

ATP-PE-MDR1 1.1 ATP-PE-MRP1 1.1 ATP-PE-MRP2 1.1 ATP-PE-MRP3 1.1 ATP-PE-ratMdr1b 1.1	 SOLVO Biotechnology	
---	---	--

PREDEASY Membrane Product Data Sheet [SB-Beta-gal-Sf9-CTRL]

Catalogue number: SB-PE-Beta-gal-Sf9-CTRL
Description: Isolated Sf9 cell membranes, purified from cells infected with a baculovirus stock containing the β -galactosidase gene.

Date of production (dd.mmm.yyyy):
Expiry date (dd.mmm.yyyy): when stored at -80°C

Packaging: 1 tube containing membrane suspended in TMEP solution.
(TMEP: 50 mM Tris, 50 mM mannitol, 2 mM EGTA, 8 $\mu\text{g/ml}$ aprotinin,
10 $\mu\text{g/ml}$ leupeptin, 50 $\mu\text{g/ml}$ PMSF, 2 mM DTT, pH 7.0.)

Total volume: 180 μl

Protein concentration: 5 mg/ml

Total protein: 0.90 mg

	Normal Range	Specific activity
Basal vanadate sensitive ATPase activity (nmol Pi/mg/min)	0-5	

Intended use: Control membrane for PREDEASY ATPase assay (for SB-MDR1-Sf9, SB-ratMdr1b-Sf9, SB-MRP1-Sf9, SB-MRP2-Sf9 and SB-MRP3-Sf9 membranes)

Methods:
Protein concentrations were determined using the BCA assay. ATPase activity was measured according to the ATPase assay protocol. Protocols are supplied with the membranes.

- Storage and handling:**
- Store at -80°C
 - Thaw membranes in a water bath at 25°C , then store on ice until use.
 - The baseline vanadate sensitive ATPase activity of the membranes does not decrease significantly after one freeze-thaw cycle. If you are planning to reuse the same vial, minimize the number of freeze-thaw cycles by making smaller aliquots.

Validated by:

Date: