

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

P450 Reductase + Cytochrome b₅ Insect Cell Control SUPERSOMES™

Catalog Number.....456244
Lot Number.....2409190

Storage Conditions. **STORE AT -80°C**
Date Released2024 September
Expiration Date.....2034 September

Package Contents.....0.5 ml
Protein Content.....5.0 mg/ml in 100mM potassium phosphate (pH 7.4)
Cytochrome c Reductase Activity.....1900nmole/(min x mg protein)
Cytochrome b₅ Content.....2500 pmol per mg protein
Cytochrome P450 Content.....Not Detectable

Insect cells (BTI-TN-5B1-4) infected with baculoviruses containing cDNA-expressed human P450 reductase and human cytochrome b₅ were used to prepare these microsomes.

ADVICE

- Thaw rapidly in a 37°C water bath. Keep on ice until use.
- Aliquot to minimize freeze-thawing cycles.
- This product is intended for use as a control for any effect of P450 reductase in the absence of P450. The OR catalytic activity in this product is lower than for some P450 + OR SUPERSOMES, but is twenty fold higher than the OR activity in the insect cell control SUPERSOMES (cat. no. 456201).

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
<https://www.dls.com>

INSECT CELL MICROSOMES

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.



01 October 2024

Quality Assurance

Date

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
<https://www.dls.com>

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

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

© DLS