

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

# Human CYP1A2 + P450 Reductase SUPERSOMES™

Catalog Number.....456203  
Lot Number.....2407262

Storage Conditions. STORE AT -80°C  
Date Released .....2024 August  
Expiration Date.....2034 July

Package Contents.....0.5 nmole cytochrome P450 in 0.5 mL

Protein Content.....7.9 mg/mL in 100 mM potassium phosphate (pH 7.4)

Cytochrome c Reductase Activity.....1500 nmole/(min x mg protein)

Cytochrome P450 Content.....1000 pmol per mL

Phenacetin Deethylase Activity.....34 pmol product/(min x pmol P450)

**PRODUCT DESCRIPTION:** This activity is catalyzed by CYP1A2 which is expressed from human CYP1A2 cDNA using a baculovirus expression system. Baculovirus infected insect cells (BTI-TN-5B1-4) were used to prepare these microsomes. A microsome preparation using wild type virus (Catalog No. 456201) should be used as a control for native activities.

## ADVICE:

- Thaw rapidly in a 37°C water bath. Keep on ice until use
- Aliquot to minimize freeze-thawing cycles. Less than 20% of the catalytic activity is lost after 6 freeze thaw cycles.
- Western immunoblotting indicates the expressed CYP1A2 has similar mobility as CYP1A2 in human liver microsomes.
- Comparison of Western immunoblotting intensity and spectral P450 contents for this product and human lymphoblast-expressed CYP1A2 indicates that a substantial amount of apoprotein is found in this product.

**HAZARD WARNING:** The product was produced using baculovirus (*Autographa californica*) infected insect cells (BTI-TN-5B1-4). This virus is not known to be pathogenic to humans or other mammals.

**SAFETY INFORMATION:** Safety assessment indicates this product is not hazardous, therefore no SDS (Safety Data Sheet) is provided. Use standard laboratory practices for the handling and disposal of Biosafety Level 1 materials.

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

© DLS

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

**PRIMARY ASSAY METHOD:** A 0.5 mL reaction mixture containing 10 pmole P450, 1.3 mM NADP<sup>+</sup>, 3.3 mM glucose-6-phosphate, 0.4 U/mL glucose-6-phosphate dehydrogenase, 3.3 mM magnesium chloride and 0.2 mM phenacetin in 100 mM potassium phosphate (pH 7.4) was incubated at 37°C for 10 minutes. After incubation, the reaction was stopped by the addition of 250 µL 10 µM acetamidophenol- <sup>13</sup>C<sub>2</sub> <sup>15</sup>N in acetonitrile with 0.1% formic acid and centrifuged (10,000 x g) for 3 minutes. The product was detected by LC-MS/MS using its Q1 mass and Q3 mass with positive polarity and quantitated by comparing the peak area ratio to a standard curve of acetamidophenol.

### ANALYTICAL METHOD:

#### Materials

Column	2.1 x 50 mm 5µm C18 HPLC
Mobile Phase A	0.1% Formic Acid in dH <sub>2</sub> O
Mobile Phase B	0.1% Formic Acid in Acetonitrile

#### Mass Transitions of MRM

Compound	Q1 Mass (amu)	Q3 Mass (amu)
Analyte- Acetamidophenol	152.0 ±0.2	110.0 ±0.2
Internal Standard- Acetamidophenol- <sup>13</sup> C <sub>2</sub> <sup>15</sup> N	155.0 ±0.2	111.0 ±0.2

#### Gradient Separation Conditions

Time (minute)	Flow Composition of Mobile Phase A (%)	Flow Composition of Mobile Phase B (%)	Flow Rate (µL /min)
0.0	98	2	750
0.4	98	2	750
1.0	60	40	750
1.1	2	98	750
1.3	0	100	750
1.4	98	2	750
1.7	98	2	750

*Alexa Silvas*

Quality Assurance

08/29/2024

Date

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

© DLS