

INSTRUCTIONS FOR USE

Product Name: INVITROGRO™ CP Rodent Medium

Product No.	Description	Size
Z990028	INVITROGRO CP Rodent Medium	250 mL
Z990027	TORPEDO Rodent Antibiotic Mix	5.5 mL

Product Description

INVITROGRO CP Rodent Medium is used for plating BioIVT rat and mouse cryoplateable hepatocytes. INVITROGRO CP Rodent Medium is formulated without antibiotics to increase its shelf life. BioIVT recommends the use of the TORPEDO™ Rodent Antibiotic Mix (Z990027) with INVITROGRO CP Rodent Medium (Z990028).

Storage

Stability: 3 months for INVITROGRO CP Rodent Medium without TORPEDO Rodent Antibiotic Mix
7 days for Complete INVITROGRO CP Rodent Medium

Storage: 4°C

Handling/Caution Statement

Caution: This product is being sold for research and/or manufacturing purposes only. The biological samples supplied by BioIVT, or any material isolated from the samples, are for in-vitro research use only and are not to be used as a source of material for clinical therapies. Human material may be used in vivo in animals. The user assumes all responsibility for its usage and disposal, in accordance with all regulations.

Related Products

Product No.	Description	Size
M00005-P	Male Sprague-Dawley Rat	5 million cells
M00025-P	Male Wistar Rat	5 million cells
M00065-P	Male Han Wistar Rat	5 million cells
M005052-P	Male ICR/CD-1 Mouse	5 million cells
M005152-P	Male C57BL/6 Mouse	5 million cells