

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG030-01	A01	HQP002833	NM_000675	ADORA2A
QG030-01	A02	HQP003089	NM_000677	ADORA3
QG030-01	A03	HQP004991	NM_001014431	AKT1
QG030-01	A04	HQP011680	NM_001030287	ATF3
QG030-01	A05	HQP016926	NM_022461	AZI2
QG030-01	A06	HQP021725	NM_003921	BCL10
QG030-01	A07	HQP016287	NM_005178	BCL3
QG030-01	A08	HQP017988	NM_000061	BTK
QG030-01	A09	HQP016832	NM_052813	CARD9
QG030-01	A10	HQP022972	NM_000610	CD44
QG030-01	A11	HQP022500	NM_001040280	CD83
QG030-01	A12	HQP000369	NM_000077	CDKN2A
QG030-01	B01	HQP001800	NM_145071	CISH
QG030-01	B02	HQP003218	NM_004383	CSK
QG030-01	B03	HQP003761	NM_001042355	CYLD
QG030-01	B04	HQP004426	NM_001381	DOK1
QG030-01	B05	HQP021953	NM_003974	DOK2
QG030-01	B06	HQP012603	NM_016581	ECSIT
QG030-01	B07	HQP004605	NM_005228	EGFR
QG030-01	B08	HQP022771	NM_004836	EIF2AK3
QG030-01	B09	HQP021526	NM_003824	FADD
QG030-01	B10	HQP018231	NM_003299	HSP90B1
QG030-01	B11	HQP009077	NM_005345	HSPA1A
QG030-01	B12	HQP009098	NM_002156	HSPD1
QG030-01	C01	HQP009102	NM_002157	HSPE1
QG030-01	C02	HQP009642	NM_000877	IL1R1
QG030-01	C03	HQP022241	NM_003856	IL1RL1
QG030-01	C04	HQP009736	NM_002164	INDO
QG030-01	C05	HQP009771	NM_001025242	IRAK1
QG030-01	C06	HQP009776	NM_001570	IRAK2
QG030-01	C07	HQP001408	NM_007199	IRAK3
QG030-01	C08	HQP012441	NM_016123	IRAK4
QG030-01	C09	HQP009778	NM_002198	IRF1
QG030-01	C10	HQP009779	NM_002199	IRF2
QG030-01	C11	HQP009780	NM_001571	IRF3
QG030-01	C12	HQP009782	NM_002200	IRF5
QG030-01	D01	HQP009785	NM_001572	IRF7
QG030-01	D02	HQP009251	NM_002163	IRF8
QG030-01	D03	HQP010033	NM_001014797	KCNMA1
QG030-01	D04	HQP010567	NM_005564	LCN2
QG030-01	D05	HQP018936	NM_024119	LGP2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG030-01	D06	HQP022851	NM_004862	LITAF
QG030-01	D07	HQP010995	NM_002371	MAL
QG030-01	D08	HQP017891	NM_003188	MAP3K7
QG030-01	D09	HQP011559	NM_001018016	MUC1
QG030-01	D10	HQP011603	NM_002468	MYD88
QG030-01	D11	HQP017417	NM_000265	NCF1
QG030-01	D12	HQP000603	NM_006540	NCOA2
QG030-01	E01	HQP007453	NM_007052	NOX1
QG030-01	E02	HQP012143	NM_016931	NOX4
QG030-01	E03	HQP015459	NM_020651	PELI1
QG030-01	E04	HQP015458	NM_021255	PELI2
QG030-01	E05	HQP006417	NM_145065	PELI3
QG030-01	E06	HQP013168	NM_006221	PIN1
QG030-01	E07	HQP008424	NM_013258	PYCARD
QG030-01	E08	HQP016063	NM_006908	RAC1
QG030-01	E09	HQP016088	NM_002880	RAF1
QG030-01	E10	HQP021524	NM_003821	RIPK2
QG030-01	E11	HQP022198	NM_080657	RSAD2
QG030-01	E12	HQP005757	NM_015077	SARM1
QG030-01	F01	HQP016150	NM_021805	SIGIRR
QG030-01	F02	HQP017463	NM_000578	SLC11A1
QG030-01	F03	HQP010965	NM_005585	SMAD6
QG030-01	F04	HQP021399	NM_003745	SOCS1
QG030-01	F05	HQP021889	NM_003955	SOCS3
QG030-01	F06	HQP017673	NM_000582	SPP1
QG030-01	F07	HQP017764	NM_007315	STAT1
QG030-01	F08	HQP017845	NM_003177	SYK
QG030-01	F09	HQP000019	NM_004180	TANK
QG030-01	F10	HQP008428	NM_013254	TBK1
QG030-01	F11	HQP023164	NM_014726	TBKBP1
QG030-01	F12	HQP009617	NM_021649	TICAM2
QG030-01	G01	HQP022424	NM_052864	TIFA
QG030-01	G02	HQP001674	NM_001039661	TIRAP
QG030-01	G03	HQP018145	NM_006290	TNFAIP3
QG030-01	G04	HQP013496	NM_019009	TOLLIP
QG030-01	G05	HQP018234	NM_003300	TRAF3
QG030-01	G06	HQP018237	NM_004620	TRAF6
QG030-01	G07	HQP018684	NM_005082	TRIM25
QG030-01	G08	HQP018383	NM_003348	UBE2N
QG030-01	G09	HQP018384	NM_001032288	UBE2V1
QG030-01	G10	HQP019940	NM_030930	UNC93B1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG030-01	G11	HQP015659	NM_020746	VISA
QG030-01	G12	HQP014742	NM_005400	PRKCE
QG030-01	H01	HGDC		
QG030-01	H02	HGDC		
QG030-01	H03	HQP006940	NM_002046	GAPDH
QG030-01	H04	HQP016381	NM_001101	ACTB
QG030-01	H05	HQP015171	NM_004048	B2M
QG030-01	H06	HQP006171	NM_012423	RPL13A
QG030-01	H07	HQP009026	NM_000194	HPRT1
QG030-01	H08	HQP054253	NR_003286	RN18S1
QG030-01	H09	RT		
QG030-01	H10	RT		
QG030-01	H11	PCR		
QG030-01	H12	PCR		