

VT-PV-ratMdr1b 1.1	 SOLVO Biotechnology	
--------------------	---	--

PREDIVEZ Membrane Product Data Sheet [ratMdr1b-Sf9-VT]

Catalogue number:	SB-PV-ratMdr1b-Sf9-VT		
Description:	Isolated Sf9 cell membranes, containing rat Mdr1b		
Date of production (dd.mmm.yyyy):	<input type="text"/>		
Expiry date (dd.mmm.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="420 <math>\mu\text{l}</math>"/>		
Protein concentration:	<input type="text" value="5 mg/ml"/>		
Total protein:	<input type="text" value="2.1 mg"/>		
		Normal range:	Specific activity:
ATP dependent $^3\text{H-NMQ}$ transport (at 2 μM) [pmol/mg/min]	<input type="text" value="30-80"/>		<input type="text"/>
Intended use:	<input type="text" value="for vesicular transport assay only"/>		

Methods:

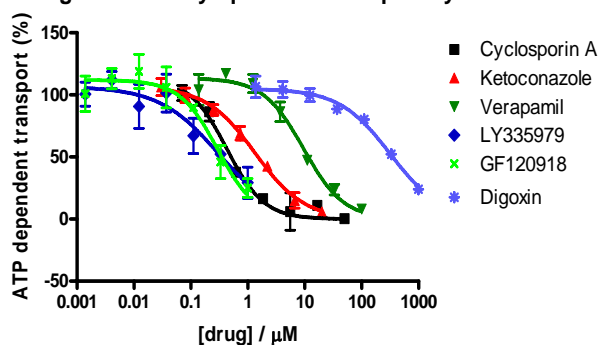
Protein concentrations were determined using the BCA assay. ATP dependent $^3\text{H-NMQ}$ transport was determined as described in the assay protocol (drug free control).

Storage and handling:

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

Effect of drugs on N-methyl-quinidine transport by rat Mdr1b



Validated by:

Date: