

Codex Cat#	Locus ID	Gene_product	Name	NA_info
CB-80100-001	134	adenosine A1 receptor [Homo sapiens] >gi 231473.	adenosine A1 receptor	XM_001687/ NM_000674
CB-80100-002	135	adenosine A2a receptor; adenosine A..	adenosine A2a receptor	NM_000675
CB-80100-003	136	adenosine A2b receptor [Homo sapiens] >gi 11425.	adenosine A2b receptor	XM_008279
CB-80100-004	140	adenosine A3 receptor [Homo sapiens] >gi 114259.	adenosine A3 receptor	XM_052155
CB-80100-005	146	adrenergic alpha-1D- receptor; ad..	alpha-1D-adrenergic receptor	NM_000678
CB-80100-006	147	adrenergic alpha-1B- receptor; ad..	alpha-1B-adrenergic receptor	XM_011248
CB-80100-007	148	Homo sapiens adrenergic, alpha-1A-, receptor (ADRA1A)	alpha-1A-adrenergic receptor, isoform 3	NM_000680
CB-80100-008	150	alpha-2A-adrenergic receptor – human >gi 178196 gb A.	alpha-2A-adrenergic receptor	NM_000681
CB-80100-009	151	(AF316895) alpha 2B adrenergic receptor.	alpha-2B-adrenergic receptor	XM_010871
CB-80100-010	154	(J02960) beta-2 adrenergic receptor [Homo sapiens.	adrenergic, beta-2-, receptor, surface	XM_004030
CB-80100-011	185	angiotensin receptor 1 [Homo sapiens] >gi 67155.	angiotensin receptor 1	XM_051470
CB-80100-012	186	angiotensin receptor 2 [Homo sapiens] >gi 17032.	angiotensin receptor 2	XM_030897
CB-80100-013	187	angiotensin receptor-like 1 [Homo s..	angiotensin receptor-like 1	XM_006349
CB-80100-014	552	arginine vasopressin receptor 1A [Homo sapiens]	arginine vasopressin receptor 1A	XM_006934
CB-80100-015	553	arginine vasopressin receptor 1B; arginine vaso.	arginine vasopressin receptor 1B	NM_000707
CB-80100-016	554	arginine vasopressin receptor 2 [Homo sapiens] .	arginine vasopressin receptor 2	NM_000054
CB-80100-017	623	bradykinin receptor B1	bradykinin receptor B1	XM_007275
CB-80100-018	624	bradykinin receptor B2 [Homo sapiens] >gi 11433.	bradykinin receptor B2	XM_040854
CB-80100-019	643	Burkitt lymphoma receptor 1 GTP-bi..	Burkitt lymphoma receptor 1, isoform 1	NM_001716

CB-80100-020	719	complement component 3a receptor 1	complement component 3a receptor 1	XM_006929
CB-80100-021	728	complement component 5 receptor 1	complement component 5 receptor 1 (C5a ligand)	XM_084025
CB-80100-022	799	(U26553) calcitonin receptor	calcitonin receptor	NM_001742
CB-80100-023	887	cholecystokinin B receptor [Homo	cholecystokinin B receptor	XM_044493
CB-80100-024	1128	cholinergic receptor muscarinic 1	cholinergic receptor, muscarinic 1	XM_057337
CB-80100-025	1129	cholinergic receptor muscarinic 2	cholinergic receptor, muscarinic 2	NM_000739
CB-80100-026	1131	cholinergic receptor muscarinic 3	cholinergic receptor, muscarinic 3	NM_000740
CB-80100-027	1132	(AF385590) m4 muscarinic cholinergic	cholinergic receptor, muscarinic 4	XM_006296
CB-80100-028	1133	cholinergic receptor muscarinic 5;..	cholinergic receptor, muscarinic 5	NM_012125
CB-80100-029	1230	Homo sapiens chemokine (C-C motif) receptor 1 (CCR1)	chemokine (C-C motif) receptor 1	NM_001295
CB-80100-030	1231	chemokine (C-C motif) receptor 2, isoform A	chemokine (C-C motif) receptor 2, isoform A	NM_000647
CB-80100-031	1232	chemokine (C-C motif) receptor 3 [Homo sapiens].	chemokine (C-C motif) receptor 3	NM_001837
CB-80100-032	1234	(AF011504) CCR5 receptor [Homo sapiens]	chemokine (C-C motif) receptor 5	XM_030397
CB-80100-033	1236	chemokine (C-C motif) receptor 7; C	chemokine (C-C motif) receptor 7	XM_049959
CB-80100-034	1237	chemokine (C-C motif) receptor 8; chemokine (C-	chemokine (C-C motif) receptor 8	XM_041048
CB-80100-035	1238	chemokine binding protein 2 [Homo sapiens] >gi .	chemokine binding protein 2	NM_001296
CB-80100-036	1240	chemokine-like receptor 1 [Homo sapiens] >gi 13.	chemokine-like receptor 1	NM_004072
CB-80100-037	1241	leukotriene b4 receptor (chemokine ..	leukotriene b4 receptor (chemokine receptor-like	XM_033240
CB-80100-038	1268	central cannabinoid receptor [Homo sapiens] >gi.	brain cannabinoid receptor 1	XM_004350
CB-80100-039	1269	cannabinoid receptor 2 (macrophage)..	cannabinoid receptor 2 (macrophage)	XM_086356
CB-80100-040	1394	(L23332) corticotropin releasing factor receptor .	corticotropin releasing hormone receptor 1	NM_004382
CB-80100-041	1524	chemokine (C-X3-C) receptor 1; chem..	G protein-coupled receptor 13, chemokine (C-C) receptor-like 1	XM_047502

CB-80100-042	1812	dopamine receptor D1 [Homo sapiens] >gi 118228 .	dopamine receptor D1	XM_084185
CB-80100-043	1813	dopamine receptor D2 [Homo sapiens] >gi 118206 .	dopamine receptor D2, isoform short	NM_016574
CB-80100-044	1814	dopamine receptor D3 [Homo sapiens]..	dopamine receptor D3, isoform b	NM_033658
CB-80100-045	1816	dopamine receptor D5; Dopamine	dopamine receptor D5	NM_000798
CB-80100-046	1880	EBV-induced G protein-coupled recep..	EBV-induced G protein-coupled receptor 2	XM_083897
CB-80100-047	1901	endothelial differentiation sphingolipid G-pro.	G protein-coupled sphingolipid receptor	XM_001499
CB-80100-048	1902	endothelial differentiation lysophosphatidic a.	lysophosphatidic acid receptor EDG2	NM_001401
CB-80100-049	1903	endothelial differentiation sphingolipid G-pro.	endothelial differentiation, sphingolipid	NM_005226
CB-80100-050	1909	endothelin receptor type A [Homo sapiens] >gi 1.	endothelin receptor type A	XM_034331
CB-80100-051	1910	(NM_000115) endothelin receptor	similar to endothelin receptor type B, isoform	NM_000115
CB-80100-052	2149	coagulation factor II receptor precursor [Homo .	coagulation factor II receptor precursor	XM_037260
CB-80100-053	2357	formyl peptide receptor 1	formyl peptide receptor 1	NM_002029
CB-80100-054	2358	formyl peptide receptor-like 1; lipoxin A4 rece.	formyl peptide receptor-like 1	XM_031082
CB-80100-055	2359	formyl peptide receptor-like 2 [Ho..	formyl peptide receptor-like 2	XM_009373
CB-80100-056	2492	follicle stimulating hormone receptor	follicle stimulating hormone receptor	XM_002212
CB-80100-057	2532	Duffy blood group [Homo sapiens]	Duffy blood group	XM_086609
CB-80100-058	2550	gamma-aminobutyric acid (GABA) B receptor 1,	gamma-aminobutyric acid (GABA) B receptor 1,	NM_021905
CB-80100-059	2587	galanin receptor 1	galanin receptor 1	XM_050937
CB-80100-060	2642	glucagon receptor [Homo sapiens] >gi.	glucagon receptor	XM_046488
CB-80100-061	2692	growth hormone releasing hormone receptor	growth hormone releasing hormone receptor	XM_030066
CB-80100-062	2693	growth hormone secretagogue receptor [Homo sapi.	growth hormone secretagogue receptor	XM_003199

CB-80100-063	2696	gastric inhibitory polypeptide recee..	gastric inhibitory polypeptide receptor	NM_000164
CB-80100-064	2740	glucagon-like peptide 1 receptor	glucagon-like peptide 1 receptor	XM_028205
CB-80100-065	2798	gonadotropin-releasing hormone receptor; gonado.	gonadotropin-releasing hormone receptor	XM_003386
CB-80100-066	2825	similar to G protein-coupled recept.	G protein-coupled receptor 1	NM_005279
CB-80100-067	2827	G protein-coupled receptor 3; adenylate cyclase.	G protein-coupled receptor 3	NM_005281
CB-80100-068	2828	G protein-coupled receptor 4	G protein-coupled receptor 4	XM_009140
CB-80100-069	2830	G protein-coupled receptor 6 [Homo sapiens] >gi.	G protein-coupled receptor 6	NM_005284
CB-80100-070	2831	G protein-coupled receptor 7	G protein-coupled receptor 7	NM_005285
CB-80100-071	2832	G protein-coupled receptor 8 [Homo sapiens] >gi.	G protein-coupled receptor 8	NM_005286
CB-80100-072	2833	G protein-coupled receptor 9;	G protein-coupled receptor 9	NM_001504
CB-80100-073	2834	prolactin-releasing hormone recepto..	prolactin-releasing hormone receptor	NM_004248
CB-80100-074	2835	G protein-coupled receptor 12 [Homo sapiens] >g.	G protein-coupled receptor 12	NM_005288
CB-80100-075	2838	G protein-coupled receptor 15 [Homo sapiens] >g.	G protein-coupled receptor 15	NM_005290
CB-80100-076	2840	G protein-coupled receptor 17	G protein-coupled receptor 17	NM_005291
CB-80100-077	2841	PROBABLE G PROTEIN-COUPLED RECE..	G protein-coupled receptor 18	NM_005292
CB-80100-078	2842	G protein-coupled receptor 19	G protein-coupled receptor 19	NM_006143
CB-80100-079	2843	G protein-coupled receptor 20	G protein-coupled receptor 20	NM_005293
CB-80100-080	2844	G protein-coupled receptor 21 [Homo..	G protein-coupled receptor 21	NM_005294
CB-80100-081	2845	G protein-coupled receptor 22	G protein-coupled receptor 22	NM_005295
CB-80100-082	2846	G protein-coupled receptor 23	G protein-coupled receptor 23	XM_018505
CB-80100-083	2847	G protein-coupled receptor 24	G protein-coupled receptor 24	NM_005297
CB-80100-084	2850	super conserved receptor expressed in brain 1 [.	super conserved receptor expressed in brain 1	NM_018971
CB-80100-085	2852	G protein-coupled receptor 30;	G protein-coupled receptor 30	NM_001505
CB-80100-086	2853	G protein-coupled receptor 31 [Homo sapiens] >g.	G protein-coupled receptor 31	NM_005299

CB-80100-087	2854	G protein-coupled receptor 32 [Homo sapiens] >g.	G protein-coupled receptor 32	NM_001506
CB-80100-088	2857	G protein-coupled receptor 34 [Homo sapiens] >g.	G protein-coupled receptor 34	XM_013100
CB-80100-089	2859	G protein-coupled receptor 35	G protein-coupled receptor 35	XM_054157
CB-80100-090	2863	G protein-coupled receptor 39	G protein-coupled receptor 39	NM_001508
CB-80100-091	2864	G protein-coupled receptor 40	G protein-coupled receptor 40	NM_005303
CB-80100-092	2865	G protein-coupled receptor 41	G protein-coupled receptor 41	NM_005304
CB-80100-093	2866	G protein-coupled receptor 42	G protein-coupled receptor 42	
CB-80100-094	2867	G protein-coupled receptor 43 [Homo sapiens] >g.	G protein-coupled receptor 43	NM_005306
CB-80100-095	2912	(AB045011) metabotropic glutamate r..	glutamate receptor, metabotropic 2 precursor	XM_003207
CB-80100-096	2913	metabotropic glutamate receptor 3 precursor	glutamate receptor, metabotropic 3 precursor	NM_000840
CB-80100-097	2914	glutamate receptor metabotropic 4 ..	glutamate receptor, metabotropic 4	NM_000841
CB-80100-098	2917	glutamate receptor metabotropic 7 ..	glutamate receptor, metabotropic 7	XM_041399
CB-80100-099	2918	glutamate receptor metabotropic 8	glutamate receptor, metabotropic 8 precursor	NM_000845
CB-80100-100	2925	gastrin-releasing peptide receptor ..	gastrin-releasing peptide receptor	NM_005314
CB-80100-101	3061	orexin receptor 1	orexin receptor 1	NM_001525
CB-80100-102	3062	orexin receptor 2	orexin receptor 2	NM_001526
CB-80100-103	3269	histamine receptor H1; histamine receptor subc.	histamine receptor H1	XM_052382
CB-80100-104	3274	histamine receptor H2; gastric receptor 1 [Homo.	histamine receptor H2	NM_022304
CB-80100-105	3350	5-hydroxytryptamine (serotonin) receptor 1A [Ho.	5-hydroxytryptamine (serotonin) receptor 1A	XM_003692
CB-80100-106	3351	5-hydroxytryptamine (serotonin) rec..	5-hydroxytryptamine (serotonin) receptor 1B	XM_004117
CB-80100-107	3352	5-hydroxytryptamine (serotonin) receptor 1D [Ho.	5-hydroxytryptamine (serotonin) receptor 1D	XM_001542
CB-80100-108	3354	5-hydroxytryptamine (serotonin) rec..	5-hydroxytryptamine (serotonin) receptor 1E	XM_004134
CB-80100-109	3355	5-hydroxytryptamine (serotonin) receptor 1F; 5-.	5-hydroxytryptamine (serotonin) receptor 1F	NM_000866

CB-80100-110	3356	5-hydroxytryptamine (serotonin) re..	5-hydroxytryptamine (serotonin) receptor 2A	XM_007123
CB-80100-111	3358	5-hydroxytryptamine (serotonin) receptor 2C	5-hydroxytryptamine (serotonin) receptor 2C	NM_000868
CB-80100-112	3360	5-hydroxytryptamine (serotonin) receptor 4 [Hom.	5-hydroxytryptamine (serotonin) receptor 4	XM_052171
CB-80100-113	3361	5-hydroxytryptamine (serotonin) re..	5-hydroxytryptamine (serotonin) receptor 5A	NM_024012
CB-80100-114	3577	interleukin 8 receptor alpha; chemokine (C-X-C.	interleukin 8 receptor, alpha	NM_000634
CB-80100-115	3579	interleukin 8 receptor beta [Homo	interleukin 8 receptor, beta	NM_001557
CB-80100-116	4142	MAS1 oncogene; Oncogene MAS1 [Homo sapiens] >gi.	MAS1 oncogene	XM_004238
CB-80100-117	4157	MELANOCYTE STIMULATING HORMONE RECEPTOR (MS.	melanocortin 1 receptor (alpha melanocyte	XM_047456
CB-80100-118	4158	melanocortin 2 receptor; Melanocortin-2 recepto.	melanocortin 2 receptor	NM_000529
CB-80100-119	4159	melanocortin 3 receptor	melanocortin 3 receptor	NM_019888
CB-80100-120	4160	melanocortin 4 receptor [Homo sapiens]	melanocortin 4 receptor	NM_005912
CB-80100-121	4161	melanocortin 5 receptor [Homo sapiens] >gi 1142.	melanocortin 5 receptor	NM_005913
CB-80100-122	4543	melatonin receptor 1A; melatonin re..	melatonin receptor 1A	XM_038446
CB-80100-123	4886	neuropeptide Y receptor Y1;	neuropeptide Y receptor Y1	XM_059645
CB-80100-124	4887	neuropeptide Y receptor Y2 [Homo sapiens] >gi 1.	neuropeptide Y receptor Y2	XM_011169
CB-80100-125	4923	neurotensin receptor 1 [Homo sapiens]	neurotensin receptor 1	XM_009612
CB-80100-126	4986	opioid receptor kappa 1; Opiate re..	opioid receptor, kappa 1	XM_051505
CB-80100-127	4987	4987_P0011_D11_T7.aln opiate receptor-like 1 [Homo sapiens]	opiate receptor-like 1	XM_047339
CB-80100-128	4989	sigma receptor (SR31747 binding protein 1) [Hom.	opioid receptor, sigma 1	ND
CB-80100-129	5021	oxytocin receptor [Homo sapiens] >..	oxytocin receptor	XM_003177
CB-80100-130	5028	purinergic receptor P2Y G-protein coupled 1 [.	purinergic receptor P2Y, G-protein coupled, 1	NM_002563
CB-80100-131	5029	purinergic receptor P2Y G-protein..	purinergic receptor P2Y, G-protein coupled, 2	XM_054215

CB-80100-132	5030	pyrimidinergic receptor P2Y G-protein coupled .	pyrimidinergic receptor P2Y, G-protein coupled,	NM_002565
CB-80100-133	5031	P2 purinoceptor - human >gi 1407633 gb AAB03572.1 (.	pyrimidinergic receptor P2Y, G-protein coupled,	NM_004154
CB-80100-134	5032	purinergic receptor P2Y G-protein coupled 11 .	purinergic receptor P2Y, G-protein coupled, 11	XM_049496
CB-80100-135	5540	pancreatic polypeptide receptor 1 [Homo sapiens.	pancreatic polypeptide receptor 1	XM_011916
CB-80100-136	5724	platelet-activating factor receptor [Homo sapie.	platelet-activating factor receptor	XM_001857
CB-80100-137	5729	prostaglandin D2 receptor (DP)	prostaglandin D2 receptor (DP)	XM_051711
CB-80100-138	5731	prostaglandin E receptor 1	prostaglandin E receptor 1 (subtype EP1), 42kD	NM_000955
CB-80100-139	5732	similar to prostaglandin E receptor 2 (subtype .	prostaglandin E receptor 2 (subtype EP2), 53kD	XM_051710
CB-80100-140	5733	(D86096) EP3-III [Homo sapiens]	prostaglandin E receptor 3 (subtype EP3)	NM_000957
CB-80100-141	5734	prostaglandin E receptor 4 (subtype EP4) [Homo .	prostaglandin E receptor 4 (subtype EP4)	NM_000958
CB-80100-142	5737	(AB041713) prostaglandin F2alpha rec..	prostaglandin F receptor (FP)	XM_001630
CB-80100-143	5739	prostaglandin I2 (prostacyclin) receptor (IP) [.	prostaglandin I2 (prostacyclin) receptor (IP)	XM_028274
CB-80100-144	5745	parathyroid hormone receptor 1 [Homo sapiens] >.	parathyroid hormone receptor 1	NM_000316
CB-80100-145	5746	parathyroid hormone receptor 2	parathyroid hormone receptor 2	NM_005048
CB-80100-146	6751	somatostatin receptor 1 [Homo sapiens] >gi 1273.	somatostatin receptor 1	NM_001049
CB-80100-147	6752	somatostatin receptor 2	somatostatin receptor 2	XM_085745
CB-80100-148	6753	somatostatin receptor 3 [Homo sapiens] >gi 1141.	somatostatin receptor 3	NM_001051
CB-80100-149	6754	somatostatin receptor 4 [Homo sapiens] >gi 5689.	somatostatin receptor 4	XM_009594
CB-80100-150	6755	somatostatin receptor 5 [Homo sapiens] >gi 1264.	somatostatin receptor 5	NM_001053
CB-80100-151	6865	neurokinin 2 receptor – human >gi 189222 g..	tachykinin receptor 2	XM_005747
CB-80100-152	6869	tachykinin receptor 1 isoform long; Tachykinin.	tachykinin receptor 1, isoform long	NM_001058

CB-80100-153	7201	(AJ011701) thyrotropin-releasing hormone recepto.	thyrotropin-releasing hormone receptor	NM_003301
CB-80100-154	7253	thyroid stimulating hormone recept..	thyroid stimulating hormone receptor	NM_000369
CB-80100-155	7433	vasoactive intestinal peptide recep.	vasoactive intestinal peptide receptor 1	XM_003226
CB-80100-156	7434	vasoactive intestinal peptide receptor 2 [Homo .	vasoactive intestinal peptide receptor 2	NM_003382
CB-80100-157	7852	chemokine (C-X-C motif) receptor 4 (fusin); Ne.	chemokine (C-X-C motif), receptor 4 (fusin)	NM_003467
CB-80100-158	8111	G protein-coupled receptor 68; Ovarian cancer G.	G protein-coupled receptor 68	XM_007383
CB-80100-159	8321	frizzled 1; Frizzled drosophila h..	frizzled 1	NM_003505
CB-80100-160	8326	frizzled 9 [Homo sapiens] >gi 11419..	frizzled 9	XM_004646
CB-80100-161	8477	G protein-coupled receptor 65 [Homo sapiens]	G protein-coupled receptor 65	XM_007392
CB-80100-162	8698	endothelial differentiation G protein coupled .	endothelial differentiation, G protein coupled	XM_009219
CB-80100-163	8834	putative receptor protein [Homo sapiens] >gi 13.	putative receptor protein	XM_008193
CB-80100-164	8843	putative chemokine receptor; GTP-binding protei.	putative chemokine receptor; GTP-binding	XM_015921
CB-80100-165	9002	(AF384819) coagulation fac..	coagulation factor II (thrombin) receptor-like	XM_008930
CB-80100-166	9034	(AF015524) putative chemokine receptor [Homo sapi.	chemokine (C-C motif) receptor-like 2	XM_002926
CB-80100-167	9038	putative neurotransmitter receptor..	putative neurotransmitter receptor	NM_003967
CB-80100-168	9052	retinoic acid induced 3; retinoic acid inducibl	retinoic acid induced 3	NM_003979
CB-80100-169	9248	G protein-coupled receptor 50	G protein-coupled receptor 50	XM_010228
CB-80100-170	9283	endothelin type b receptor-like protein 2 [Homo.	endothelin type b receptor-like protein 2	NM_004767
CB-80100-171	9287	G protein-coupled receptor 58	G protein-coupled receptor 58	NM_014626
CB-80100-172	9288	GPR57 pseudogene	G protein-coupled receptor 57(psuedogene)	NM_014627
CB-80100-173	9289	(BC008770) G protein-coupled..	G protein-coupled receptor 56	XM_048049
CB-80100-174	9290	G protein-coupled receptor 55 [Homo sapiens]	G protein-coupled receptor 55	NM_005683
CB-80100-175	9293	G protein-coupled receptor 52	G protein-coupled receptor 52	XM_001543

CB-80100-176	9294	endothelial differentiation sphingolipid G-pro.	endothelial differentiation, sphingolipid	NM_004230
CB-80100-177	9340	glucagon-like peptide 2 receptor pr..	glucagon-like peptide 2 receptor precursor	NM_004246
CB-80100-178	9934	KIAA0001 gene product; putative G-protein-coupl.	G protein-coupled receptor 105	XM_003091
CB-80100-179	10149	G protein-coupled receptor 64;	G protein-coupled receptor 64	XM_096288
CB-80100-180	10161	purinergic receptor (family A group..	purinergic receptor (family A group 5)	NM_005767
CB-80100-181	10203	calcitonin receptor-like	calcitonin receptor-like	XM_002673
CB-80100-182	10316	G protein-coupled receptor 66 [Homo sapiens] >g.	G protein-coupled receptor 66	XM_087047
CB-80100-183	10430	(BC001118) Similar to seven	Homo sapiens seven transmembrane domain protein (NIFIE14)	NM_032635
CB-80100-184	10663	Homo sapiens chemokine (C-X-C motif) receptor 6 (CXCR6)	chemokine (C-X-C motif) receptor 6	NM_006564
CB-80100-185	10800	cysteinyl leukotriene receptor 1 [Homo sapiens].	cysteinyl leukotriene receptor 1	NM_006639
CB-80100-186	10803	chemokine (C-C motif) receptor 9	chemokine (C-C motif) receptor 9, isoform A	NM_031200
CB-80100-187	10888	(AB040973) KIAA1540 protein [Homo sa..	G protein-coupled receptor 72	XM_045812
CB-80100-188	10936	G protein-coupled receptor 75 [Homo..	G protein-coupled receptor 75	XM_002624
CB-80100-189	11245	putative G protein coupled receptor	putative G protein coupled receptor	NM_007223
CB-80100-190	11250	G protein-coupled receptor 45	G protein-coupled receptor 45	XM_086954
CB-80100-191	11251	G protein-coupled receptor 44	G protein-coupled receptor 44	NM_004778
CB-80100-192	11255	histamine receptor H3	histamine receptor H3	NM_007232
CB-80100-193	11318	Homo sapiens adrenomedullin receptor (ADMR),	adrenomedullin receptor	NM_007264
CB-80100-194	23432	G-protein coupled receptor RE2	G-protein coupled receptor RE2	NM_007369
CB-80100-195	23566	endothelial cell differentiation gene 7; calciu.	endothelial cell differentiation gene 7	NM_012152
CB-80100-196	23620	neurotensin receptor 2; neurotensin receptor t.	neurotensin receptor 2	NM_012344
CB-80100-197	26996	Homo sapiens putative G protein-coupled receptor (GPCR150),	Putative G protein-coupled receptor (GPCR150),	NM_014373
CB-80100-198	27197	G protein-coupled receptor 82	G protein-coupled receptor 82	NM_080817
CB-80100-199	27198	G protein-coupled receptor 81 putative chemokine receptor	G protein-coupled receptor 81	NM_032554

CB-80100-200	27334	putative purinergic receptor [Homo sapiens] >gi.	putative purinergic receptor	NM_014499
CB-80100-201	29909	platelet activating receptor homolog [Homo sapi.	platelet activating receptor homolog	NM_013308
CB-80100-202	29933	G protein-coupled receptor (G2A)	G protein-coupled receptor (G2A)	XM_085103
CB-80100-203	51289	G-protein coupled receptor SALPR; s..	G-protein coupled receptor SALPR; somatostatin	NM_016568
CB-80100-204	51463	putative G-protein coupled receptor (SH120)	putative G-protein coupled receptor (SH120)	NM_016334
CB-80100-205	51554	Homo sapiens chemokine (C-C motif) receptor-like 1 (CCRL1)	chemokine (C-C motif) receptor-like 1	NM_016557
CB-80100-206	51704	G protein-coupled receptor, family C, group 5, member B	G protein-coupled receptor, family C, group 5, member B	NM_016235
CB-80100-207	53637	Homo sapiens endothelial differentiation, sphingolipid G-protein-coupled receptor, 8 (EDG8)	endothelial differentiation, sphingolipid G-protein-coupled receptor, 8	NM_030760
CB-80100-208	53829	G protein-coupled receptor 86	G protein-coupled receptor 86	NM_023914
CB-80100-209	53831	inflammation-related G protein-coupl.	inflammation-related G protein-coupled receptor	NM_020370
CB-80100-210	53836	G protein-coupled receptor 87	G protein-coupled receptor 87	NM_023915
CB-80100-211	54328	super conserved receptor expressed in brain 3 [.	super conserved receptor expressed in brain 3	NM_018969
CB-80100-212	54329	G protein-coupled receptor 85; super conserved .	G protein-coupled receptor 85	XM_033529
CB-80100-213	54982	Homo sapiens ceroid-lipofuscinosis, neuronal 6, late infantile, variant (CLN6)	Ceroid-lipofuscinosis, neuronal 6 (CLN6)	NM_017882
CB-80100-214	55507	Homo sapiens G protein-coupled receptor, family C, group 5, member D (GPRC5D),	G protein-coupled receptor, family C, group 5, member D	NM_018654
CB-80100-215	55868	G protein-coupled receptor C5L2	G protein-coupled receptor C5L2	NM_018485
CB-80100-216	55890	G protein-coupled receptor, family C, group 5,	G protein-coupled receptor, family C, group 5,	NM_018653
CB-80100-217	56413	seven transmembrane receptor BLTR2;..	seven transmembrane receptor BLTR2; leukotriene	NM_019839
CB-80100-218	56670	G protein-coupled receptor 91	G protein-coupled receptor 91	NM_033050

CB-80100-219	56923	neuromedin U receptor 2 [Homo sapie..	neuromedin U receptor 2	NM_020167
CB-80100-220	57007	G PROTEIN-COUPLED RECEPTOR RDC1 HOMOLOG >gi.	G protein-coupled receptor RDC1	XM_051522
CB-80100-221	57105	cysteinyl leukotriene CysLT2 receptor; cDNA: PS.	cysteinyl leukotriene CysLT2 receptor; cDNA:	NM_020377
CB-80100-222	57121	putative G protein-coupled receptor 92 [Homo sa.	putative G protein-coupled receptor 92	NM_020400
CB-80100-223	59340	histamine H4 receptor	histamine H4 receptor	XM_033995
CB-80100-224	59350	leucine-rich repeat-containing G p	leucine-rich repeat-containing G protein-coupled	NM_021634
CB-80100-225	64582	putative leukocyte platelet-activa..	putative leukocyte platelet-activating factor	XM_096782
CB-80100-226	64805	platelet ADP receptor [Homo sapiens] >gi 120839.	platelet ADP receptor	NM_022788
CB-80100-227	81491	GPR63/PSP24 short isoform	brain expressed G-protein-coupled receptor PSP24	NM_030784
CB-80100-228	81491			
CB-80100-229	83550	G protein-coupled receptor 101	G protein-coupled receptor 101	NM_054021
CB-80100-230	83551	G protein-coupled receptor 102	G protein-coupled receptor 102, trace amine receptor 5	NM_053278
CB-80100-231	83873	G protein-coupled receptor 61	G protein-coupled receptor 61	XM_086232
CB-80100-232	84539	G protein-coupled receptor slt	G protein-coupled receptor slt	XM_046588
CB-80100-233	84636	FKSG79 putative purinergic receptor	putative purinergic receptor (FKSG79)	NM_032553
CB-80100-234	117194	G protein-coupled receptor MRGX2	G protein-coupled receptor MRGX2	NM_054030
CB-80100-235	117195	G protein-coupled receptor MRGX3	G protein-coupled receptor MRGX3	NM_054031
CB-80100-236	117196	G protein-coupled receptor MRGX4	G protein-coupled receptor MRGX4	NM_054032
CB-80100-237	119586	Homo sapiens similar to G protein-coupled receptor 26	similar to G protein-coupled receptor 26	XM_061555
CB-80100-238	119685	similar to olfactory receptor-like protein	similar to olfactory receptor-like protein	XM_061617
CB-80100-239	119686	PSGR-like	similar to prostate specific G-protein coupled	XM_061618
CB-80100-240	119769	RA1c		
CB-80100-241	120052	LOC120052		
CB-80100-242	120795	similar to odorant receptor S19	similar to odorant receptor S19	XM_062271
CB-80100-243	120806	similar to olfactory receptor S83	similar to olfactory receptor S83	XM_062282
CB-80100-244	121901	(P2YR1-like)	similar to purinergic receptor P2Y, G-protein	XM_062888
CB-80100-245	124601	Chemokine-like	similar to putative transmembrane receptor	XM_064220
CB-80100-246	128227	similar to G-protein coupled receptor SALPR;	similar to G-protein coupled receptor SALPR;	XM_060898

CB-80100-247	134864	Homo sapiens trace amine receptor 1 (Tar1)	Trace amine receptor 1 (Tar1)	NM_138327
CB-80100-248	139760	Homo sapiens G-protein coupled receptor 2 (GPCR2),	G-protein coupled receptor 2 (GPCR2),	XM_066873
CB-80100-249	143475	similar to MrgX1 G protein-coupled receptor	similar to MrgX1 G protein-coupled receptor	XM_089843
CB-80100-250	143502	Homo sapiens similar to olfactory receptor MOR41-1	Similar to olfactory receptor MOR41-1	XM_089863
CB-80100-251	146127	similar to serpentine receptor (secretin)	similar to serpentine receptor (secretin)	XM_085330
CB-80100-252	159948	MAS1-like(GPR90-like)	Homo sapiens similar to putative G-protein coupled receptor	XM_089955
CB-80100-253	163535	A-like	similar to gene rich cluster, A gene	XM_088965
CB-80100-254	165140	Homo sapiens G-protein coupled receptor TG1019	G-protein coupled receptor TG1019	XM_092406
CB-80100-255	166071	R35-like	hypothetical protein XP_093638	XM_093638
CB-80100-256	167417	LOC167417	Oxytocin-like	XM_094471.2
CB-80100-257	2840V2	Homo sapiens G protein-coupled receptor 17	Homo sapiens G protein-coupled receptor 17	NM_005291
CB-80100-258	AAE27643	Chemokine-like		
CB-80100-259	AAH22401	prostate specific G-protein coupled receptor [H.	PSGR-like3	ND
CB-80100-260	AF380189	trace amine receptor 3 TA3	trace amine receptor 3	AF380189
CB-80100-261	AF380192	trace amine receptor 4 TA4	trace amine receptor 4	AF380192
CB-80100-262	AF380193	Homo sapiens trace amine receptor 5 (TA5) gene	Trace amine receptor 5 (TA5) gene	AF380193
CB-80100-263	AGPCR15	FPRL1-like	similar to circumsporozoite protein	XM_061833
CB-80100-264	AGPCR16	similar to putative G-protein coupled receptor RA1c	similar to putative G-protein coupled receptor RA1c	XM_061671
CB-80100-265	AGPCR35	PSGR-like	PSGR-like	ND
CB-80100-266	AGPCR38	PSGR-like	PSGR-like	ND
CB-80100-267	AGPCR50	Chemokine-like	5-HT1B-like	ND
CB-80100-268	AX136399	RTA-like2	RTA-like2	AX136399