

CERTIFICATE OF ANALYSIS
GENTEST® METMAX® MOUSE HEPATOCYTES

Catalog Number	4.82224
Lot Number	2405220

Storage Conditions	Store at -80°C
Date Released	2024 September

Number of Donors	>20
Strain	CD-1
Gender	Male
Volume	2.5 mL
Cell Concentration	2.0 x 10 ⁶ /mL

Drug Metabolism Activity

Metabolic Pathway	Substrate	Substrate Conc. (µM)	Marker Metabolite	Metabolic Activity (pmol/million cells/min)
ECOD	7-Ethoxycoumarin	100	7-OH Coumarin	480
		100	7-Hydroxycoumarin Glucuronide	41
		100	7-Hydroxycoumarin Sulfate	120
UGT	7-Hydroxycoumarin	100	7-Hydroxycoumarin Glucuronide	1800
SULT	7-Hydroxycoumarin	100	7-Hydroxycoumarin Sulfate	13

Drug Metabolism Activity Assessment: Gentest® MetMax® hepatocytes were thawed in a 37°C water bath. A 1000 µL pipette (with tip) was used to pipet up and down three times to achieve homogeneity and then transferred to 1 vial of Cofactor N10/N10+, catalog 4.82212. Contents pipette mixed 2-3 times to ensure homogeneity. Substrate(s) prepared at 2X the desired final concentration. Each substrate added at a volume of 0.05 mL to a 96-well plate. Incubation initiated with addition of 0.05 mL (2x10⁶ cells/mL) Gentest® MetMax® hepatocytes combined with Cofactor N10/N10+ (100,000 hepatocytes/well) and incubated for 30 min in a 37°C, 5% CO₂ incubator and stopped with addition of 0.1 mL of acetonitrile + 0.1% Formic Acid containing internal standard. Samples centrifuged at 4000 rpm at 4°C for 20 minutes and supernatant added to injection plate. Metabolites identified using LC-MS/MS.

SAFETY INFORMATION:

This product is non-hazardous according to US OSHA hazard communication/GHS 29CFR1910.1200. Therefore, a SDS (Safety Data Sheet) is not required. Handle in accordance with good industrial hygiene and laboratory safety practices.



12 September 2024

Quality Assurance

Date