

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.....456200
Lot Number.....2405242

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

Package Contents.....1.5 mL
Protein Content.....5.0 mg/mL in 100mM potassium phosphate (pH 7.4)
Cytochrome c Reductase Activity.....38 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.

© 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.



31 May 2024

Quality Assurance

Date

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

© DLS