

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_01	A01	HmiRQP2171	MIMAT0019064	hsa-miR-4525	
QM047_01	A02	HmiRQP0509	MIMAT0001631	hsa-miR-451a	
QM047_01	A03	HmiRQP0316	MIMAT0000076	hsa-miR-21-5p	
QM047_01	A04	HmiRQP0315	MIMAT0004494	hsa-miR-21-3p	
QM047_01	A05	HmiRQP0192	MIMAT0000437	hsa-miR-145-5p	
QM047_01	A06	HmiRQP0309	MIMAT0000241	hsa-miR-208a	
QM047_01	A07	HmiRQP0857	MIMAT0000689	hsa-miR-99b-5p	
QM047_01	A08	HmiRQP0184	MIMAT0000432	hsa-miR-141-3p	
QM047_01	A09	HmiRQP0002	MIMAT0000062	hsa-let-7a-5p	
QM047_01	A10	HmiRQP2458	MIMAT0019927	hsa-miR-4772-3p	
QM047_01	A11	HmiRQP0221	MIMAT0000646	hsa-miR-155-5p	
QM047_01	A12	HmiRQP4680	MIMAT0037306	hsa-miR-103a-1-5p	
QM047_01	B01	HmiRQP0157	MIMAT0004593	hsa-miR-130a-5p	
QM047_01	B02	HmiRQP0220	MIMAT0004658	hsa-miR-155-3p	
QM047_01	B03	HmiRQP0022	MIMAT0000101	hsa-miR-103a-3p	
QM047_01	B04	HmiRQP0156	MIMAT0000425	hsa-miR-130a-3p	
QM047_01	B05	HmiRQP0196	MIMAT0000449	hsa-miR-146a-5p	
QM047_01	B06	HmiRQP1298	MIMAT0009196	hsa-miR-103a-2-5p	
QM047_01	B07	HmiRQP0195	MIMAT0004608	hsa-miR-146a-3p	
QM047_01	B08	HmiRQP0023	MIMAT0007402	hsa-miR-103b	
QM047_01	B09	HmiRQP0191	MIMAT0004601	hsa-miR-145-3p	
QM047_01	B10	HmiRQP0263	MIMAT0000440	hsa-miR-191-5p	
QM047_01	B11	HmiRQP0264	MIMAT0001618	hsa-miR-191-3p	
QM047_01	B12	HmiRQP0283	MIMAT0000461	hsa-miR-195-5p	
QM047_01	C01	HmiRQP0282	MIMAT0004615	hsa-miR-195-3p	
QM047_01	C02	HmiRQP0341	MIMAT0004570	hsa-miR-223-5p	
QM047_01	C03	HmiRQP0342	MIMAT0000280	hsa-miR-223-3p	
QM047_01	C04	HmiRQP0393	MIMAT0000420	hsa-miR-30b-5p	
QM047_01	C05	HmiRQP0486	MIMAT0003329	hsa-miR-411-5p	
QM047_01	C06	HmiRQP0244	MIMAT0000261	hsa-miR-183-5p	
QM047_01	C07	HmiRQP0392	MIMAT0004589	hsa-miR-30b-3p	
QM047_01	C08	HmiRQP0243	MIMAT0004560	hsa-miR-183-3p	
QM047_01	C09	HmiRQP0396	MIMAT0000244	hsa-miR-30c-5p	
QM047_01	C10	HmiRQP0394	MIMAT0004674	hsa-miR-30c-1-3p	
QM047_01	C11	HmiRQP0395	MIMAT0004550	hsa-miR-30c-2-3p	
QM047_01	C12	HmiRQP0056	MIMAT0000421	hsa-miR-122-5p	
QM047_01	D01	HmiRQP0055	MIMAT0004590	hsa-miR-122-3p	
QM047_01	D02	HmiRQP0305	MIMAT0000264	hsa-miR-203a-3p	
QM047_01	D03	HmiRQP2343	MIMAT0019813	hsa-miR-203b-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_01	D04	HmiRQP2344	MIMAT0019814	hsa-miR-203b-3p	
QM047_01	D05	HmiRQP0337	MIMAT0004568	hsa-miR-221-5p	
QM047_01	D06	HmiRQP0338	MIMAT0000278	hsa-miR-221-3p	
QM047_01	D07	HmiRQP0340	MIMAT0004569	hsa-miR-222-5p	
QM047_01	D08	HmiRQP0339	MIMAT0000279	hsa-miR-222-3p	
QM047_01	D09	HmiRQP0347	MIMAT0004587	hsa-miR-23b-5p	
QM047_01	D10	HmiRQP0852	MIMAT0000095	hsa-miR-96-5p	
QM047_01	D11	HmiRQP0325	MIMAT0000274	hsa-miR-217-5p	
QM047_01	D12	HmiRQP0346	MIMAT0000418	hsa-miR-23b-3p	
QM047_01	E01	HmiRQP0289	MIMAT0000232	hsa-miR-199a-3p	
QM047_01	E02	HmiRQP0274	MIMAT0000222	hsa-miR-192-5p	
QM047_01	E03	HmiRQP0425	MIMAT0004701	hsa-miR-338-5p	
QM047_01	E04	HmiRQP0093	MIMAT0004602	hsa-miR-125a-3p	
QM047_01	E05	HmiRQP0495	MIMAT0003393	hsa-miR-425-5p	
QM047_01	E06	HmiRQP0398	MIMAT0000245	hsa-miR-30d-5p	
QM047_01	E07	HmiRQP0208	MIMAT0000450	hsa-miR-149-5p	
QM047_01	E08	HmiRQP4726	MIMAT0037308	hsa-miR-217-3p	
QM047_01	E09	HmiRQP0034	MIMAT0000254	hsa-miR-10b-5p	
QM047_01	E10	HmiRQP0275	MIMAT0004543	hsa-miR-192-3p	
QM047_01	E11	HmiRQP0496	MIMAT0001343	hsa-miR-425-3p	
QM047_01	E12	HmiRQP0008	MIMAT0004484	hsa-let-7d-3p	
QM047_01	F01	HmiRQP1574	MIMAT0016878	hsa-miR-4257	
QM047_01	F02	HmiRQP0207	MIMAT0004609	hsa-miR-149-3p	
QM047_01	F03	HmiRQP0427	MIMAT0000764	hsa-miR-339-5p	
QM047_01	F04	HmiRQP0790	MIMAT0004945	hsa-miR-744-5p	
QM047_01	F05	HmiRQP0428	MIMAT0004702	hsa-miR-339-3p	
QM047_01	F06	HmiRQP0789	MIMAT0004946	hsa-miR-744-3p	
QM047_01	F07	HmiRQP0391	MIMAT0000087	hsa-miR-30a-5p	
QM047_01	F08	HmiRQP0390	MIMAT0000088	hsa-miR-30a-3p	
QM047_01	F09	HmiRQP0183	MIMAT0004598	hsa-miR-141	
QM047_01	F10	HmiRQP0486	MIMAT0003329	hsa-miR-411	
QM047_01	F11	HmiRQP0523	MIMAT0002177	hsa-miR-486-5p	
QM047_01	F12	HmiRQP0832	MIMAT0000092	hsa-miR-92a-3p	
QM047_01	G01	HmiRQP0435	MIMAT0004694	hsa-miR-342-5p	
QM047_01	G02	HmiRQP0315	MIMAT0004494	hsa-miR-21	
QM047_01	G03	HmiRQP0225	MIMAT0000417	hsa-miR-15b-5p	
QM047_01	G04	HmiRQP0301	MIMAT0004657	hsa-miR-200c	
QM047_01	G05	HmiRQP0466	MIMAT0000728	hsa-miR-375	
QM047_01	G06	HmiRQP0099	MIMAT0000445	hsa-miR-126-3p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_01	G07	HmiRQP0313	MIMAT0004752	hsa-miR-20b	
QM047_01	G08	HmiRQP0399	MIMAT0000693	hsa-miR-30e-3p	
QM047_01	G09	HmiRQP0406	MIMAT0005792	hsa-miR-320b	
QM047_01	G10	HmiRQP0302	MIMAT0000617	hsa-miR-200c-3p	
QM047_01	G11	HmiRQP0446	MIMAT0000703	hsa-miR-361-5p	
QM047_01	G12	HmiRQP0290	MIMAT0000231	hsa-miR-199a-5p	
QM047_01	H01	NC	NC	NC	
QM047_01	H02	NC	NC	NC	
QM047_01	H03	hsnRNA U6	NR_002752	HK1	
QM047_01	H04	hsnoRNA U44	NR_002750	HK2	
QM047_01	H05	hsnoRNA U48	NR_002745	HK3	
QM047_01	H06	HsnoRNA U47	NR_002746	HK4	
QM047_01	H07	HsnoRNA U49	NR_002744	HK5	
QM047_01	H08	HsnRNA U68	NR_002450	HK6	
QM047_01	H09	RT	RT	RT	
QM047_01	H10	RT	RT	RT	
QM047_01	H11	PCR	PCR	PCR	
QM047_01	H12	PCR	PCR	PCR	
QM047_02	A01	HmiRQP3046	MIMAT0000250	hsa-miR-139-5p	
QM047_02	A02	HmiRQP0354	MIMAT0000082	hsa-miR-26a-5p	
QM047_02	A03	HmiRQP0007	MIMAT0000065	hsa-let-7d-5p	
QM047_02	A04	HmiRQP0343	MIMAT0000281	hsa-miR-224-5p	
QM047_02	A05	HmiRQP0308	MIMAT0000462	hsa-miR-206	
QM047_02	A06	HmiRQP0237	MIMAT0002821	hsa-miR-181d-5p	
QM047_02	A07	HmiRQP0782	MIMAT0004284	hsa-miR-675-5p	
QM047_02	A08	HmiRQP0837	MIMAT0000093	hsa-miR-93-5p	
QM047_02	A09	HmiRQP0078	MIMAT0005898	hsa-miR-1246	
QM047_02	A10	HmiRQP0291	MIMAT0000263	hsa-miR-199b-5p	
QM047_02	A11	HmiRQP1295	MIMAT0009198	hsa-miR-224-3p	
QM047_02	A12	HmiRQP1310	MIMAT0006790	hsa-miR-675-3p	
QM047_02	B01	HmiRQP0836	MIMAT0004509	hsa-miR-93-3p	
QM047_02	B02	HmiRQP2236	MIMAT0019704	hsa-miR-4644	
QM047_02	B03	HmiRQP0096	MIMAT0000423	hsa-miR-125b-5p	
QM047_02	B04	HmiRQP0357	MIMAT0000083	hsa-miR-26b-5p	
QM047_02	B05	HmiRQP1554	MIMAT0016858	hsa-miR-4306	
QM047_02	B06	HmiRQP0361	MIMAT0000419	hsa-miR-27b-3p	
QM047_02	B07	HmiRQP0301	MIMAT0004657	hsa-miR-200c-5p	
QM047_02	B08	HmiRQP0351	MIMAT0004498	hsa-miR-25-5p	
QM047_02	B09	HmiRQP0440	MIMAT0000255	hsa-miR-34a-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_02	B10	HmiRQP0314	MIMAT0001413	hsa-miR-20b-5p	
QM047_02	B11	HmiRQP0070	MIMAT0005591	hsa-miR-1236	
QM047_02	B12	HmiRQP4732	MIMAT0037313	hsa-miR-375-5p	
QM047_02	C01	HmiRQP0010	MIMAT0000066	hsa-let-7e-5p	
QM047_02	C02	HmiRQP0292	MIMAT0004490	hsa-miR-19a-5p	
QM047_02	C03	HmiRQP0630	MIMAT0004806	hsa-miR-548c-5p	
QM047_02	C04	HmiRQP3416	MIMAT0022727	hsa-miR-1307-5p	
QM047_02	C05	HmiRQP0434	MIMAT0004692	hsa-miR-340-5p	
QM047_02	C06	HmiRQP0400	MIMAT0000692	hsa-miR-30e-5p	
QM047_02	C07	HmiRQP0834	MIMAT0004792	hsa-miR-92b-5p	
QM047_02	C08	HmiRQP0359	MIMAT0000084	hsa-miR-27a-3p	
QM047_02	C09	HmiRQP0020	MIMAT0004513	hsa-miR-101-5p	
QM047_02	C10	HmiRQP0284	MIMAT0000226	hsa-miR-196a-5p	
QM047_02	C11	HmiRQP0854	MIMAT0000097	hsa-miR-99a-5p	
QM047_02	C12	HmiRQP0230	MIMAT0000070	hsa-miR-17-5p	
QM047_02	D01	HmiRQP0312	MIMAT0000075	hsa-miR-20a-5p	
QM047_02	D02	HmiRQP0352	MIMAT0000081	hsa-miR-25-3p	
QM047_02	D03	HmiRQP0439	MIMAT0004557	hsa-miR-34a-3p	
QM047_02	D04	HmiRQP0378	MIMAT0000688	hsa-miR-301a-3p	
QM047_02	D05	HmiRQP0095	MIMAT0004592	hsa-miR-125b-1-3p	
QM047_02	D06	HmiRQP0293	MIMAT0000073	hsa-miR-19a-3p	
QM047_02	D07	HmiRQP0286	MIMAT0001080	hsa-miR-196b-5p	
QM047_02	D08	HmiRQP0154	MIMAT0005951	hsa-miR-1307-3p	
QM047_02	D09	HmiRQP0021	MIMAT0000099	hsa-miR-101-3p	
QM047_02	D10	HmiRQP0415	MIMAT0000752	hsa-miR-328-3p	
QM047_02	D11	HmiRQP0442	MIMAT0000685	hsa-miR-34b-5p	
QM047_02	D12	HmiRQP0345	MIMAT0004496	hsa-miR-23a-5p	
QM047_02	E01	HmiRQP0097	MIMAT0004603	hsa-miR-125b-2-3p	
QM047_02	E02	HmiRQP0506	MIMAT0001545	hsa-miR-450a-5p	
QM047_02	E03	HmiRQP0831	MIMAT0004507	hsa-miR-92a-1-5p	
QM047_02	E04	HmiRQP0285	MIMAT0004562	hsa-miR-196a-3p	
QM047_02	E05	HmiRQP0168	MIMAT0000447	hsa-miR-134-5p	
QM047_02	E06	HmiRQP2147	MIMAT0019041	hsa-miR-4505	
QM047_02	E07	HmiRQP0441	MIMAT0004676	hsa-miR-34b-3p	
QM047_02	E08	HmiRQP0344	MIMAT0000078	hsa-miR-23a-3p	
QM047_02	E09	HmiRQP0615	MIMAT0002888	hsa-miR-532-5p	
QM047_02	E10	HmiRQP3060	MIMAT0026481	hsa-miR-134-3p	
QM047_02	E11	HmiRQP0247	MIMAT0000455	hsa-miR-185-5p	
QM047_02	E12	HmiRQP0444	MIMAT0000686	hsa-miR-34c-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_02	F01	HmiRQP0674	MIMAT0004795	hsa-miR-574-5p	
QM047_02	F02	HmiRQP0181	MIMAT0000431	hsa-miR-140-5p	
QM047_02	F03	HmiRQP0780	MIMAT0003880	hsa-miR-671-5p	
QM047_02	F04	HmiRQP2402	MIMAT0019871	hsa-miR-4741	
QM047_02	F05	HmiRQP0443	MIMAT0004677	hsa-miR-34c-3p	
QM047_02	F06	HmiRQP0673	MIMAT0003239	hsa-miR-574-3p	
QM047_02	F07	HmiRQP0006	MIMAT0000064	hsa-let-7c-5p	
QM047_02	F08	HmiRQP1333	MIMAT0010133	hsa-miR-2110	
QM047_02	F09	HmiRQP0456	MIMAT0000722	hsa-miR-370-3p	
QM047_02	F10	HmiRQP0363	MIMAT0004502	hsa-miR-28-3p	
QM047_02	F11	HmiRQP0279	MIMAT0004767	hsa-miR-193b-5p	
QM047_02	F12	HmiRQP0220	MIMAT0004658	hsa-miR-155	
QM047_02	G01	HmiRQP0003	MIMAT0004482	hsa-let-7b	
QM047_02	G02	HmiRQP0171	MIMAT0004698	hsa-miR-135b	
QM047_02	G03	HmiRQP0344	MIMAT0000078	hsa-miR-23a	
QM047_02	G04	HmiRQP0209	MIMAT0004610	hsa-miR-150	
QM047_02	G05	HmiRQP0211	MIMAT0000757	hsa-miR-151a	
QM047_02	G06	HmiRQP0254	MIMAT0002891	hsa-miR-18a	
QM047_02	G07	HmiRQP0203	MIMAT0004549	hsa-miR-148a	
QM047_02	G08	HmiRQP0844	MIMAT0004982	hsa-miR-939-5p	
QM047_02	G09	HmiRQP0161	MIMAT0000426	hsa-miR-132-3p	
QM047_02	G10	HmiRQP0358	MIMAT0004501	hsa-miR-27a	
QM047_02	G11	HmiRQP0844	MIMAT0004982	hsa-miR-939	
QM047_02	G12	HmiRQP0311	MIMAT0004493	hsa-miR-20a-3p	
QM047_02	H01	NC	NC	NC	
QM047_02	H02	NC	NC	NC	
QM047_02	H03	hsnRNA U6	NR_002752	HK1	
QM047_02	H04	hsnoRNA U44	NR_002750	HK2	
QM047_02	H05	hsnoRNA U48	NR_002745	HK3	
QM047_02	H06	HsnoRNA U47	NR_002746	HK4	
QM047_02	H07	HsnoRNA U49	NR_002744	HK5	
QM047_02	H08	HsnRNA U68	NR_002450	HK6	
QM047_02	H09	RT	RT	RT	
QM047_02	H10	RT	RT	RT	
QM047_02	H11	PCR	PCR	PCR	
QM047_02	H12	PCR	PCR	PCR	
QM047_03	A01	HmiRQP0364	MIMAT0000690	hsa-miR-296-5p	
QM047_03	A02	HmiRQP0295	MIMAT0000074	hsa-miR-19b-3p	
QM047_03	A03	HmiRQP0033	MIMAT0004556	hsa-miR-10b	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_03	A04	HmiRQP0276	MIMAT0000459	hsa-miR-193a	
QM047_03	A05	HmiRQP0016	MIMAT0004585	hsa-let-7i	
QM047_03	A06	HmiRQP0544	MIMAT0002871	hsa-miR-500a-3P	
QM047_03	A07	HmiRQP0229	MIMAT0000071	hsa-miR-17	
QM047_03	A08	HmiRQP0311	MIMAT0004493	hsa-miR-20a	
QM047_03	A09	HmiRQP0186	MIMAT0000434	hsa-miR-142-3p	
QM047_03	A10	HmiRQP0185	MIMAT0000433	hsa-miR-142-5p	
QM047_03	A11	HmiRQP0026	MIMAT0000103	hsa-miR-106a	
QM047_03	A12	HmiRQP0263	MIMAT0000440	hsa-miR-191	
QM047_03	B01	HmiRQP0125	MIMAT0000424	hsa-miR-128-3p	
QM047_03	B02	HmiRQP0228	MIMAT0004518	hsa-miR-16-2-3p	
QM047_03	B03	HmiRQP3040	MIMAT0022693	hsa-miR-204-3p	
QM047_03	B04	HmiRQP0330	MIMAT0000276	hsa-miR-219a-5p	
QM047_03	B05	HmiRQP0414	MIMAT0000756	hsa-miR-326	
QM047_03	B06	HmiRQP0519	MIMAT0002174	hsa-miR-484	
QM047_03	B07	HmiRQP2007	MIMAT0018204	hsa-miR-676-3p	
QM047_03	B08	HmiRQP3342	MIMAT0022717	hsa-miR-873-3p	
QM047_03	B09	HmiRQP0094	MIMAT0000443	hsa-miR-125a-5p	
QM047_03	B10	HmiRQP0139	MIMAT0004605	hsa-miR-129-2-3p	
QM047_03	B11	HmiRQP0160	MIMAT0004594	hsa-miR-132-5p	
QM047_03	B12	HmiRQP0190	MIMAT0000436	hsa-miR-144-3p	
QM047_03	C01	HmiRQP0227	MIMAT0000069	hsa-miR-16-5p	
QM047_03	C02	HmiRQP0306	MIMAT0000265	hsa-miR-204-5p	
QM047_03	C03	HmiRQP0332	MIMAT0000077	hsa-miR-22-3p	
QM047_03	C04	HmiRQP0501	MIMAT0002814	hsa-miR-432-5p	
QM047_03	C05	HmiRQP0784	MIMAT0004927	hsa-miR-708-3p	
QM047_03	C06	HmiRQP0804	MIMAT0004911	hsa-miR-874-3p	
QM047_03	C07	HmiRQP0025	MIMAT0000102	hsa-miR-105-5p	
QM047_03	C08	HmiRQP0137	MIMAT0000242	hsa-miR-129-5p	
QM047_03	C09	HmiRQP0174	MIMAT0004606	hsa-miR-136-3p	
QM047_03	C10	HmiRQP0189	MIMAT0004600	hsa-miR-144-5p	
QM047_03	C11	HmiRQP0232	MIMAT0000256	hsa-miR-181a-5p	
QM047_03	C12	HmiRQP0416	MIMAT0001629	hsa-miR-329-3p	
QM047_03	D01	HmiRQP0502	MIMAT0001627	hsa-miR-433-3p	
QM047_03	D02	HmiRQP0548	MIMAT0004775	hsa-miR-502-3p	
QM047_03	D03	HmiRQP0812	MIMAT0004947	hsa-miR-885-5p	
QM047_03	D04	HmiRQP0058	MIMAT0005458	hsa-miR-1224-5p	
QM047_03	D05	HmiRQP0145	MIMAT0005800	hsa-miR-1298-5p	
QM047_03	D06	HmiRQP0210	MIMAT0000451	hsa-miR-150-5p	

Plate	Position	Catalog No.	Mature miRNA Accession No.	Mature miRNA ID	miRNAs also Amplified
QM047_03	D07	HmiRQP0234	MIMAT0000257	hsa-miR-181b-5p	
QM047_03	D08	HmiRQP0327	MIMAT0000275	hsa-miR-218-5p	
QM047_03	D09	HmiRQP0417	MIMAT0000751	hsa-miR-330-3p	
QM047_03	D10	HmiRQP2076	MIMAT0018967	hsa-miR-4448	
QM047_03	D11	HmiRQP0622	MIMAT0004954	hsa-miR-543	
QM047_03	D12	HmiRQP0792	MIMAT0004957	hsa-miR-760	
QM047_03	E01	HmiRQP0824	MIMAT0000442	hsa-miR-9-3p	
QM047_03	E02	HmiRQP0074	MIMAT0000422	hsa-miR-124-3p	
QM047_03	E03	HmiRQP0098	MIMAT0000444	hsa-miR-126-5p	
QM047_03	E04	HmiRQP0182	MIMAT0004597	hsa-miR-140-3p	
QM047_03	E05	HmiRQP0223	MIMAT0000068	hsa-miR-15a-5p	
QM047_03	E06	HmiRQP0410	MIMAT0000755	hsa-miR-323a-3p	
QM047_03	E07	HmiRQP0418	MIMAT0004693	hsa-miR-330-5p	
QM047_03	E08	HmiRQP0767	MIMAT0003332	hsa-miR-656-3p	
QM047_03	E09	HmiRQP0796	MIMAT0003882	hsa-miR-767-5p	
QM047_03	E10	HmiRQP0835	MIMAT0003218	hsa-miR-92b-3p	
QM047_03	F01	NC	NC	NC	
QM047_03	F02	NC	NC	NC	
QM047_03	F03	hsnRNA U6	NR_002752	HK1	
QM047_03	F04	hsnoRNA U44	NR_002750	HK2	
QM047_03	F05	hsnoRNA U48	NR_002745	HK3	
QM047_03	F06	HsnoRNA U47	NR_002746	HK4	
QM047_03	F07	HsnoRNA U49	NR_002744	HK5	
QM047_03	F08	HsnRNA U68	NR_002450	HK6	
QM047_03	F09	RT	RT	RT	
QM047_03	F10	RT	RT	RT	
QM047_03	F11	PCR	PCR	PCR	
QM047_03	F12	PCR	PCR	PCR	