

VT-HTS-MRP1-2-Sf9-LTC4 1.0 VT-HTS-MRP2-Sf9-E217bG 1.0 VT-nHTS-MRP1-2-Sf9-LTC4 1.0 VT-nHTS-MRP2-Sf9-E217bG 1.0	 SOLVO Biotechnology	
------------------------------------------------------------------------------------------------------------------------	-----------------------------------------------------------------------------------------------------------------	--

Membrane Product Data Sheet [MRP2-Sf9-VT]

Catalogue number:	SB-MRP2-Sf9-VT	
Description:	Isolated Sf9 cell membranes, containing human MRP2	
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:	500 µl	
Protein concentration:	5 mg/ml	
Total protein:	2.5 mg	
	Normal range:	Specific activity:
ATP dependent ³H-E₂17βG transport (at 50 µM) [pmol/mg/min]	500-1000	
ATP dependent ³H-LTC₄ transport (at 50 nM) [pmol/mg/min]	6-20	
Intended use:	for vesicular transport assay only	

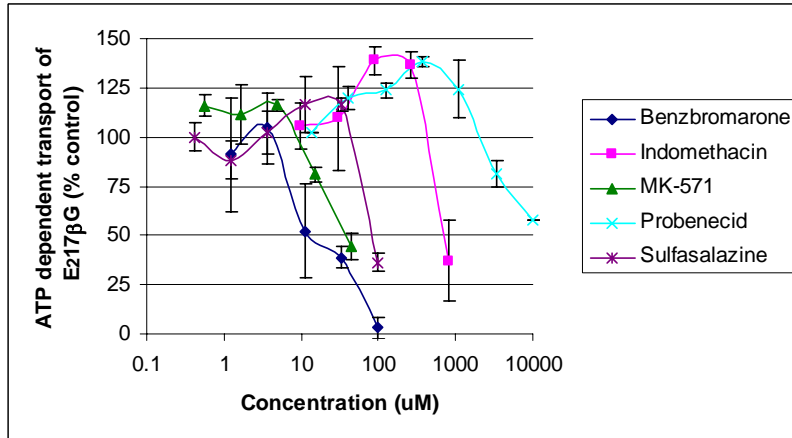
Methods:

Protein concentrations were determined using the BCA assay. ATP dependent ³H-E₂17βG 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-defMRP-Sf9-CTRL as a transporter negative control.



Validated by:

Date: