

VT-PV-BCRP-M 1.1	 SOLVO Biotechnology	
------------------	-----------------------------------------------------------------------------------------------------------------	--

PREDIV EZ Membrane Product Data Sheet [BCRP-M-VT]

Catalogue number:	SB-PV-BCRP-M-VT	
Description:	Isolated mammalian (M) cell membranes, containing human BCRP	
Date of production (dd.mmm.yyyy):	<input type="text"/>	when stored at -80°C
Expiry date (dd.mmm.yyyy):	<input type="text"/>	
Packaging:	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.)	
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 Lucifer Yellow transport (at 50 μM) [pmol/mg/min]	<input type="text" value="20-35"/>	<input type="text"/>
Intended use:	<input type="text" value="for vesicular transport assay only"/>	

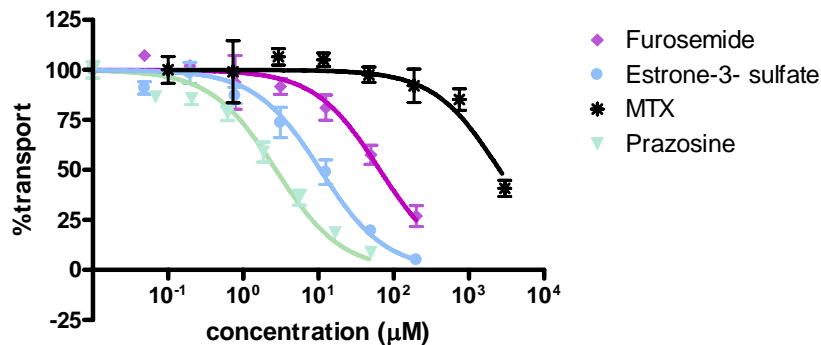
Methods:

Protein concentrations were determined using the BCA assay. ATP dependent Lucifer Yellow 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-M-CTRL as a transporter negative control (included in the Kit).



Validated by:

Date: