

H0610.M Lot No. 2310303

Human Liver Mitochondria

Mixed Gender, Pool of 5

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Specific Activities

MAO-A

Substrate

5-Hydroxytryptamine

[S] (μM)

500

5-Hydroxytryptophol formation

10500 ± 200 pmol/mg protein/min

MAO-B

Substrate

4-(Dimethylamino)benzylamine

[S] (μM)

300

4-(Dimethylamino)benzaldehyde formation

670 ± 16 pmol/mg protein/min

Values for enzyme activities are mean ± standard deviation of three or more determinations.

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
332	F	67	Caucasian	Cerebrovascular accident
347	F	49	Caucasian	Cerebrovascular accident
415	M	56	Caucasian	Anoxia
432	M	60	African American	Cerebrovascular accident
469	F	56	Hispanic	Cerebrovascular accident

Serology information

- Cytomegalovirus: 4 donors tested positive.
- RPR*: 5 donors tested negative.
- HIV, HTLV, HbsAg, and HCV**: All donors tested negative.

* Rapid Plasma Reagin

** Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic Virus, Hepatitis B Surface Antigen, Antibody to Hepatitis C Virus, respectively.



Store at -80°C

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.

This data sheet serves as a Certificate of Analysis and has been approved by Stephanie Helmstetter, Assistant Director.

Signature and Date: Stephanie Helmstetter 13 November 2023