



# Known G-LISA® Compatible Cells and Tissues By Species

Species	Cell/Tissue Type	Citation
Bovine	Bovine brain microvascular endothelial cells	Higashibata et al., 2006. BMC Biochem. 7, 19
Canine	MDCK cells	Chastre et al., 2009. FASEB J. 23, 916
Chick	Embryonic neural plate tissue	Kinoshita et al., 2008. Mol Biol Cell. 19, 2289
Hamster	Chinese hamster ovary cells	Abu-Elneel et al., 2008. J Biol Chem. 283, 32781
Human	PC-3 human prostate cancer cells	Sequeira et al., 2008. Clin Exp Metastasis. 25, 569
Human	Bone marrow mesenchymal stromal cells	Musso et al., 2010. Haematologica. DOI: 10.3324/haematol.2010.031633
Human	Rectus abdominus muscle tissue	Jin et al., 2011. J Clin Invest. 121, 918
Human	Human microvascular endothelial cells	Schlegel and Waschke, 2010. Microcirculation. 16, 521
Human	Human umbilical vein endothelial cells	Stankiewicz et al., 2010. Biochem Biophys Res Commun. 391, 443
Human	Brain Metastatic Melanoma Cells	Ridgway et al., 2010. J Cell Biochem. 111, 1299
Human	HEp-2 cells	Lesato et al., 2008. Respiration. 76, 434
Human	Human Alveolar Epithelial Type II Cells	Fang et al., 2010. J Biol Chem. 285, 26211
Human	U87 glioblastoma/astrocytoma cells	Pontow et al., 2007. Virology. 368, 1
Human	Primary dermal microvascular endothelial cells	Baumer et al., 2009. J Cell Physiol. 220, 716
Human	Caco-2 E cells, a subclone of Caco-2 BBE	Rapier et al., 2010. Cancer Cell Int. 10, 24
Human	HeLa cells	Mercer and Helenius, 2008. Science. 320, 531
Human	HEK293 cells	Chastre et al., 2009. FASEB J. 23, 916
Human	MDA-MB-231 cells	Zuo et al., 2006. Biochem Biophys Res Commun. 351, 361
Human	Hs578T cells	Zuo et al., 2006. Biochem Biophys Res Commun. 351, 361
Human	Primary brain microvascular endothelial cells	Ramirez et al., 2008. J Immunol. 180, 1854
Mouse	Primary keratinocytes	Nini and Dagnino, 2010. Anal Biochem. 398, 135
Mouse	Frozen tumors from p190B and rtTA/Neu mice	McHenry et al., 2010. Breast Cancer Res. 12, R73
Mouse	Astrocytes in primary culture	Lichtenstein et al., 2010. J Alz Dis. 22, 1135
Mouse	Mouse bone marrow-derived monocytes	Murthy et al., 2009. Am J Physiol Lung Cell Mol Physiol. 297, L846



# Known G-LISA® Compatible Cells and Tissues By Species

Species	Cell/Tissue Type	Citation
Mouse	Primary myocardial endothelial cells	Schlegel et al., 2008. Am J Physiol Cell Physiol. 294, C178
Mouse	Osteoclast cells	Vives et al., 2011. J Bone Miner Res. 26, 1099
Mouse	ATDC5 cells	Woods and Beier, 2006. J Biol Chem. 281, 13134
Mouse	Fibroblast cells	Bradley et al., 2006. Mol Biol Cell. 17, 4827
Rabbit	Renal and femoral artery rings	Alvarez et al., 2010. Br J Pharmacol. 160, 1326
Rat	Hippocampal tissue lysate	Johanna et al., 2010. Neurochem Int. 57, 762
Rat	Aortic rings	Yang et al., 2010. Environ Toxicol Pharmacol. 29, 290
Rat	GP8/3.9 cells	Schreibelt et al., 2007. FASEB J. 21, 3666
Rat	Spinal commissural neurons	Moore et al., 2008. Development. 135, 2855
Rat	Primary pancreatic beta-cells from islet cells	Hammar et al., 2009. Endocrinology. 150, 2072
Rat	B35 neuroblastoma cells	Seifert et al., 2008. Neurosci Lett. 450, 176
Rat	Hippocampal primary neurons	Romero et al., 2010. Toxicol Sci. 118, 602
Rat	Cerebellar primary neurons	Tanaka et al., 2007. J Biol Chem. 282, 10506
Rat	Primary cardiomyocytes	Ganguly et al., 2011. Endocrinology. 152, 247