

VT-nonHTS-BCRP-M E3S 1.0 VT-HTS-BCRP-M E3S 1.0 VT-nonHTS-BCRP-M MTX 1.0 VT-HTS-BCRP-M MTX 1.0	 SOLVO Biotechnology	
--	---	--

Membrane Product Data Sheet [BCRP-M-VT]

Catalogue number:	SB-BCRP-M-VT
Description:	Isolated mammalian cell membranes, containing human BCRP (ABCG2-482R wild type variant)
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 µg/ml aprotinin, 10 µg/ml leupeptin, 50 µg/ml PMSF, 2 mM DTT, pH 7.0.)
Total volume:	<input type="text" value="500 µl"/>
Protein concentration:	<input type="text" value="5 mg/ml"/>
Total protein:	<input type="text" value="2,5 mg"/>

	Normal range:	Specific activity:
ATP dependent ³ H-estrone-3-sulfate transport (pmol/mg/min)	40-100	
ATP dependent ³ H-methotrexate transport (pmol/mg/min)	40-90	
Intended use:	<input type="text" value="for vesicular transport assay only"/>	

Methods:

Protein concentrations were determined using the BCA assay. ATP dependent ³H-methotrexate and ³H-estrone-3-sulfate 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-M-CTRL as a transporter negative control.

Validated by:

Date: