

VT-PV-MDR1-K 1.1	 <b>SOLVO</b> Biotechnology	
------------------	---	--

## PREDIV EZ Membrane Product Data Sheet [K-CTRL]

<b>Catalogue number:</b>	SB-PV-K-CTRL		
<b>Description:</b>	Isolated mammalian (K) cell membranes		
<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 50 mM TrisHCl, 250 mM sucrose (pH 7.4), 8 $\mu\text{g/ml}$ aprotinin, 10 $\mu\text{g/ml}$ leupeptin, 50 $\mu\text{g/ml}$ PMSF		
<b>Total volume:</b>	<input type="text"/>	200 $\mu\text{l}$	
<b>Protein concentration:</b>	<input type="text"/>	5 mg/ml	
<b>Total protein:</b>	<input type="text"/>	1.0 mg	
		<b>Normal range:</b>	<b>Specific activity:</b>
<b>ATP dependent <math>^3\text{H-NMQ}</math> transport (at 2 <math>\mu\text{M}</math>) [pmol/mg/min]</b>	<input type="text"/>	0-20	<input type="text"/>
<b>Intended use:</b>	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  $-80^{\circ}\text{C}$ .
- 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).

Validated by:

Date: