

ATP-PE-MRP2 1.1	 SOLVO Biotechnology	
-----------------	---	--

PREDEASY Membrane Product Data Sheet [MRP2- Sf9-ATPase]

Catalogue number: SB-PE-MRP2-Sf9-ATPase

Description: Isolated Sf9 cell membranes, containing human MRP2

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: 300 μl

Protein concentration: 5 mg/ml

Total protein: 1.50 mg

	Normal range:	Specific activity:
Basal vanadate-sensitive ATPase activity [nmol Pi/mg/min]	5-12	
Modified vanadate-sensitive ATPase activity in the presence of reference substrate [nmol Pi/mg/min]	15-25	
-fold activation vs. control	2-3	

Reference substrate [Concentration]: Sulfasalazine (100 μM)

Intended use: for PREDEASY ATPase only

Methods:

Protein concentrations were determined using the BCA assay. Vanadate sensitive background ATPase activity was determined as the difference between Pi liberated in the assay buffer and in the presence of 1.2 mM Na_3VO_4 . Total ATPase activity was determined as the difference between the Pi liberated in the presence of the reference substrate and in the presence of 1.2 mM Na_3VO_4 . See assay protocol for further details.

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 MRP2-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.

Note: We strongly recommend using SB-defMRP-Sf9-CTRL as a transporter negative control.

Validated by:

Date: