

ATP-PE-BCRP-HAM 1.1	 SOLVO Biotechnology	
---------------------	---	--

PREDEASY Membrane Product Data Sheet [BCRP-HAM-Sf9-ATPase]

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:
Basal vanadate-sensitive ATPase activity [nmol Pi/mg/min]	10-30	
Modified vanadate-sensitive ATPase activity in the presence of reference substrate [nmol Pi/mg/min]	40-60	
-fold activation vs. control	2-4	

Reference substrate [Concentration]:

Intended use:

Methods:

Protein concentrations were determined using the BCA assay. Vanadate sensitive background ATPase activity was determined as the difference between Pi liberated in the assay buffer and in the presence of 1.2 mM Na_3VO_4 . Modified ATPase activity was determined as the difference between the Pi liberated in the presence of the reference substrates and in the presence of 1.2 mM Na_3VO_4 . See assay protocol for further details.

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 BCRP-HAM-ATPase activity of the membranes does not decrease significantly after one freeze-thaw cycle. If you are planning to reuse the same vial, minimize the number of freeze-thaw cycles by making smaller aliquots.

Note: We strongly recommend using SB-defBCRP-HAM-Sf9-CTRL as a transporter negative control.

Validated by:

Date: