

VT-PV-MRP5-HEK293 1.1	 SOLVO Biotechnology	
-----------------------	---	--

PREDIVEZ Membrane Product Data Sheet [HEK293-CTRL]

Catalogue number:
Description:

Date of production (dd.mmm.yyyy):
Expiry date (dd.mmm.yyyy): when stored at -80°C

Packaging:

Total volume:

Protein concentration:

Total protein:

	Normal Range	Specific activity
ATP dependent CDCF transport (at 5 µM) [pmol/mg/min]	0-0.5	

Intended use:

Methods:

Protein concentrations were determined using the BCA assay. ATP dependent CDCF transport was determined as described in the assay protocol (see negative 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).

Validated by:

Date: