

M1500.H15 Lot No. 0810436

Cryopreserved CD-1 Mouse Hepatocytes

Animal Information

Tissue was obtained from female CD-1 mice (n=8), between 8-12 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: 2.0×10^6 per vial**Viability:** 85.8%

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

Marker substrate reaction (pmol/cells per inc/min)

7-Ethoxycoumarin <i>O</i> -dealkylation	228 ± 5
7-Hydroxycoumarin glucuronidation	146 ± 0
7-Hydroxycoumarin sulfonation	335 ± 6

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 2/4/09