

Gs-coupled receptor cell lines

Cat: CB-80200-209	Vasoactive Intestinal Peptide Receptor 2 (VIPR2)
Cat: CB-80200-211	Melanocortin 4 Receptor (MC4R)
Cat: CB-80200-215	Melanocortin 5 Receptor (MC5R)
Cat: CB-80200-216	Parathyroid Hormone Receptor 1 (PTH1R)
Cat: CB-80200-217	Glucagon Receptor (GCGR)
Cat: CB-80200-218	Dopamine Receptor 1 (DRD1)
Cat: CB-80200-219	Prostaglandin E Receptor 4 (PTGER4)
Cat: CB-80200-220	Vasoactive Intestinal Peptide Receptor 1 (VIPR1)
Cat: CB-80200-221	Gastric Inhibitor Peptide Receptor (GIPR)
Cat: CB-80200-222	Dopamine Receptor 5 (DRD5)
Cat: CB-80200-228	Parathyroid Hormone Receptor 2 (PTH2R)
Cat: CB-80200-231	Serotonin Receptor 4 (HTR4)
Cat: CB-80200-233	Corticotropin Releasing Hormone Receptor 2 (CRHR2)
Cat: CB-80200-234	Adenylate Cyclase Activating Polypeptide 1 Receptor (ADCYAP1R1)
Cat: CB-80200-235	Secretin Receptor (SCTR)
Cat: CB-80200-236	Follicle Stimulating Hormone Receptor (FSHR)
Cat: CB-80200-237	Serotonin Receptor 6 (HTR6)
Cat: CB-80200-240	Serotonin Receptor Receptor 7 (HTR7)
Cat: CB-80200-241	Glucagon-like Receptor 2 (GLP2R)
Cat: CB-80200-242	Thyroid Stimulating Hormone Receptor (TSHR)
Cat: CB-80200-243	Melanocortin 3 Receptor (MC3R)
Cat: CB-80200-246	Arginine Vasopressin 2 Receptor (AVPR2)
Cat: CB-80200-247	Prostaglandin I2 Receptor (PTGIR)
Cat: CB-80200-248	Glucagon-like Receptor 1 (GLP1R)
Cat: CB-80200-249	Corticotropin Releasing Hormone Receptor 1 (CRHR1)
Cat: CB-80200-251	Calcitonin-like Receptor (CALCRL)
Cat: CB-80200-253	Beta-2 Adrenergic Receptor
Cat: CB-80200-255	Prostaglandin D Receptor 2 (PTGDR2)
Cat: CB-80200-256	PTGER2 (EP2)
Cat: CB-80200-257	Melanocortin 1 Receptor (MC1R)
Cat: CB-80200-258	Calcitonin Receptor (CALCR)
Cat: CB-80200-260	Adenosine A2b Receptor
Cat: CB-80200-261	Adenosine A2a Receptor
Cat: CB-80200-271	Amylin 3 Receptor
Cat: CB-80200-272	Adrenergic receptor beta 1 (ADRB1)
Cat: CB-80200-274	GPR119
Cat: CB-80200-275	TGR5
Cat: CB-80200-275	Luteinizing hormone/Choriogonadotropin receptor (LHCGR)