



## Membrane Product Data Sheet [BCRP-HAM-Sf9-VT]

<b>Catalogue number:</b>	SB-BCRP-HAM-Sf9-VT		
<b>Description:</b>	Isolated Sf9 cell membranes, containing wild type human BCRP		
<b>Date of production (dd.mmm.yyyy):</b>	<input type="text"/>		
<b>Expiry date (dd.mmm.yyyy):</b>	<input type="text"/> when stored at $-80^{\circ}\text{C}$		
<b>Packaging:</b>	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.)		
<b>Total volume:</b>	<input type="text"/> 500 $\mu\text{l}$		
<b>Protein concentration:</b>	<input type="text"/> 5 mg/ml		
<b>Total protein:</b>	<input type="text"/> 2.5 mg		
	<b>Normal range:</b>	<b>Specific activity:</b>	
ATP dependent $^3\text{H}$ -estrone-3-sulfate transport (pmol/mg/min)	100-200	<input type="text"/>	
ATP dependent $^3\text{H}$ -methotrexate transport (pmol/mg/min)	200-350	<input type="text"/>	
<b>Intended use:</b>	<input type="text"/> for vesicular transport assay only		

### Methods:

Protein concentrations were determined using the BCA assay. ATP dependent  $^3\text{H}$ -methotrexate and  $^3\text{H}$ -estrone-3-sulfate transport was determined as described in the assay protocol (drug free control).

### Storage and handling:

- Store at  $-80^{\circ}\text{C}$ .
- Thaw membranes in a water bath at  $25^{\circ}\text{C}$ , 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-defBCRP-HAM-Sf9-CTRL as a transporter negative control.

Validated by:

Date: