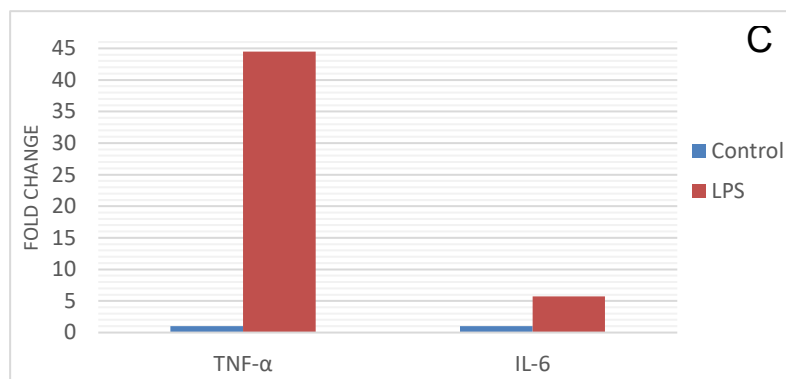
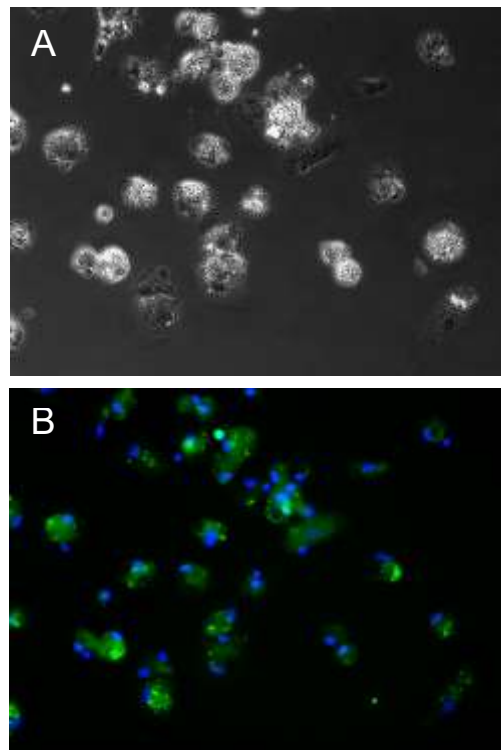


HK1000.H15 Lot No. H1294

Cryopreserved Human Kupffer Cells
Human, Male, Individual

Assured Minimum Yield: 1.0×10^6 cells per vial
Average Viability: 89%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol.



- A) Photomicrograph of cryopreserved Kupffer cells at day 7 in culture (200x).
B) Immuno-fluorescent staining of CD163 expression in cryopreserved Kupffer cells with DAPI nuclear counterstain at day 7 in culture (200x).
C) Expression fold change of TNF-α and IL-6 in cryopreserved Kupffer cells in response to 24 hour 1μg/mL lipopolysaccharide (LPS) treatment.

Donor Information

Gender:	Male
Age:	51 years of age
Race:	Caucasian
Cause of Death:	Cerebrovascular accident
Antibody to Cytomegalovirus (CMV):	Negative
Antibody to Human Immunodeficiency Virus (HIV):	Negative
Hepatitis B Surface Antigen (HbsAg):	Negative
Antibody to Hepatitis C Virus (HCV):	Negative



Store in liquid nitrogen, vapor phase

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for *in vitro* use only.

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 03 November 2016