

INVITROGRO™ KHB

Product No.	Description	Size
Z99074	INVITROGRO KHB	250 mL

Product Description:

INVITROGRO KHB is a modified Krebs-Henseleit Buffer for use with hepatocytes. BioIVT recommends the use of the INVITROGRO KHB along with our INVITROGRO hepatocyte media when using LIVERPOOL® 5-, 10-, 20- and 50-Donor Pooled Cryopreserved Human Hepatocytes and regular Cryopreserved Human and Animal Hepatocytes.

Stability: 6 months

Storage: 4°C

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.