

ExProfile™ Human Brain Cancer Gene qPCR Array

For focused group profiling of human brain cancer related genes expression

Cat. No. QG062-A (3 x 96-well plate, Format A)

Cat. No. QG062-B (3 x 96-well plate, Format B)

Cat. No. QG062-C (3 x 96-well plate, Format C)

Cat. No. QG062-D (3 x 96-well plate, Format D)

Cat. No. QG062-E (3 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 252 unique gene primers deposited in three 96-well plates.

Introduction

The ExProfile human brain cancer gene qPCR array profiles 252 human genes to aberrantly expressed human genes involved in human brain cancer. These genes are carefully chosen for their close cancer correlation based on a thorough literature search of peer-reviewed publications. Abnormal gene expression is often observed in cancer development and progression. The ExProfile human brain cancer array allows researchers to study the cancer-related genes to gain understanding of their roles in brain cancer pathogenesis.

- QG062 plate 01: 84 unique gene PCR primer pairs
- QG062 plate 02: 84 unique gene PCR primer pairs
- QG062 plate 03: 84 unique gene PCR primer pairs

Shipping and storage condition

Shipped at room temperate

Stable for at least 6 months when stored at -20°C

Array format

GeneCopeia provides five qPCR array formats (A, B, C, D, and E) suitable for use with the following real-time cyclers.

Important note: Upon receiving, please check to make sure that the correct array format was ordered to ensure the compatibility with your qPCR instrument.

Plate format	Instrument provider	qPCR instrument model
A (96-well)	Applied Biosystems	5700, 7000, 7300, 7500, 7700, 7900HT (Standard 96-well block), ViiA™7 (Standard 96-well block)
B (96-well)	Applied Biosystems	7500 (Fast block), 7900HT (Fast block), StepOnePlus™, ViiA™7 (Fast block)
C (96-well)	Bio-Rad Laboratories	iCycler iQ®, MyiQ™, iQ™5
D (96-well)	Bio-Rad Laboratories	CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™
E (96-well)	Roche Applied Science	LightCycler® 480 (96-well block)

Quality control

1. Each pair of primers in the ExProfile gene qPCR array has been experimentally validated to yield a single dissociation curve peak and to generate a single amplicon of the correct size for the targeted gene.
2. The positive PCR controls (PCR) have been verified to amplify a single amplicon of the correct size with Ct values around **20±2**.
3. The Spike-in reverse transcription controls (RT) have been verified to amplify a single amplicon of the correct size with Ct values around **20±3**.
4. $R^2 > 0.99$ was observed for high inter/ intra-array reproducibility.

Materials required but not provided

All-in-One™ First-Strand cDNA Synthesis Kit
 All-in-One™ qPCR Mix
 Total RNA extraction kit (RNAzol® RT RNA extraction reagent is recommended)
 DNase/RNase free tips, PCR reaction tubes, 1.5 ml microcentrifuge tubes
 5 ml and 10 ml graduated pipettes, beakers, flasks, and cylinders
 10 µl to 1,000 µl adjustable single channel micropipettes with disposable tips
 5 µl to 20 µl adjustable multichannel micropipette, disposable tips, and reservoir
 qPCR instrument, compatible with gene qPCR arrays ordered

Array layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	TP53	IDH1	EGF	IDH2	MDM2	EGFR	CDKN2A	XRCC1	XRCC3	PTEN	HLA-A	HLA-DRB1
B	RTEL1	MTHFR	ERCC2	PON1	IL13	IL4R	CASP8	XRCC5	XRCC4	SOD1	NFKB1	NCF2
C	NBN	MTR	LIG4	IL6	IL4	GSTM3	PHLDB1	NAT2	ALAD	NQO1	RAD54L	XRCC2
D	XDH	TCN2	STAT1	BRAF	SOD2	SELP	CCL2	CCND1	RAD52	RAC1	PTGS2	ABCB1
E	NOS3	NOS2A	NF2	NCF4	MPO	MMP3	MMP1	LIG1	KRAS	KIR3DL1	KIR2DL3	KIR2DL1
F	APEX1	GLTSCR1	XRCC6	ALOX5	FCGR2A	ERCC1	CYP1A1	CTNBN1	PARP1	IRAK3	CDK4	SOD3
G	NOS1	LPO	ICHTHYIN	EXO1	NAT1	CDH1	LRIG2	CLCA2	KLK4	SLIT2	CD14	NUMBL
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG062 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	DLG5	AURKB	CCNH	SOCS3	TNFSF18	STK19	SOCS2	CBS	ADAM19	SOCS1	TSLP	LMO4
B	CNTNAP1	LNX1	ACAD11	TLR10	DEFB126	SHFM1	C7orf24	XPC	WRN	VHL	VEGFA	VCAM1
C	SCGB1A1	C8G	C8B	C7	TYK2	C5AR1	C4BPB	C4BPA	C2	C1QC	TNF	TLR3
D	TLR1	C1QBP	TGFB1	SULT1A1	STAT4	STAT3	STAT2	BRCA2	SSTR2	BPI	SPP1	SOS1
E	SLC2A11	SMARCB1	DCLRE1C	NOD2	SELPLG	SELE	SECTM1	CCL24	CCL20	CCL18	CCL13	CCL11
F	CCL5	CCL1	SERPINB3	CFB	SERPIN1	RFC1	ZNF350	RB1	BCHE	RAF1	RAD51L1	RAD23B
G	BARD1	PTPRD	SFRS15	PTGDR	GPR108	DHX33	KLK6	KLK7	MAP2K2	MAP2K1	FOXJ2	BTBD2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG062 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	KLK15	TOLLIP	POLB	TLR9	PIK3CA	SERPINI1	PFKFB2	REV1	C1RL	IRAK4	PCNA	DEF6
B	SERPINB2	SERPINE1	OPRM1	NTRK1	NFKBIA	NDUFS2	NAB2	SERPINC1	MYD88	MTRR	MSH2	MRE11A
C	MMP9	MIF	MEN1	CD46	MBP	MAL	CD180	LMO1	LMAN1	LBP	KLKB1	KLK1
D	KIR2DS3	KDR	JAK3	IRAK2	IL12B	IL8RB	IL8RA	IL8	APOE	IGFBP3	FLJ40869	IGF2
E	IGF1R	IGF1	ICAM4	APOA2	HSPA6	HRAS	APC	ACAA1	CD209	HGF	ATG9B	GNB3
F	GLS	SENP3	RAD54B	ALOX15B	ALOX15	IL27	DEFB119	ALOX5AP	ALOX12	LY96	TRAM1	FMO3
G	FMO1	CD93	FGF2	FCER2	FCER1G	FCER1A	ALDH3A1	F5	ERCC4	ERBB3	EPHX1	DRD1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG062 plate 03

Product Data Sheet

- **Gene primer pairs:** 84 wells (A row to G row) are designated for a real-time PCR assay for genes (see the primer list).
- **HK1-6:** Six pre-deposited housekeeping gene (HK1-6) primer pairs, which can be used as endogenous positive controls as well as for array normalization.
- **GDC:** Genomic DNA controls, which can be used to specifically detect genomic DNA contamination with a high level of sensitivity.
- **RT:** Spike-in reverse transcription controls, which can be used to monitor the efficiency of the RT reactions. These pre-deposited primer pairs specifically amplify the cDNA template reversed transcribed from the spike-in control RNA in the sample.
- **PCR:** Positive PCR controls, which are used to verify the PCR efficiency by amplifying the pre-deposited DNA template with its specific pre-deposited primer pairs.

Gene primer list

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-01	A01	HQP018175	NM_000546	TP53
QG062-01	A02	HQP009351	NM_005896	IDH1
QG062-01	A03	HQP004599	NM_001963	EGF
QG062-01	A04	HQP009353	NM_002168	IDH2
QG062-01	A05	HQP011135	NM_002392	MDM2
QG062-01	A06	HQP004605	NM_005228	EGFR
QG062-01	A07	HQP000370	NM_058195	CDKN2A
QG062-01	A08	HQP018562	NM_006297	XRCC1
QG062-01	A09	HQP018564	NM_005432	XRCC3
QG062-01	A10	HQP015535	NM_000314	PTEN
QG062-01	A11	HQP008849	NM_002116	HLA-A
QG062-01	A12	HQP054047	BC008403	HLA-DRB1
QG062-01	B01	HQP013006	NM_016434	RTEL1
QG062-01	B02	HQP011547	NM_005957	MTHFR
QG062-01	B03	HQP004976	NM_000400	ERCC2
QG062-01	B04	HQP013473	NM_000446	PON1
QG062-01	B05	HQP009697	NM_002188	IL13
QG062-01	B06	HQP009664	NM_000418	IL4R
QG062-01	B07	HQP018966	NM_001080124	CASP8
QG062-01	B08	HQP018568	NM_021141	XRCC5
QG062-01	B09	HQP018565	NM_003401	XRCC4
QG062-01	B10	HQP017615	NM_000454	SOD1
QG062-01	B11	HQP011807	NM_003998	NFKB1
QG062-01	B12	HQP011693	NM_000433	NCF2
QG062-01	C01	HQP011687	NM_002485	NBN
QG062-01	C02	HQP011554	NM_000254	MTR
QG062-01	C03	HQP010613	NM_002312	LIG4
QG062-01	C04	HQP009670	NM_000600	IL6
QG062-01	C05	HQP009662	NM_000589	IL4
QG062-01	C06	HQP008483	NM_000849	GSTM3
QG062-01	C07	HQP005843	NM_015157	PHLDB1

Product Data Sheet

QG062-01	C08	HQP001136	NM_000015	NAT2
QG062-01	C09	HQP005011	NM_000031	ALAD
QG062-01	C10	HQP004317	NM_000903	NQO1
QG062-01	C11	HQP020707	NM_003579	RAD54L
QG062-01	C12	HQP018563	NM_005431	XRCC2
QG062-01	D01	HQP018552	NM_000379	XDH
QG062-01	D02	HQP017979	NM_000355	TCN2
QG062-01	D03	HQP017764	NM_007315	STAT1
QG062-01	D04	HQP017733	NM_004333	BRAF
QG062-01	D05	HQP017616	NM_000636	SOD2
QG062-01	D06	HQP016746	NM_003005	SELP
QG062-01	D07	HQP016621	NM_002982	CCL2
QG062-01	D08	HQP016204	NM_053056	CCND1
QG062-01	D09	HQP016087	NM_134424	RAD52
QG062-01	D10	HQP016064	NM_018890	RAC1
QG062-01	D11	HQP015598	NM_000963	PTGS2
QG062-01	D12	HQP013100	NM_000927	ABCB1
QG062-01	E01	HQP011868	NM_000603	NOS3
QG062-01	E02	HQP011866	NM_000625	NOS2A
QG062-01	E03	HQP011778	NM_000268	NF2
QG062-01	E04	HQP011694	NM_000631	NCF4
QG062-01	E05	HQP011309	NM_000250	MPO
QG062-01	E06	HQP011257	NM_002422	MMP3
QG062-01	E07	HQP011255	NM_002421	MMP1
QG062-01	E08	HQP010609	NM_000234	LIG1
QG062-01	E09	HQP010133	NM_004985	KRAS
QG062-01	E10	HQP010095	NM_013289	KIR3DL1
QG062-01	E11	HQP010086	NM_015868	KIR2DL3
QG062-01	E12	HQP010083	NM_014218	KIR2DL1
QG062-01	F01	HQP009061	NM_001641	APEX1
QG062-01	F02	HQP008673	NM_015711	GLTSCR1
QG062-01	F03	HQP006547	NM_001469	XRCC6
QG062-01	F04	HQP006359	NM_000698	ALOX5
QG062-01	F05	HQP005280	NM_021642	FCGR2A
QG062-01	F06	HQP004975	NM_202001	ERCC1
QG062-01	F07	HQP003772	NM_000499	CYP1A1
QG062-01	F08	HQP003539	NM_001904	CTNNB1
QG062-01	F09	HQP003120	NM_001618	PARP1
QG062-01	F10	HQP001408	NM_007199	IRAK3
QG062-01	F11	HQP000245	NM_000075	CDK4
QG062-01	F12	HQP017618	NM_003102	SOD3
QG062-01	G01	HQP011865	NM_000620	NOS1
QG062-01	G02	HQP010851	NM_006151	LPO
QG062-01	G03	HQP054326	NM_001099287	ICHTHYIN
QG062-01	G04	HQP054030	NM_130398	EXO1
QG062-01	G05	HQP023467	NM_000662	NAT1
QG062-01	G06	HQP023466	NM_004360	CDH1
QG062-01	G07	HQP023300	NM_014813	LRIG2

Product Data Sheet

QG062-01	G08	HQP023012	NM_006536	CLCA2
QG062-01	G09	HQP022998	NM_004917	KLK4
QG062-01	G10	HQP022584	NM_004787	SLIT2
QG062-01	G11	HQP022490	NM_000591	CD14
QG062-01	G12	HQP022406	NM_004756	NUMBL
QG062-01	H01	HGDC		
QG062-01	H02	HGDC		
QG062-01	H03	HQP006940	NM_002046	GAPDH
QG062-01	H04	HQP016381	NM_001101	ACTB
QG062-01	H05	HQP015171	NM_004048	B2M
QG062-01	H06	HQP006171	NM_012423	RPL13A
QG062-01	H07	HQP009026	NM_000194	HPRT1
QG062-01	H08	HQP054253	NR_003286	RN18S1
QG062-01	H09	RT		
QG062-01	H10	RT		
QG062-01	H11	PCR		
QG062-01	H12	PCR		
QG062-02	A01	HQP022361	NM_004747	DLG5
QG062-02	A02	HQP022323	NM_004217	AURKB
QG062-02	A03	HQP021906	NM_001239	CCNH
QG062-02	A04	HQP021889	NM_003955	SOCS3
QG062-02	A05	HQP021834	NM_005092	TNFSF18
QG062-02	A06	HQP021636	NM_032454	STK19
QG062-02	A07	HQP021602	NM_003877	SOCS2
QG062-02	A08	HQP021518	NM_000071	CBS
QG062-02	A09	HQP021484	NM_033274	ADAM19
QG062-02	A10	HQP021399	NM_003745	SOCS1
QG062-02	A11	HQP021260	NM_033035	TSLP
QG062-02	A12	HQP021225	NM_006769	LMO4
QG062-02	B01	HQP021130	NM_003632	CNTNAP1
QG062-02	B02	HQP020906	NM_032622	LNX1
QG062-02	B03	HQP020499	NM_032169	ACAD11
QG062-02	B04	HQP019982	NM_001017388	TLR10
QG062-02	B05	HQP019941	NM_030931	DEFB126
QG062-02	B06	HQP019268	NM_006304	SHFM1
QG062-02	B07	HQP018854	NM_024051	C7orf24
QG062-02	B08	HQP018556	NM_004628	XPC
QG062-02	B09	HQP018545	NM_000553	WRN
QG062-02	B10	HQP018485	NM_000551	VHL
QG062-02	B11	HQP018475	NM_001025366	VEGFA
QG062-02	B12	HQP018466	NM_080682	VCAM1
QG062-02	C01	HQP018411	NM_003357	SCGB1A1
QG062-02	C02	HQP018391	NM_000606	C8G
QG062-02	C03	HQP018380	NM_000066	C8B
QG062-02	C04	HQP018353	NM_000587	C7
QG062-02	C05	HQP018340	NM_003331	TYK2
QG062-02	C06	HQP018326	NM_001736	C5AR1
QG062-02	C07	HQP018276	NM_000716	C4BPB

Product Data Sheet

QG062-02	C08	HQP018259	NM_000715	C4BPA
QG062-02	C09	HQP018226	NM_000063	C2
QG062-02	C10	HQP018170	NM_172369	C1QC
QG062-02	C11	HQP018141	NM_000594	TNF
QG062-02	C12	HQP018115	NM_003265	TLR3
QG062-02	D01	HQP018113	NM_003263	TLR1
QG062-02	D02	HQP018106	NM_001212	C1QBP
QG062-02	D03	HQP018044	NM_000660	TGFB1
QG062-02	D04	HQP017811	NM_177536	SULT1A1
QG062-02	D05	HQP017770	NM_003151	STAT4
QG062-02	D06	HQP017767	NM_003150	STAT3
QG062-02	D07	HQP017766	NM_005419	STAT2
QG062-02	D08	HQP017753	NM_000059	BRCA2
QG062-02	D09	HQP017744	NM_001050	SSTR2
QG062-02	D10	HQP017702	NM_001725	BPI
QG062-02	D11	HQP017673	NM_000582	SPP1
QG062-02	D12	HQP017625	NM_005633	SOS1
QG062-02	E01	HQP017549	NM_030807	SLC2A11
QG062-02	E02	HQP017526	NM_003073	SMARCB1
QG062-02	E03	HQP016982	NM_001033855	DCLRE1C
QG062-02	E04	HQP016801	NM_022162	NOD2
QG062-02	E05	HQP016747	NM_003006	SELPLG
QG062-02	E06	HQP016744	NM_000450	SELE
QG062-02	E07	HQP016738	NM_003004	SECTM1
QG062-02	E08	HQP016644	NM_002991	CCL24
QG062-02	E09	HQP016639	NM_004591	CCL20
QG062-02	E10	HQP016637	NM_002988	CCL18
QG062-02	E11	HQP016630	NM_005408	CCL13
QG062-02	E12	HQP016629	NM_002986	CCL11
QG062-02	F01	HQP016626	NM_002985	CCL5
QG062-02	F02	HQP016620	NM_002981	CCL1
QG062-02	F03	HQP016581	NM_006919	SERPINB3
QG062-02	F04	HQP016565	NM_001710	CFB
QG062-02	F05	HQP016500	NM_001032295	SERPING1
QG062-02	F06	HQP016224	NM_002913	RFC1
QG062-02	F07	HQP016170	NM_021632	ZNF350
QG062-02	F08	HQP016131	NM_000321	RB1
QG062-02	F09	HQP016108	NM_000055	BCHE
QG062-02	F10	HQP016088	NM_002880	RAF1
QG062-02	F11	HQP016082	NM_133509	RAD51L1
QG062-02	F12	HQP016076	NM_002874	RAD23B
QG062-02	G01	HQP015946	NM_000465	BARD1
QG062-02	G02	HQP015912	NM_001040712	PTPRD
QG062-02	G03	HQP015625	NM_020706	SFRS15
QG062-02	G04	HQP015537	NM_000953	PTGDR
QG062-02	G05	HQP015248	NM_001080452	GPR108
QG062-02	G06	HQP015240	NM_020162	DHX33
QG062-02	G07	HQP015102	NM_001012965	KLK6

Product Data Sheet

QG062-02	G08	HQP015096	NM_005046	KLK7
QG062-02	G09	HQP014909	NM_030662	MAP2K2
QG062-02	G10	HQP014907	NM_002755	MAP2K1
QG062-02	G11	HQP014732	NM_018416	FOXJ2
QG062-02	G12	HQP014543	NM_017797	BTBD2
QG062-02	H01	HGDC		
QG062-02	H02	HGDC		
QG062-02	H03	HQP006940	NM_002046	GAPDH
QG062-02	H04	HQP016381	NM_001101	ACTB
QG062-02	H05	HQP015171	NM_004048	B2M
QG062-02	H06	HQP006171	NM_012423	RPL13A
QG062-02	H07	HQP009026	NM_000194	HPRT1
QG062-02	H08	HQP054253	NR_003286	RN18S1
QG062-02	H09	RT		
QG062-02	H10	RT		
QG062-02	H11	PCR		
QG062-02	H12	PCR		
QG062-03	A01	HQP014462	NM_017509	KLK15
QG062-03	A02	HQP013496	NM_019009	TOLLIP
QG062-03	A03	HQP013422	NM_002690	POLB
QG062-03	A04	HQP013388	NM_017442	TLR9
QG062-03	A05	HQP013150	NM_006218	PIK3CA
QG062-03	A06	HQP013133	NM_005025	SERPINI1
QG062-03	A07	HQP013071	NM_006212	PFKFB2
QG062-03	A08	HQP012755	NM_001037872	REV1
QG062-03	A09	HQP012584	NM_016546	C1RL
QG062-03	A10	HQP012441	NM_016123	IRAK4
QG062-03	A11	HQP012420	NM_002592	PCNA
QG062-03	A12	HQP012166	NM_022047	DEF6
QG062-03	B01	HQP012155	NM_002575	SERPINB2
QG062-03	B02	HQP012154	NM_000602	SERPINE1
QG062-03	B03	HQP012062	NM_001008505	OPRM1
QG062-03	B04	HQP011938	NM_001007792	NTRK1
QG062-03	B05	HQP011810	NM_020529	NFKBIA
QG062-03	B06	HQP011728	NM_004550	NDUFS2
QG062-03	B07	HQP011663	NM_005967	NAB2
QG062-03	B08	HQP011620	NM_000488	SERPINC1
QG062-03	B09	HQP011603	NM_002468	MYD88
QG062-03	B10	HQP011555	NM_002454	MTRR
QG062-03	B11	HQP011491	NM_000251	MSH2
QG062-03	B12	HQP011320	NM_005590	MRE11A
QG062-03	C01	HQP011263	NM_004994	MMP9
QG062-03	C02	HQP011219	NM_002415	MIF
QG062-03	C03	HQP011171	NM_000244	MEN1
QG062-03	C04	HQP011113	NM_002389	CD46
QG062-03	C05	HQP011085	NM_001025081	MBP
QG062-03	C06	HQP010996	NM_022438	MAL
QG062-03	C07	HQP010932	NM_005582	CD180

Product Data Sheet

QG062-03	C08	HQP010690	NM_002315	LMO1
QG062-03	C09	HQP010660	NM_005570	LMAN1
QG062-03	C10	HQP010560	NM_004139	LBP
QG062-03	C11	HQP010104	NM_000892	KLKB1
QG062-03	C12	HQP010100	NM_002257	KLK1
QG062-03	D01	HQP010092	NM_012313	KIR2DS3
QG062-03	D02	HQP010070	NM_002253	KDR
QG062-03	D03	HQP009851	NM_000215	JAK3
QG062-03	D04	HQP009776	NM_001570	IRAK2
QG062-03	D05	HQP009693	NM_002187	IL12B
QG062-03	D06	HQP009681	NM_001557	IL8RB
QG062-03	D07	HQP009679	NM_000634	IL8RA
QG062-03	D08	HQP009678	NM_000584	IL8
QG062-03	D09	HQP009556	NM_000041	APOE
QG062-03	D10	HQP009544	NM_000598	IGFBP3
QG062-03	D11	HQP009543	NM_182625	FLJ40869
QG062-03	D12	HQP009529	NM_000612	IGF2
QG062-03	E01	HQP009523	NM_000875	IGF1R
QG062-03	E02	HQP009518	NM_000618	IGF1
QG062-03	E03	HQP009201	NM_001039132	ICAM4
QG062-03	E04	HQP009139	NM_001643	APOA2
QG062-03	E05	HQP009085	NM_002155	HSPA6
QG062-03	E06	HQP009036	NM_005343	HRAS
QG062-03	E07	HQP009024	NM_000038	APC
QG062-03	E08	HQP008846	NM_001607	ACAA1
QG062-03	E09	HQP008808	NM_021155	CD209
QG062-03	E10	HQP008801	NM_001010931	HGF
QG062-03	E11	HQP008198	NM_173681	ATG9B
QG062-03	E12	HQP007767	NM_002075	GNB3
QG062-03	F01	HQP007722	NM_014905	GLS
QG062-03	F02	HQP007100	NM_015670	SENP3
QG062-03	F03	HQP006715	NM_012415	RAD54B
QG062-03	F04	HQP006427	NM_001039130	ALOX15B
QG062-03	F05	HQP006425	NM_001140	ALOX15
QG062-03	F06	HQP006424	NM_145659	IL27
QG062-03	F07	HQP006396	NM_153289	DEFB119
QG062-03	F08	HQP006375	NM_001629	ALOX5AP
QG062-03	F09	HQP006356	NM_000697	ALOX12
QG062-03	F10	HQP006296	NM_015364	LY96
QG062-03	F11	HQP006131	NM_014294	TRAM1
QG062-03	F12	HQP005959	NM_001002294	FMO3
QG062-03	G01	HQP005939	NM_002021	FMO1
QG062-03	G02	HQP005589	NM_012072	CD93
QG062-03	G03	HQP005403	NM_002006	FGF2
QG062-03	G04	HQP005242	NM_002002	FCER2
QG062-03	G05	HQP005240	NM_004106	FCER1G
QG062-03	G06	HQP005238	NM_002001	FCER1A
QG062-03	G07	HQP005100	NM_000691	ALDH3A1

Product Data Sheet

QG062-03	G08	HQP005058	NM_000130	F5
QG062-03	G09	HQP004984	NM_005236	ERCC4
QG062-03	G10	HQP004971	NM_001982	ERBB3
QG062-03	G11	HQP004948	NM_000120	EPHX1
QG062-03	G12	HQP004445	NM_000794	DRD1
QG062-03	H01	HGDC		
QG062-03	H02	HGDC		
QG062-03	H03	HQP006940	NM_002046	GAPDH
QG062-03	H04	HQP016381	NM_001101	ACTB
QG062-03	H05	HQP015171	NM_004048	B2M
QG062-03	H06	HQP006171	NM_012423	RPL13A
QG062-03	H07	HQP009026	NM_000194	HPRT1
QG062-03	H08	HQP054253	NR_003286	RN18S1
QG062-03	H09	RT		
QG062-03	H10	RT		
QG062-03	H11	PCR		
QG062-03	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Brain Cancer Gene qPCR Array (the Product). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products or deliver information obtained in service without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund limited to the actual purchase price.

GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2016 GeneCopoeia, Inc.

GeneCopoeia, Inc.
9620 Medical Center Drive, Suite 101
Rockville, MD 20850
+1 (301) 762-0888
+1 (866) 360-9531
inquiry@genecopoeia.com