

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

PREDEASY Membrane Product Data Sheet [defMRP-Sf9-CTRL]

Catalogue number: SB-PE-defMRP-Sf9-CTRL

Description: Isolated Sf9 cell membranes, purified from cells infected with a baculovirus stock containing a defective MRP1 gene.

Date of production (dd.mmm.yyyy):

Expiry date (dd.mmm.yyyy): when stored at $-80\text{ }^{\circ}\text{C}$

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:

Protein concentration:

Total protein:

	Normal range	Specific activity
Basal vanadate sensitive ATPase activity (nmol Pi/mg/min)	0-10	
Modified vanadate sensitive ATPase activity in the presence of the reference substrate (nmol Pi/mg/min)	0-10 no activation	

Intended use: control membrane for PREDEASY ATPase assay (for 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. Vesicular transport was measured according to the corresponding vesicular transport protocol. Protocols are supplied with the membranes.

Storage and handling:

- Store at $-80\text{ }^{\circ}\text{C}$
- Thaw membranes in a water bath at $25\text{ }^{\circ}\text{C}$, 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.
- The vesicular structure of the membrane preparation might be destroyed upon freezing and thawing. If you are using a membrane stock that has been thawed and frozen always include membrane validation in your assay (drug free control – see assay protocol for details).

Validated by:

Date: