

FRESH PLATED HEPATOCYTES

Human Plated Hepatocytes (Collagen coated with collagen overlay)		
Product No.	Description	Size
M91537	Male Human	6-well culture plate
M91538	Male Human	12-well culture plate
M91539	Male Human	24-well culture plate
M91540	Male Human	48-well culture plate
M91541	Male Human	96-well culture plate
F91537	Female Human	6-well culture plate
F91538	Female Human	12-well culture plate
F91539	Female Human	24-well culture plate
F91540	Female Human	48-well culture plate
F91541	Female Human	96-well culture plate

Male Cynomolgus Monkey Plated Hepatocytes (Collagen coated with collagen overlay)		
Product No.	Description	Size
M91569	Male Monkey	6-well culture plate
M91570	Male Monkey	12-well culture plate
M91571	Male Monkey	24-well culture plate
M91572	Male Monkey	48-well culture plate
M91573	Male Monkey	96-well culture plate

Male Beagle Dog Plated Hepatocytes (Collagen coated with collagen overlay)		
Product No.	Description	Size
M91601	Male Beagle	6-well culture plate
M91602	Male Beagle	12-well culture plate
M91603	Male Beagle	24-well culture plate
M91604	Male Beagle	48-well culture plate
M91605	Male Beagle	96-well culture plate

Male Sprague-Dawley Rat Plated Hepatocytes (Collagen coated with collagen overlay)		
Product No.	Description	Size
M91633	Male Rat	6-well culture plate
M91634	Male Rat	12-well culture plate
M91635	Male Rat	24-well culture plate
M91636	Male Rat	48-well culture plate
M91637	Male Rat	96-well culture plate

Male ICR/CD-1 Mouse Plated Hepatocytes (Collagen coated with collagen overlay)		
Product No.	Description	Size
M91665	Male Mouse	6-well culture plate
M91666	Male Mouse	12-well culture plate
M91667	Male Mouse	24-well culture plate
M91668	Male Mouse	48-well culture plate
M91669	Male Mouse	96-well culture plate

**Formats listed above are hepatocytes on collagen coated plates with a collagen overlay. Other matrix configurations are available including collagen coated plate with Matrigel™ overlay and collagen coated plate with no overlay. Inquire with customer service for more information on these and other matrix configurations.*

Product Description:

Freshly isolated hepatocytes represent a living biological system in which coupled Phase I and II chemical metabolism can be evaluated.¹ The hepatocytes are plated on collagen in multi-well plates filled with shipping medium and are available in a number of formats. Plates should be unpacked immediately upon receipt and prepared for later use. The isolation schedule for our fresh animal hepatocytes can be found on the BioIVT animal isolation calendar at can be found online under the resources tab. Our hepatocytes perform the best when used with BioIVT's INVITROGRO hepatocyte media.

Stability: Unpack immediately upon receipt

Storage: 5% CO₂, 37 °C incubator

Procedure:

Medium Preparation

1. Prepare complete INVITROGRO™ HI Medium
 - Thaw the TORPEDO Antibiotic Mix in a 37° C water bath for 3 to 5 minutes.
 - Add 1.0 mL TORPEDO Antibiotic Mix per 45 mL INVITROGRO HI medium.
 - Note: Following the addition of TORPEDO Antibiotic Mix, the shelf life for the complete medium is 7 days.

Unpacking multi-well plates

1. Remove plate from cardboard sleeve.
2. Place plate in a sterile environment (biosafety cabinet).
3. Remove lid and gently peel away sealing film, starting at a corner.
4. Aspirate shipping medium carefully to avoid disrupting cell monolayer. Use a sterile glass Pasteur pipette attached to a vacuum source.
5. Immediately after aspirating, replace shipping medium with an appropriate volume of room-temperature complete INVITROGRO HI medium. Replace medium one plate at a time to avoid drying out the cells. Appropriate volumes are:
 - 100 µL per well for 96-Well plates.
 - 200 µL per well for 48-Well plates.
 - 500 µL per well for 24-Well plates.
 - 1.0 mL per well for 12-Well plates.
 - 2.5 mL per well for 6-Well plates.
6. Cover plate with lid and incubate cells in 5% CO₂, 37 °C incubator at saturating humidity for no less than 2 hours and no longer than 24 hours to allow cells to resume normal physiological function. (If using INVITROGRO HI Medium, it is important to use a 5% CO₂ incubator since this medium is buffered with bicarbonate.)

Related Products:

Product No.	Description	Size
Z99009	INVITROGRO HI Medium	250 mL
Z990012	INVITROGRO HI Medium	500 mL
Z990002	INVITROGRO HI Medium	1 L
Z99000	TORPEDO Antibiotic Mix	5.5 mL
Z990007	TORPEDO Antibiotic Mix	11 mL
Z990008	TORPEDO Antibiotic Mix	22 mL

References:

1. Li, A. P. Primary hepatocyte cultures as an in vitro experimental model for the evaluation of pharmacokinetic drug-drug interactions. *Adv. Pharmacol. Series* **1997**, 43, 103–130.

Caution: Treat all products containing human and monkey-derived materials a potentially infectious, as no known test methods can offer assurance that products derived from human or monkey tissues will not transmit infectious agents.

All products are for research use only. Do not use in animals or humans. These products have not been approved for any diagnostic or clinical procedures.