

## INVITROGRO™ HI MEDIUM

Product No.	Description	Size
Z99009	INVITROGRO HI Medium	250 mL
Z990012	INVITROGRO HI Medium	500 mL
Z990002	INVITROGRO HI Medium	1 L

### Product Description:

INVITROGRO HI Medium is a hepatocyte incubation medium used for culturing primary hepatocytes<sup>1</sup>. INVITROGRO HI Medium is formulated without antibiotics to increase its shelf life. BioIVT recommends the use of the TORPEDO™ Antibiotic Mix with INVITROGRO HI Medium.

**Stability:** 6 months for INVITROGRO HI Medium without TORPEDO Antibiotic Mix  
7 days for Complete INVITROGRO HI Medium

**Storage:** 4°C

### Related Products:

Product No.	Size
Z99000	TORPEDO Antibiotic Mix 5.5 ml
Z990007	TORPEDO Antibiotic Mix 11 mL
Z990008	TORPEDO Antibiotic Mix 22 mL

### Reference:

1. Roymans, D.; Van Looveren, C.; Leone, A.; Parker, J. B.; McMillan, M.; Johnson, M. D.; Koganti, A.; Gilissen, R.; Silber, P.; Mannens, G.; Meuldermans, W. Determination of cytochrome P450 1A2 and cytochrome P450 3A4 induction in cryopreserved human hepatocytes. *Biochem. Pharmacol.* **2004**, *67*(3), 427–437.

**Caution:** 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.