

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

PREDIVEZ Membrane Product Data Sheet [defPgp-Sf9-CTRL]

Catalogue number:	SB-PV-defPgp-Sf9-CTRL		
Description:	Isolated Sf9 cell membranes, purified from cells infected with a baculovirus stock containing a defective MDR1/Pgp gene.		
Date of production (dd.mmm.yyyy):			
Expiry date (dd.mmm.yyyy):		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 $\mu\text{g}/\text{ml}$ aprotinin, 10 $\mu\text{g}/\text{ml}$ leupeptin, 50 $\mu\text{g}/\text{ml}$ PMSF, 2 mM DTT, pH 7.0.)		
Total volume:	200 μl		
Protein concentration:	5 mg/ml		
Total protein:	1.0 mg		
		Normal range:	Specific activity:
ATP dependent $^3\text{H-NMQ}$ transport (at 2 μM) [pmol/mg/min]	0-5		
Intended use:	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 (see negative 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).

Validated by:

Date: