

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG089-01	A01	HQP020342	NM_031915	SETDB2
QG089-01	A02	HQP019798	NM_030648	SETD7
QG089-01	A03	HQP019797	NM_030647	JHDM1D
QG089-01	A04	HQP019656	NM_025209	EPC1
QG089-01	A05	HQP019357	NM_024827	HDAC11
QG089-01	A06	HQP019284	NM_024757	EHMT1
QG089-01	A07	HQP019201	NM_024670	SUV39H2
QG089-01	A08	HQP019083	NM_024545	SAP130
QG089-01	A09	HQP017124	NM_022743	SMYD3
QG089-01	A10	HQP016908	NM_022455	NSD1
QG089-01	A11	HQP014823	NM_018682	MLL5
QG089-01	A12	HQP014792	NM_018489	ASH1L
QG089-01	B01	HQP014790	NM_018486	HDAC8
QG089-01	B02	HQP014740	NM_018433	JMJD1A
QG089-01	B03	HQP014599	NM_018039	JMJD2D
QG089-01	B04	HQP014093	NM_018091	ELP3
QG089-01	B05	HQP013847	NM_017778	WHSC1L1
QG089-01	B06	HQP013570	NM_019071	ING3
QG089-01	B07	HQP013031	NM_016604	JMJD1B
QG089-01	B08	HQP012861	NM_015401	HDAC7A
QG089-01	B09	HQP012846	NM_016539	SIRT6
QG089-01	B10	HQP008383	NM_014159	SETD2
QG089-01	B11	HQP007502	NM_001025247	TAF5L
QG089-01	B12	HQP007324	NM_019040	ELP4
QG089-01	C01	HQP006172	NM_012330	MYST4
QG089-01	C02	HQP006080	NM_012238	SIRT1
QG089-01	C03	HQP005736	NM_015061	JMJD2C
QG089-01	C04	HQP005682	NM_015013	AOF2

QG089-01	C05	HQP005645	NM_012308	FBXL11
QG089-01	C06	HQP005604	NM_012237	SIRT2
QG089-01	C07	HQP000969	NM_006662	SRCAP
QG089-01	C08	HQP000900	NM_006618	JARID1B
QG089-01	C09	HQP000857	NM_012215	MGEA5
QG089-01	C10	HQP000758	NM_006473	TAF6L
QG089-01	C11	HQP000630	NM_006388	HTATIP
QG089-01	C12	HQP000572	NM_006354	TADA3L
QG089-01	D01	HQP000500	NM_001039619	PRMT5
QG089-01	D02	HQP000023	NM_001015053	HDAC5
QG089-01	D03	HQP000022	NM_006044	HDAC6
QG089-01	D04	HQP023359	NM_014860	SUPT7L
QG089-01	D05	HQP023308	NM_012432	SETDB1
QG089-01	D06	HQP023167	NM_006037	HDAC4
QG089-01	D07	HQP023143	NM_014712	SETD1A
QG089-01	D08	HQP023134	NM_014707	HDAC9
QG089-01	D09	HQP023073	NM_014663	JMJD2A
QG089-01	D10	HQP022927	NM_004898	CLOCK
QG089-01	D11	HQP022336	NM_004739	MTA2
QG089-01	D12	HQP021621	NM_003884	PCAF
QG089-01	E01	HQP021612	NM_003883	HDAC3
QG089-01	E02	HQP020068	NM_004187	JARID1C
QG089-01	E03	HQP020040	NM_006534	NCOA3
QG089-01	E04	HQP019800	NM_003482	MLL2
QG089-01	E05	HQP019414	NM_006766	MYST3
QG089-01	E06	HQP018755	NM_001007257	PRDM2
QG089-01	E07	HQP018521	NM_001042424	WHSC1
QG089-01	E08	HQP018453	NM_021140	UTX
QG089-01	E09	HQP017889	NM_005644	TAF12
QG089-01	E10	HQP017887	NM_006284	TAF10

QG089-01	E11	HQP017884	NM_001015891	TAF9
QG089-01	E12	HQP017873	NM_004606	TAF1
QG089-01	F01	HQP017871	NM_001488	TADA2L
QG089-01	F02	HQP011171	NM_000244	MEN1
QG089-01	F03	HQP009852	NM_004973	JARID2
QG089-01	F04	HQP008746	NM_001527	HDAC2
QG089-01	F05	HQP008745	NM_004964	HDAC1
QG089-01	F06	HQP007253	NM_021078	GCN5L2
QG089-01	F07	HQP005050	NM_004456	EZH2
QG089-01	F08	HQP005049	NM_001991	EZH1
QG089-01	F09	HQP004897	NM_001429	EP300
QG089-01	F10	HQP002920	NM_001079846	CREBBP
QG089-01	F11	HQP017753	NM_000059	BRCA2
QG089-01	F12	HQP020087	NM_004653	JARID1D
QG089-01	G01	HQP005787	NM_001080424	JMJD3
QG089-01	G02	HQP017833	NM_003173	SUV39H1
QG089-01	G03	HQP020846	NM_003599	SUPT3H
QG089-01	G04	HQP021143	NM_001033085	HAT1
QG089-01	G05	HQP021389	NM_003743	NCOA1
QG089-01	G06	HQP021473	NM_003792	EDF1
QG089-01	G07	HQP022740	NM_004824	CDYL
QG089-01	G08	HQP053932	NM_004825	CDY2A
QG089-01	G09	HQP001050	NM_006709	EHMT2
QG089-01	G10	HQP001272	NM_052996	PRDM7
QG089-01	G11	HQP001313	NM_007067	MYST2
QG089-01	G12	HQP005685	NM_015015	JMJD2B
QG089-01	H01	HGDC		
QG089-01	H02	HGDC		
QG089-01	H03	HQP006940	NM_002046	GAPDH
QG089-01	H04	HQP016381	NM_001101	ACTB

QG089-01	H05	HQP015171	NM_004048	B2M
QG089-01	H06	HQP006171	NM_012423	RPL13A
QG089-01	H07	HQP009026	NM_000194	HPRT1
QG089-01	H08	HQP054253	NR_003286	RN18S1
QG089-01	H09	RT		
QG089-01	H10	RT		
QG089-01	H11	PCR		
QG089-01	H12	PCR		