

L1000.H15 Lot No. 1110135

Cryopreserved New Zealand White Rabbit Hepatocytes

Animal Information

Tissue was obtained from male New Zealand white rabbits (n=3), approximately 6 months of age.
(Covance Research Products Inc., Denver, PA)

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

Assured Minimum Yield: 5.0×10^6 per vial**Viability:** 83.0%

(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	952 ± 72
7-Hydroxycoumarin glucuronidation	156 ± 12
7-Hydroxycoumarin sulfonation	46.6 ± 2.9

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 4-May-2011