

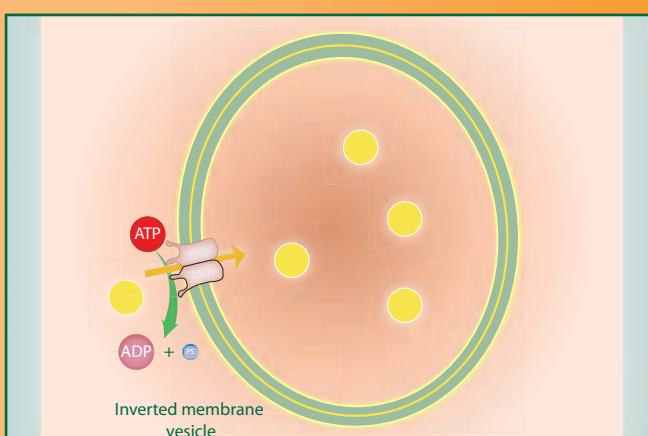
PREDIVEZ™ Vesicular Transport kit

The new **PREDIVEZ™ Vesicular Transport Kit** is a unique product for measuring the drug-transporter interaction in a convenient, “ready to assay” package. The **PREDIVEZ™** kit operates with fluorescent reporter substrate, and contains all you need to conduct vesicular transport assays, including all necessary reagents and solutions.

SOLVO’s PREDIVEZ™ Kit is available in **3-, 6- and 9-compound** sizes.

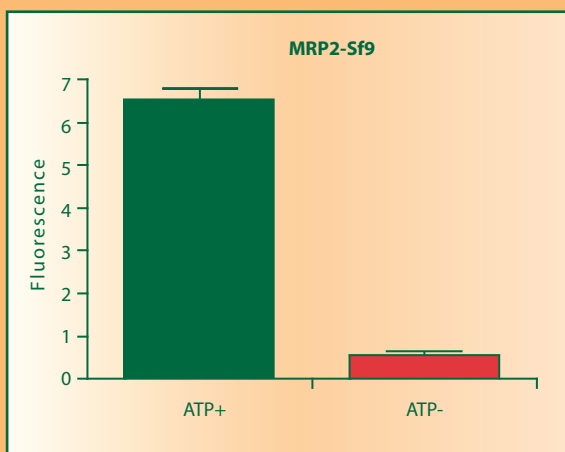
The **PREDIVEZ™ Vesicular Transport kit** is an inhibitory assay. It provides information on any interaction between the ABC transporter and the test drug that would affect the transport of the fluorescent reporter substrate.

The **Fluorescent PREDIVEZ™ Kit** delivers major advantages over the conventional radioactive Vesicular Transport assay



Representation of the vesicular transport assay

ADVANTAGES:



Sensitive and quick fluorescent measurement

With the use of a fluorescent reporter substrate you can avoid radioactive materials and lengthy scintillation counting.

Larger dynamic range / signal-to-background ratio compared to the radioactive Vesicular transport assay

The reporter substrate provides low non-specific binding, so you can get a larger signal and smaller standard deviations for your test compounds.

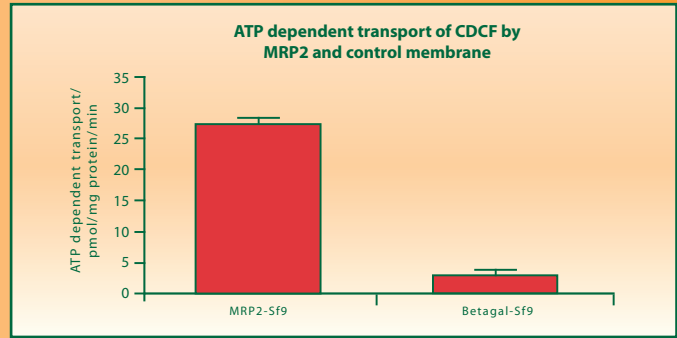
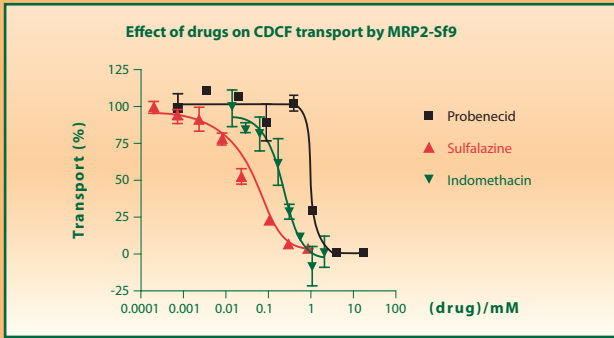
Instant results

Data can be presented in transport values (with the help of the calibration curve, included), or in percentages. Negative control vesicles are also included in the kit for performance monitoring.

3 compounds in 7 concentrations can be screened using one 3-compound Kit.

	MRP2
CDCF	10-15 - fold
E217βG	4-6 - fold

Fold difference of ATP+ and ATP- signal in case of two different reporter substrates.



Reproducible results

- Low inter-assay variability
- Z' factor for MRP2 is ~ 0.74
- Increased automated system compatibility

Recommended equipment:

- 96-well filter plates from Millipore (Cat No. MSFB6B) or equivalent
- Filtration manifold from Millipore (Cat No. MSVMHTS00) or equivalent
- Fluorescent plate reader

Easy to use

All reagents to perform the assay are included!
Every kit comes with a CD containing important information about the PREDIVEZ assay:

- Assay protocol
- Excel data evaluation sheet
- MSDSs
- Reference curves with known interactors



PREDIVEZ Kits are now available for the following transporters

MDR1 PREDIVEZ™ Kit

ratMdr1b PREDIVEZ™ Kit

MRP1 PREDIVEZ™ Kit

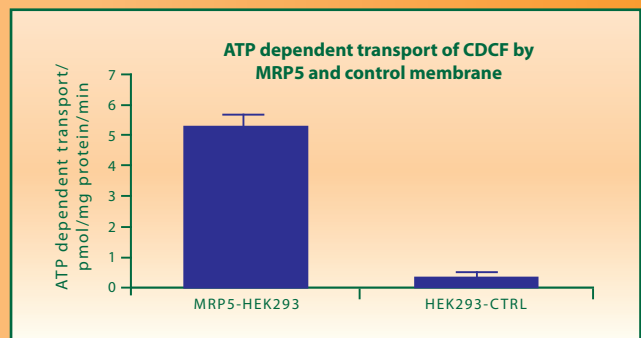
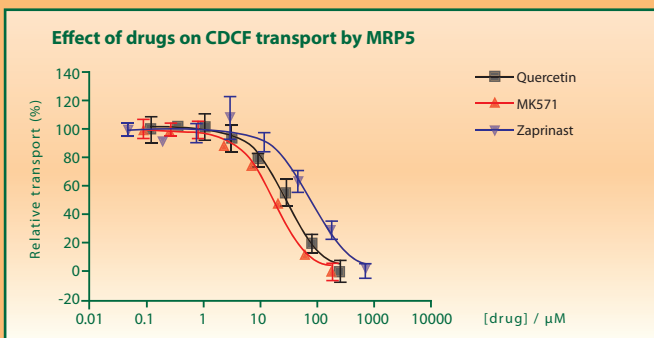
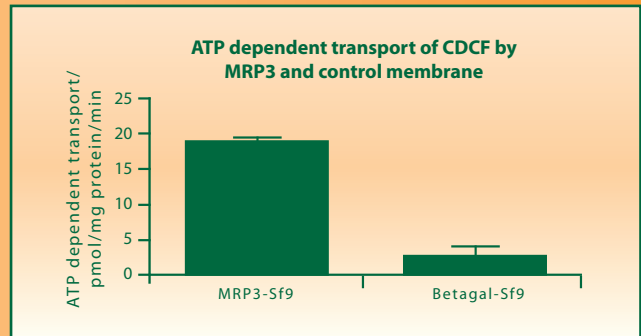
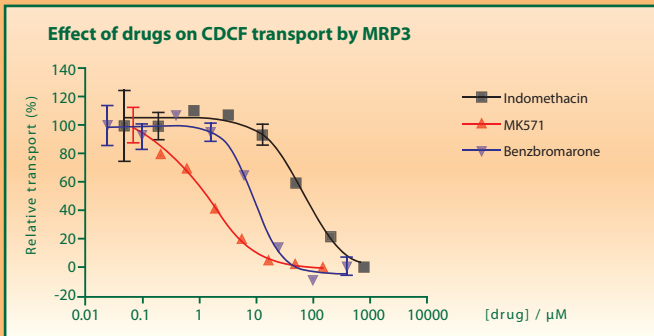
MRP2 PREDIVEZ™ Kit

ratMrp2 PREDIVEZ™ Kit

MRP3 PREDIVEZ™ Kit

MRP5 PREDIVEZ™ Kit

BCRP-M PREDIVEZ™ Kit



CONTACT INFORMATION:

SOLVO Biotechnology

E-mail: sales@solvo.com web: www.solvo.com

Phone: +36-23/503-940 Fax: +36-23/503-941 2 Gyár St. H-2040 Budaörs, HUNGARY

SOLVO distributors: In US and Canada: XenoTech LLC www.xenotechllc.com, in Europe: tebu-bio www.tebu-bio.com, in Japan: Nosan www.nosan.co.jp, in India: www.krishngen.com

