

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

PREDIVEZ Membrane Product Data Sheet [SB-M-CTRL]

Catalogue number:	SB-PV-M-CTRL	
Description:	Isolated M cell membranes	
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 50 mM TrisHCl, 250 mM sucrose (pH 7.4)	
Total volume:	<input type="text" value="200 μl"/>	
Protein concentration:	<input type="text" value="5 mg/ml"/>	
Total protein:	<input type="text" value="1.0 mg"/>	
ATP dependent Lucifer Yellow transport (at 50 μM) [pmol/mg/min]	Normal Range	Specific activity
	0.0-1.0	<input type="text"/>
Intended use:	Control membrane for BCRP-VT assay	

Methods:

Protein concentrations were determined using the BCA assay. ATP dependent Lucifer Yellow 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 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).

Validated by:

Date: