

Gi-coupled receptor cell lines

Cat: CB-80300-205	Cannabinoid Receptor 1
Cat: CB-80300-206	Dopamine Receptor 2 (DRD2)
Cat: CB-80300-207	Somatostatin Receptor 5 (SSTR5)
Cat: CB-80300-212	Adenosine A1 Receptor
Cat: CB-80300-213	Chemokine (C-C motif) Receptor 5
Cat: CB-80300-214	Melanin Concentrating Hormone Receptor 1 (MCHR1)
Cat: CB-80300-225	Cannabinoid Receptor 2
Cat: CB-80300-230	Opioid receptor, kappa 1
Cat: CB-80300-232	Adenosine A3 Receptor
Cat: CB-80300-239	Glutamate Receptor, Metabotropic 8
Cat: CB-80300-244	Neuropeptide Y Receptor Y1 (NPY1R)
Cat: CB-80300-245	Neuropeptide Y Receptor Y2 (NPY2R)
Cat: CB-80300-254	Sphingolipid G-protein-coupled receptor 1 (EDG1)
Cat: CB-80300-259	Glutamate Receptor, Metabotropic 7
Cat: CB-80300-262	Glutamate Receptor, Metabotropic 2
Cat: CB-80300-265	Chemokine (C-X-C motif) Receptor 4
Cat: CB-80300-266	Opiate Receptor-like 1 (ORL1)
Cat: CB-80300-267	Angiotensin Receptor-like 1
Cat: CB-80300-268	Opioid receptor, Mu
Cat: CB-80300-269	Glutamate Receptor, Metabotropic 4
Cat: CB-80300-270	Neuropeptide Y Receptor Y4 (NPY4R)
Cat: CB-80300-276	Serotonin Receptor 1F (HTR1F)
Cat: CB-80300-277	Serotonin Receptor 1A (HTR1A)
Cat: CB-80300-279	Opioid receptor, delta 1
Cat: CB-80300-280	Neuropeptide Y Receptor Y5 (NPY5R)