

VT-PV-MRP2-Sf9 1.1	 SOLVO Biotechnology	
--------------------	---	--

PREDIVAZ Membrane Product Data Sheet [MRP2-Sf9-VT]

Catalogue number:	SB-PV-MRP2-Sf9-VT	
Description:	Isolated Sf9 cell membranes, containing human MRP2	
Date of production (mm.dd.yyyy):	<input type="text"/>	
Expiry date (mm.dd.yyyy):	<input 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 type="text" value="320 <math>\mu\text{l}</math>"/>	
Protein concentration:	<input type="text" value="5 mg/ml"/>	
Total protein:	<input type="text" value="1.6 mg"/>	
	Normal range:	Specific activity:
ATP dependent CDCF transport (at 5 μM) [pmol/mg/min]	<input type="text" value="10-40"/>	<input type="text"/>
Intended use:	<input type="text" value="for vesicular transport assay only"/>	

Methods:

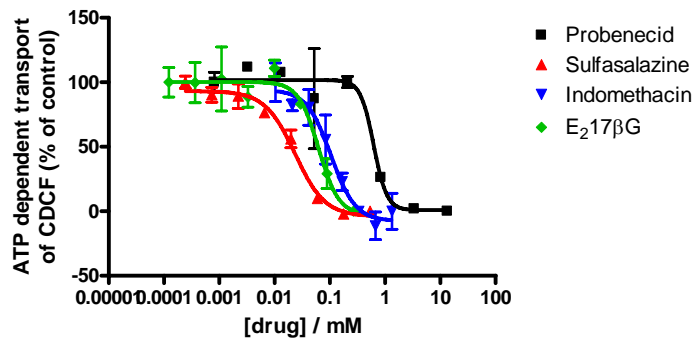
Protein concentrations were determined using the BCA assay. ATP dependent CDCF transport was determined as described in the assay protocol (drug free control).

Storage and handling:

- Store at -80°C .
- Thaw membranes in a water bath at 25°C than 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-Beta-gal-Sf9-CTRL as a transporter negative control (included in the Kit).

Effect of drugs on CDCF transport by MRP2



Validated by:

Date: