

P2500.H15A Lot No. 1210031

Cryopreserved Cynomolgus Monkey Hepatocytes

Animal Information

Tissue was obtained from female Cynomolgus monkeys (n=3), 2-4 years of age. (Sierra Biomedical, Sparks, NV)

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

Assured Minimum Yield: 4.0×10^6 per vial**Viability:** 86.2%

(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 O-dealkylation	487 ± 42
7-Hydroxycoumarin glucuronidation	390 ± 5
7-Hydroxycoumarin sulfonation	92.1 ± 3.1

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 9 Mar 2012