

## INVITROGRO™ HT MEDIUM

Product No.	Description	Size
Z99019	INVITROGRO HT Medium	250 mL
Z990005	INVITROGRO HT Medium	500 mL
Z990006	INVITROGRO HT Medium	1 L

### Product Description:

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

**Stability:** 6 months for INVITROGRO HT Medium without TORPEDO Antibiotic Mix  
7 days for Complete INVITROGRO HT 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. Prabhu, S.; Fackett, A.; Lloyd, S.; McClellan, H. A.; Terrell, C. M.; Silber, P. M.; Li, A. P. Identification of glutathione conjugates of troglitazone in human hepatocytes. *Chem. Biol. Interact.* **2002**, *142*(1-2), 83-97.

**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.