

VT-HTS-MRP3-Sf9-E217bG 1.0 VT-nHTS-MRP3-Sf9-E217bG 1.0	 <b>SOLVO</b> Biotechnology	
---	---	--

## Membrane Product Data Sheet [MRP3-Sf9-VT]

**Catalogue number:**

**Description:**

**Date of production (dd.mmm.yyyy):**

**Expiry date (dd.mmm.yyyy):**  when stored at  $-80\text{ }^{\circ}\text{C}$

**Packaging:**

**Total volume:**

**Protein concentration:**

**Total protein:**

	Normal range:	Specific activity:
<b>ATP dependent <math>^3\text{H-E}_217\beta\text{G}</math> transport [pmol/mg/min]</b>	20-35	<input type="text"/>

**Intended use:**

**Methods:**

Protein concentrations were determined using the BCA assay. ATP dependent  $^3\text{H-E}_217\beta\text{G}$  transport was determined as described in the assay protocol (drug free control).

**Storage and handling:**

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

Validated by:

Date: