

VT-PV-MRP1-Sf9 1.1	 SOLVO Biotechnology	
--------------------	---	--

PREDIV EZ Membrane Product Data Sheet [SB-defMRP-Sf9-CTRL]

Catalogue number:	SB-PV-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):	<input style="width: 100%;" type="text"/>		
Expiry date (dd.mmm.yyyy):	<input style="width: 100%;" type="text"/>	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:	<input style="width: 100%;" type="text" value="200 <math>\mu\text{l}</math>"/>		
Protein concentration:	<input style="width: 100%;" type="text" value="5 mg/ml"/>		
Total protein:	<input style="width: 100%;" type="text" value="1.0 mg"/>		
		Normal Range	Specific activity
ATP dependent B-GS transport (at 5 μM) [pmol/mg/min]	<input style="width: 100%;" type="text" value="5.0-20.0"/>		<input style="width: 100%;" type="text"/>
Intended use:	Control membrane for MRP1 VT assays		

Methods:

Protein concentrations were determined using the BCA assay. ATP dependent B-GS transport was determined as described in the assay protocol (see negative control).

Storage and handling:

- Store at -80°C
- Thaw membranes in a water bath at 25°C , then store on ice until use.
- 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: