

ATP-PE-MDR1 1.1 ATP-PE-ratMdr1b 1.1	 SOLVO Biotechnology	
--	---	--

PREDEASY Membrane Product Data Sheet [SB-defPgp-Sf9-CTRL]

Catalogue number: SB-PE-defPgp-Sf9-CTRL
Description: Isolated Sf9 cell membranes, purified from cells infected with a baculovirus stock containing a defective MDR1/Pgp 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 µg/ml aprotinin, 10 µg/ml leupeptin, 50 µg/ml PMSF, 2 mM DTT, pH 7.0.)

Total volume:

Protein concentration:

Total protein:

	Normal Range	Specific activity
Basal vanadate sensitive ATPase activity (nmol Pi/mg/min)	0-8	<input type="text"/>

Intended use: Control membrane for PREDEASY ATPase assay (for SB-MDR1-Sf9 and SB-ratMdr1b-Sf9)

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: