

INSTRUCTIONS FOR USE

Product Name: INVITROGRO™ CP Dog Medium

Product No.	Description	Size
Z990025	INVITROGRO CP Dog Medium	250 mL
Z990000	TORPEDO Antibiotic Mix	5.5 mL

Product Description

INVITROGRO CP Dog Medium is used for plating BioIVT dog cryoplateable hepatocytes. INVITROGRO CP Dog Medium is formulated without antibiotics to increase its shelf life. BioIVT recommends the use of the TORPEDO™ Antibiotic Mix (Z990000) with INVITROGRO CP Medium (Z990025).

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
M0002055-P	Male Beagle Dog Cryoplateable Hepatocytes, Individual	5 million cells