

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

## POOLED MALE MOUSE LIVER MICROSOMES (CD-1)

Catalog Number.....452701  
Lot Number.....2308020

Storage Conditions.....STORE AT -80°C  
Date Released.....2023 November.

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

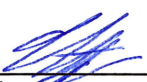
### Enzyme Assay Results

Enzyme Measured	Assay	Enzyme Activity [in pmole/(mg x min)]
Total P450	Omura and Sato	540 pmole/mg
OR	Cytochrome c Reductase	140 nmole/ (mg x min)
Cytochrome b <sub>5</sub>	Spectrophotometric	250 pmole/mg
CYP3A	Testosterone 6 $\beta$ -hydroxylase	2100
CYP1A	7-Ethoxyresorufin O-deethylase	400

- Cytochrome P450 catalytic activity assays were conducted at 0.2 mg/mL protein (testosterone 6 $\beta$ -hydroxylase) or 0.25 mg/mL (7-ethoxyresorufin O-deethylase) in 0.1 M Potassium Phosphate buffer (pH 7.4) with a NADPH regenerating system (1.3 mM NADP<sup>+</sup>, 3.3 mM glucose 6-phosphate, 3.3 mM MgCl<sub>2</sub>, and 0.4 U/mL glucose 6-phosphate dehydrogenase). Assays were conducted in 0.1 M potassium phosphate buffer (pH 7.4). Incubation times for the assays were 4 minutes (7-ethoxyresorufin O-deethylase) and 20 minutes (testosterone 6 $\beta$ -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 unbleached cellulose bedding for at least 7 days prior to use.
- Pool is comprised of livers from approx. 67 mice, 11 weeks of age.

### Safety Recommendations:

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

  
Quality Assurance

14 November 2023  
Date

*For research use only. Not for use in diagnostic or therapeutic procedures.*

© 2013 DLS