

R1000.H15+ Lot No. 1110457

Cryopreserved Sprague Dawley Rat Hepatocytes

Animal InformationTissue was obtained from male Sprague Dawley rats (n=4), between 8 and 12 weeks of age.
(Charles Rivers Laboratories, Raleigh, NC)

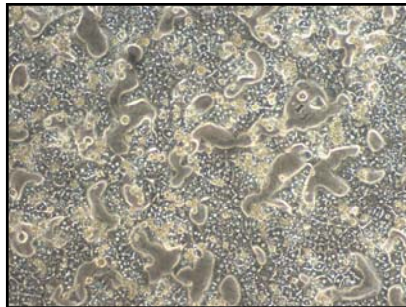
Livers were perfused and subjected to collagenase digestion for the purpose of hepatocyte isolation.

Assured Minimum Yield: 7.0×10^6 per vial**Viability:** 74.8%

(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	390 ± 57
7-Hydroxycoumarin glucuronidation	295 ± 37
7-Hydroxycoumarin sulfonation	181 ± 14



Lot 1110457 Day 3 of culture (1.2 million cells/mL)

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 27-Jan-2012