

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

POOLED MALE MOUSE LIVER MICROSOMES (CD-1)

Catalog Number.....452701 Storage Conditions.....**STORE AT -80°C**
Lot Number.....2402228 Date Released.....2024 March
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	480 pmole/mg
OR	Cytochrome c Reductase	300 nmole/(mg x min)
Cytochrome b ₅	Spectrophotometric	540 pmole/mg
CYP3A	Testosterone 6β-hydroxylase	3000
CYP1A	7-Ethoxyresorufin O-deethylase	400

- Cytochrome P450 catalytic activity assays were conducted at 0.2 mg/mL protein (testosterone 6β-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⁺, 3.3 mM glucose 6-phosphate, 3.3 mM MgCl₂, 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 6β-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. 662 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.



08 March 2024

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

© 2013 DLS

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

Quality Assurance

Date

Revision History				
Rev.	Change No.	Description of Change	Revised By	Revised Date
1	CC-02544	Updated to Corning branding, protected variable fields, replace safety section with safety statement and clarify SDS [Safety Data Sheet] information, addition of QA signature and date	L. Brown	7/23/14
2	CC 06346	Updated the safety statement	A Fernando	05/26/2017
3	CC 07052	Changed 7-Ethoxyresorufin O-Deethylase protein incubation concentration from 0.5mg/mL to 0.25mg/mL and incubation time from 2 minutes to 4 minutes.	C. McCullough	01/22/2018
4	CC-17526	The following statement was added to the product notes: "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."	C. McCullough	11/19/2020

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

© 2013 DLS