxycoumarin Sulfation	72.1

**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 1/16/07