

VT-PV-MRP1-Sf9 1.1	 <b>SOLVO</b> Biotechnology	
--------------------	---	--

## PREDIVIZ Membrane Product Data Sheet [MRP1-Sf9-VT]

<b>Catalogue number:</b>	SB-PV-MRP1-Sf9-VT	
<b>Description:</b>	Isolated Sf9 cell membranes, containing human MRP1	
<b>Date of production (dd.mmm.yyyy):</b>	<input type="text"/>	
<b>Expiry date (dd.mmm.yyyy):</b>	<input type="text"/>	when stored at $-80^{\circ}\text{C}$
<b>Packaging:</b>	1 tube containing membrane suspended in TMEP solution. (TMEP: 50 mM Tris, 50 mM mannitol, 2 mM EGTA, 8 $\mu\text{g}/\text{ml}$ aprotinin, 10 $\mu\text{g}/\text{ml}$ leupeptin, 50 $\mu\text{g}/\text{ml}$ PMSF, 2 mM DTT, pH 7.0.)	
<b>Total volume:</b>	<input type="text" value="420 &lt;math&gt;\mu\text{l}&lt;/math&gt;"/>	
<b>Protein concentration:</b>	<input type="text" value="5 mg/ml"/>	
<b>Total protein:</b>	<input type="text" value="2.1 mg"/>	
	<b>Normal range:</b>	<b>Specific activity:</b>
<b>ATP dependent B-GS transport (at 5 <math>\mu\text{M}</math>) [pmol/mg/min]</b>	<input type="text" value="180-300"/>	<input type="text"/>
<b>Intended use:</b>	<input type="text" value="for vesicular transport assay only"/>	

### Methods:

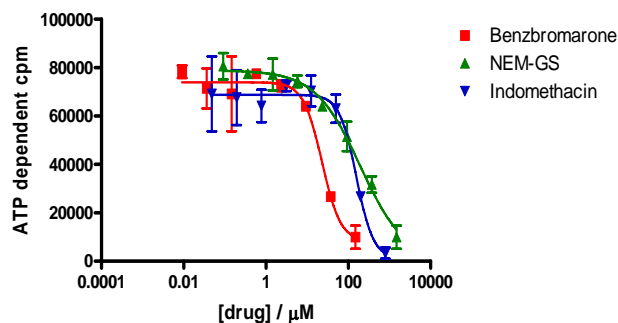
Protein concentrations were determined using the BCA assay. ATP dependent B-GS transport was determined as described in the assay protocol (drug free control).

### Storage and handling:

- Store at  $-80^{\circ}\text{C}$ .
- Thaw membranes in a water bath at  $25^{\circ}\text{C}$ , 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).

**Note:** We strongly recommend using SB-defMRP-Sf9-CTRL as a transporter negative control (included in the Kit).

### Effect of drugs on MRP1-mediated B-GS transport



Validated by:

Date: