

MEMBRANE TRANSPORTERS

	Gene symbol	Protein name/alias	Sub-family	Chrom.	LocusID	SOLVO membrane products' catalogue#	Assay	Membrane/Cell
HUMAN	ABCB1	PGY1	MDR/TAP	7q21.1	5243	SB-MDR1-Sf9-ATPase-act	ATPase Assay activation	SB-MDR1-Sf9-ATPase
		MDR1				SB-MDR1-Sf9-ATPase-inhib-I	ATPase Assay inhibition I	SB-MDR1-Sf9-ATPase
		P-GP				K562MDR1 cells-I	Dye Transport Assay I	K562MDR1 cells
		ABC20						
		HGNC:40						
	ABCC1	MRP1	CFTR/MRP	16p13.1	4363	SB-MRP1-Sf9-ATPase-act	ATPase Assay activation	SB-MRP1-Sf9-ATPase
		MRP				SB-MRP1-Sf9-ATPase-inhib-I	ATPase Assay inhibition I	SB-MRP1-Sf9-ATPase
		GS-X				SB-MRP1-Sf9-VT-inhib-I	Vesicular Transport Assay inhibition by test drug I	SB-MRP1-Sf9-VT
		ABC29				SB-MRP1-Sf9-VT-dirtsp	Vesicular Transport Assay direct transport of test drug	SB-MRP1-Sf9-VT
		MRP				HL60MRP1 cells-I	Dye Transport Assay I	HL60MRP1 cells
	ABCC2	DKZp78	CFTR/MRP	10q24	1244	SB-MRP2-Sf9-ATPase-act	ATPase Assay activation	SB-MRP2-Sf9-ATPase
		ABCC				SB-MRP2-Sf9-ATPase-inhib-I	ATPase Assay inhibition I	SB-MRP2-Sf9-ATPase
		1G125				SB-MRP2-Sf9-VT-inhib-I	Vesicular Transport Assay inhibition by test drug I	SB-MRP2-Sf9-VT
		HGNC:51				SB-MRP2-Sf9-VT-dirtsp	Vesicular Transport Assay direct transport of test drug	SB-MRP2-Sf9-VT
	ABCC3	MRP3	CFTR/MRP	17q22	8714	SB-MRP3-Sf9-ATPase-act	ATPase Assay activation	SB-MRP3-Sf9-ATPase
		MLP2				SB-MRP3-Sf9-ATPas-inhib-I	ATPase Assay inhibition I	SB-MRP3-Sf9-ATPase
		ABC31				SB-MRP3-Sf9-VT-inhib-I	Vesicular Transport Assay inhibition by test drug I	SB-MRP3-Sf9-VT
		CMOAT2				SB-MRP3-Sf9-VT-dirtsp	Vesicular Transport Assay direct transport of test drug	SB-MRP3-Sf9-VT
		EST90757						
	ABCC4	MOAT-D	CFTR/MRP	13q32	10257	SB-MRP4-Sf9-VI	Vesicular Transport Assay	SB-MRP4-Sf9-VI
MRP4		SB-MRP4-Sf9-ATPase				ATPase Assay	SB-MRP4-Sf9-ATPase	
RP11-74A12.1								
ABCG2	EST170205	WHITE	4q22	9429	SB-MXR-M-ATPase-act	ATPase Assay activation	SB-MXR-M-ATPase	
	MOAT-B				SB-MXR-M-ATPase-inhib-I	ATPase Assay inhibition I	SB-MXR-M-ATPase	
	BCRP				SB-MXR-Sf9-VT-inhib-I	Vesicular Transport Assay inhibition by test drug I	SB-MXR-Sf9-VT	
	MXR				SB-MXR-Sf9-VT-dirtsp	Vesicular Transport Assay direct transport of test drug	SB-MXR-Sf9-VT	
	MXR1				Mammalian cells-I	Dye Transport Assay I	Mammalian cells	
	ABC1				SB-MXR-R482G-Sf9-ATPase-act	ATPase Assay activation	SB-MXR-R482G-Sf9-ATPase	
	ABCP				SB-MXR-R482G-Sf9-ATPase-inhib-I	ATPase Assay inhibition I	SB-MXR-R482G-Sf9-ATPase	
	ABC15							
	BCRP1							
	HGNC:74							
ABCB11	MGC102821	MDR/TAP	2q24	8647	SB-BSEP-Sf9-VI-inhib-I	Vesicular Transport Assay inhibition by test drug I	SB-BSEP-Sf9-VI	
	EST157481				SB-BSEP-Sf9-VT-dirtsp	Vesicular Transport Assay direct transport of test drug	SB-BSEP-Sf9-VT	
	BDMP							
Abcb1	SPGP	MDR/TAP	4q12	24646	SB-ratMdr1b-Sf9-ATPase-act	ATPase Assay activation	SB-ratMdr1b-Sf9-ATPase	
	PFIC2				SB-ratMdr1b-Sf9-ATPase-inhib-I	ATPase Assay inhibition I	SB-ratMdr1b-Sf9-ATPase	
	PGY4							
	ABC16							
	HGNC: 42							
Abcc2	BSEP	CFTR/MRP	1q51	25303	SB-ratMrp2-Sf9-ATPase-act	ATPase Assay activation	SB-ratMrp2-Sf9-ATPase	
	MRP2				SB-ratMrp2-Sf9-ATPase-inhib-I	ATPase Assay inhibition I	SB-ratMrp2-Sf9-ATPase	
	CMOAT							
	MGI:1352447							
	Abc30							