

CERTIFICATE OF ANALYSIS

GENTEST® METMAX® DOG HEPATOCYTES

Catalog Number	4.82225
Lot Number	2405170

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

Number of Donors	4
Strain	Beagle
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)
CYP3A	Midazolam	20	1-Hydroxymidazolam	62
	Testosterone	200	6β-Hydroxytestosterone	120
ECOD	7-Ethoxycoumarin	100	7-OH Coumarin	110
		100	7-Hydroxycoumarin Glucuronide	31
		100	7-Hydroxycoumarin Sulfate	69
UGT	7-Hydroxycoumarin	100	7-Hydroxycoumarin Glucuronide	520
SULT	7-Hydroxycoumarin	100	7-Hydroxycoumarin Sulfate	55

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.



13 September 2024

Quality Assurance

Date

For product inquiries, service requests, or technical support, please contact us at (866) 458-1094 or info@dls.com

For Research Use Only. Not for use in diagnostic or therapeutic procedures.