

VT-PV-MDR1-K 1.1	 <b>SOLVO</b> Biotechnology	
------------------	---	--

## PREDIVAZ Membrane Product Data Sheet [MDR1-K-VT]

<b>Catalogue number:</b>	SB-PV-MDR1-K-VT	
<b>Description:</b>	Isolated mammalian (K) cell membranes, containing human MDR1/Pgp	
<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 50 mM TrisHCl, 250 mM sucrose (pH 7.4), 8 $\mu\text{g/ml}$ aprotinin, 10 $\mu\text{g/ml}$ leupeptin, 50 $\mu\text{g/ml}$ PMSF	
<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 <math>^3\text{H-NMQ}</math> transport (at 2 <math>\mu\text{M}</math>) [pmol/mg/min]</b>	<input type="text" value="350-500"/>	<input type="text"/>
<b>Intended use:</b>	<input type="text" value="for vesicular transport assay only"/>	

### Methods:

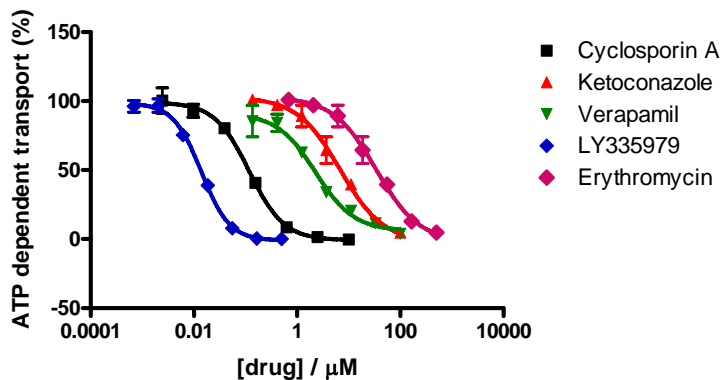
Protein concentrations were determined using the BCA assay. ATP dependent  $^3\text{H-NMQ}$  transport was determined as described in the assay protocol (drug free control).

### 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-K-CTRL as a transporter negative control (included in the Kit).

### Effect of drugs on N-methyl-quinidine transport by MDR1



Validated by:

Date: