

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

POOLED MALE RAT LIVER MICROSOMES (Sprague Dawley)

Catalog Number.....452501
Lot Number.....2409271

Storage Conditions.....**STORE AT -80°C**
Date Released.....2024 October

Package Content.....0.5 mL
Protein Content.....20 mg/mL in 250 mM Sucrose

Enzyme Assay Results

Enzyme Measured	Assay	Enzyme Activity [in pmole/(mg x min)]
Total P450	Omura and Sato	890 pmole/mg
OR	Cytochrome c Reductase	400 nmole/(mg x min)
Cytochrome b ₅	Spectrophotometric	380 pmole/mg
CYP3A	Testosterone 6 β -hydroxylase	3600
CYP2C	Testosterone 16 α -hydroxylase	4400
CYP1A	7-Ethoxyresorufin O-Deethylase	210

- Cytochrome P450 assays were conducted at 0.2 mg/mL protein (except 7-ethoxyresorufin O-deethylase assays which was conducted at 0.25 mg/mL) in 0.1 M Potassium Phosphate buffer (pH 7.4) with an NADPH generating system (1.3 mM NADP, 3.3 mM glucose 6-phosphate, 3.3 mM MgCl₂, and 0.4 U/mL glucose 6-phosphate dehydrogenase). Incubation times for the assays were 4 minutes (7-ethoxyresorufin O-deethylase) and 5 minutes (testosterone 6 β -hydroxylase, testosterone 16 α -hydroxylase). Enzyme activities are expressed as pmol product per (mg protein x minute) except cytochrome c reductase which is expressed as nmol product per (mg protein x minute).
- The protein concentration reported has been demonstrated to be within an acceptable range of 18-24 mg/mL. Range reported is due to intrinsic variability in the assay.
- Animals were acclimated on gypsum treated paper bedding for at least seven days prior to use.
- Pool is comprised of livers from 80 rats, 8-10 weeks of age.

Safety Recommendations:

Safety assessment indicates that this product is non-hazardous: therefore no SDS [Safety Data Sheet] is provided. Handle in accordance with good industrial hygiene and laboratory safety practices.



28 October 2024

Quality Assurance

Date