

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

Insect Cell Control SUPERSOMES™

Catalog Number	457238	LT Catalog Number	P2315
Lot Number	2890313	LT Lot Number	2890313
Date Released	2024 July		
Storage Conditions ...	STORE AT -80°C		

Package Contents.....0.5 ml
Protein Content.....5.0 mg/ml in 100mM potassium phosphate (pH 7.4)
Cytochrome c Reductase Activity.....41 nmole/(min x mg protein)
Cytochrome P450 Content.....Not Detectable

Insect cells (BTI-TN-5B1-4) infected with wild type baculovirus (*Autographa californica*) 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.
- When used as a control for native activities, use protein concentrations comparable to those for active microsome preparations.

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

INSECT CELL MICROSOMES

SAFETY INFORMATION

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 RECOMMENDATIONS:

When using this product, follow good laboratory safety procedures:

Do not eat, drink or smoke.

Avoid contact with skin or eyes.

Do not inhale aerosols.

Do not pipette by mouth.

Wear suitable protective clothing, gloves and eye protection.

Steam sterilize product or treat product with a 1% solution of sodium hypochlorite prior to disposal.



17 July 2024

Quality Assurance

Date

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