

SOLVO PREDISCREEN SERVICE CATALOGUE

ABC EFFLUX TRANSPORTER ASSAYS

ATPase Assays			
Transporter	Membrane used	Probe (reference substrate)	Positive control activation / inhibition
<i>MDR1</i> /P-gp (ABCB1)	SB-MDR1-Sf9-ATPase	Verapamil	Digoxin / Cyclosporine A
<i>ratMdr1b</i>	SB-ratMdr1b-Sf9-ATPase	Verapamil	Quinidine / Cyclosporine A
<i>MRP1</i> (ABCC1)	SB-MRP1-Sf9-ATPase	NEM-GS	Indomethacin / Benzbromarone
<i>MRP2</i> (ABCC2)	SB-MRP2-Sf9-ATPase	Probenecid	Sulfasalazine / Benzbromarone
<i>ratMrp2</i>	SB-ratMrp2-Sf9-ATPase	Probenecid	Sulfasalazine / Benzbromarone
<i>MRP3</i> (ABCC3)	SB-MRP3-Sf9-ATPase	Benzbromarone	E ₂ 17βG / Cyclosporine A
<i>BCRP</i> (ABCG2)	SB-BRCP-M-ATPase or SB-BCRP-HAM-Sf9-ATPase	Sulfasalazine, Hoechst 33342, Ko134	Topotecan / Ko134
<i>BCRP</i> -R482G (ABCG2-R482G)	SB-BCRP-R482G-Sf9-ATPase	Prazosin, Hoechst 33342, Ko134	Mitoxantrone / Ko134
<i>mouseBsep</i>	SB-mouseBsep-HAM-Sf9-ATPase	TCDC	

NEM-GS: N-Ethylmaleimide S-glutathione, TCDC: taurochenodeoxy-cholate, E₂17βG: estradiol-17-beta-glucuronide

Deliverables: EC₅₀, IC₅₀ • Suggested concentration range of test drug: 137 nM–300 μM (8 concentrations)

Vesicular Transport Assays			
Transporter	Membrane used	Probe (reference substrate)	Positive control
<i>MDR1</i> /P-gp (ABCB1)	SB-MDR1-K-VT	³ H-NMQ	Verapamil
<i>ratMdr1b</i>	SB-ratMdr1b-K-VT	³ H-NMQ	Verapamil
<i>MRP1</i> (ABCC1)	SB-MRP1-Sf9-VT	³ H-LTC4 or radioactive form of test drug (TD)	MK571
<i>MRP2</i> (ABCC2)	SB-MRP2-Sf9-VT	³ H-E ₂ 17βG or radioactive form of TD	Benzbromarone
<i>ratMrp2</i>	SB-ratMrp2-MDCKII-VT	fluorescent CDCF	Benzbromarone
<i>MRP3</i> (ABCC3)	SB-MRP3-Sf9-VT	³ H-E ₂ 17βG or radioactive form of TD	Benzbromarone
<i>MRP5</i> (ABCC5)	SB-MRP5-HEK293-VT	³ H-cGMP (1 μM)	Benzbromarone
<i>BCRP</i> (ABCG2)	SB-BCRP-M-VT	³ H-Estrone-3-sulfate or radioactive form of TD	Methotrexate
<i>BCRP</i> (ABCG2)	SB-BCRP-HAM-Sf9-VT	³ H-Estrone-3-sulfate or ³ H-Methotrexate	Methotrexate
<i>BCRP</i> (ABCG2)	SB-BCRP-Sf9-VT	³ H-Methotrexate or radioactive form of TD	Estrone-3-sulfate
<i>BSEP</i> (ABCB11)	SB-BSEP-Sf9-VT	³ H-Taurocholate or radioactive form of TD	Cyclosporine A
<i>mouseBsep</i>	SB-mouseBsep-Sf9-VT	³ H-Taurocholate	Cyclosporine A

³H-NMQ: ³H-N-methyl-quinidine, ³H-LTC4: ³H-leukotriene C4, ³H-cGMP: ³H-Guanosine 3',5'-cyclic phosphate, CDCF: 5(6)-Carboxy-2',7'-dichlorofluorescein

Deliverables: IC₅₀ – for interacting compounds • (Follow-up studies are available, which provide kinetic data: K_m, K_i and V_{max})
Suggested concentration range of test drug: 0,41 μM–300 μM (7 concentrations)

Dye Transport Assays			
Transporter	Cell line used	Dye used	Positive control
<i>MDR1</i> /P-gp (ABCB1)	HL60-MDR or K562-MDR	Calcein AM	Verapamil
<i>MRP1</i> (ABCC1)	HL60-MRP	Calcein AM	MK571
<i>BCRP</i> (ABCG2)	MXR-M	Hoechst 33342	Proprietary

Deliverables: IC₅₀ • Suggested concentration range of test drug: 0,07 μM–150 μM (8 concentrations)

UPTAKE TRANSPORTER ASSAYS

Transporter	Probe (reference substrate)	Reference inhibitor	Cell type	Negative control
<i>OATP1B1</i> (SLCO1B1)	³ H-Estrone-3-Sulfate	Cerivastatin	CHO	Parental cells
<i>OATP1B3</i> (SLCO1B3)	Fluo-3	Fluvastatin	CHO	Parental cells
<i>OATP2B1</i> (SLCO2B1)	³ H-Estrone-3-Sulfate	Fluvastatin	MDCKII	Parental cells
<i>ratOatp1a1</i> (Slco1a1)	³ H-Estrone-3-Sulfate	Ketoconazole	CHO	Parental cells
<i>NTCP</i> (SLC10A1)	³ H-Taurocholate	TCDC	CHO	Na ⁺ free buffer
<i>ratNtcp</i> (Slc10a1)	³ H-Taurocholate	TCDC	CHO	Na ⁺ free buffer
<i>PEPT1</i> (SLC15A1)	Gly-Sar	Tyr-Phe	CHO	Parental cells
<i>PEPT2</i> (SLC15A2)	Gly-Sar	Cefadroxil	CHO	Parental cells
<i>OAT1</i>	³ H-p-Aminohippuric acid (PAH)	Benzbromarone	CHO	Parental cells
<i>OCT1</i>	¹⁴ C-Tetraethylammonium-chloride (TEA)	Verapamil	CHO	Parental cells
<i>OCT2</i>	¹⁴ C-Tetraethylammonium-chloride (TEA)	Verapamil	CHO	Parental cells

Gly-Sar: Glycylsarcosine, Tyr-Phe: Tyrosine-Phenylalanine

Deliverables: IC₅₀ – for interacting compounds • (Follow-up studies are available, which provide kinetic data: K_m, K_i and V_{max})
Suggested concentration range of test drug: 0,41 μM–300 μM (7 concentrations)

