

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

# POOLED MALE MOUSE LIVER MICROSOMES (C57BL/6)

Catalog Number.....457145

Storage Conditions.....**STORE AT -80°C**

Lot Number.....2405080

Date Released.....2024 June

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	730 pmole/mg
OR	Cytochrome c Reductase	380 nmole/(mg x min)
cytochrome b <sub>5</sub>	Spectrophotometric	480 pmole/mg

- Enzyme activities are expressed as pmole product per (mg protein x minute) except cytochrome c reductase which is expressed as nmole 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.
- Pool is comprised of livers from approx. 261 mice, 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.



21 June 2024

\_\_\_\_\_  
Quality Assurance

\_\_\_\_\_  
Date

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

© 2013 DLS