

H0606.C(AX) Lot No. 2310090

Human Liver Cytosol Fraction - High AO/XO Activity

Mixed Gender, Pool of 20

1.0 mL at 10 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 2 mM EDTA

Enzyme Activities		Rate
Phthalazine Oxidation	(pmol/mg protein/min)	3470 ± 140
p-Vanillin Oxidation	(pmol/mg protein/min)	1320 ± 30
Zaleplon Oxidation	(pmol/mg protein/min)	164 ± 3
6-Nitroquinazoline Oxidation	(pmol/mg protein/min)	743 ± 21

Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

To measure aldehyde oxidase (AO) activity, liver cytosol samples (0.05 mg/mL) were incubated in triplicate at 37 ± 2°C for 1 minute in potassium phosphate buffer (50 mM, pH 7.4) and phthalazine (25 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure aldehyde oxidase (AO) activity, liver cytosol samples (0.05 mg/mL) were incubated in triplicate at 37 ± 2°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4) and p-Vanillin (20 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure aldehyde oxidase (AO) activity, liver cytosol samples (0.25 mg/mL) were incubated in triplicate at 37 ± 2°C for 30 minutes in potassium phosphate buffer (50 mM, pH 7.4) and Zaleplon (20 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure xanthine oxidase (XO) activity, liver cytosol samples (0.05 mg/mL) were incubated in triplicate at 37 ± 2°C for 5 minutes in potassium phosphate buffer (50 mM, pH 7.4) and 6-Nitroquinazoline (1 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.



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 26 April 2023

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
732	M	54	Caucasian	Head trauma
736	F	46	Caucasian	Anoxia
737	F	49	African American	Anoxia
766	F	38	Caucasian	Anoxia
772	M	46	Caucasian	Anoxia
780	F	48	African American	Cerebrovascular accident
795	M	69	Caucasian	Cerebrovascular accident
803	M	23	African American	Head trauma
805	F	24	African American	Anoxia
814	F	65	Caucasian	Cerebrovascular accident
815	M	52	Caucasian	Anoxia
819	F	49	Caucasian	Head trauma
838	M	51	Caucasian	Cerebrovascular accident
843	M	31	African American	Head trauma
861	F	61	Caucasian	Head trauma
871	M	30	Caucasian	Cerebrovascular accident
873	M	67	Caucasian	Anoxia
886	F	28	African American	Head trauma
887	F	32	African American	Cerebrovascular accident
899	F	49	African American	Anoxia

Serology information

- Cytomegalovirus: 15 donors tested positive
- RPR*: All 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.

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.