

Fusion transcript assays for *BCR-ABL* translocations

TaqMan Gene Expression Assays for fusion transcripts



HEMATOPATHOLOGY UNIT, HOSPITAL CLINIC, BARCELONA

Dolors Colomer, PhD, senior consultant responsible for the molecular biology area at the Hematopathology Unit and Spain Representative for the Molecular Monitoring EUTOS Program (middle), with technicians Sandra Cabezas (right) and Sandra Martinez (left)

APPLICATION

Fusion transcript expression

TECHNOLOGIES

TaqMan Gene Expression Assays for Fusion Transcripts

7900HT Fast Real-Time PCR System

Detecting fusion transcripts caused by chromosome translocations

Chromosome translocations are abnormalities caused by rearrangement of chromosome sections between nonhomologous chromosomes. These rearrangements may result in chimeric genes that express fusion transcripts. Some of these transcripts can be translated into fusion proteins that affect normal regulatory pathways and stimulate abnormal cell growth. A well-known example is the *BCR-ABL* chimeric mRNA (Philadelphia translocation), which is the result of a translocation of the *ABL* gene on chromosome 9 to the *BCR* gene breakpoint cluster on chromosome 22.

In this study, quantitative real-time PCR assays—Applied Biosystems™

TaqMan™ Gene Expression Assays for Fusion Transcripts, for detection of different *BCR-ABL* transcripts (targeting p210 and p190 isoforms)—were compared to primers and probes for the same targets recommended and standardized by the Europe Against Cancer (EAC) program and currently in widespread use [1–3]. The data presented here indicate that TaqMan Gene Expression Assays for Fusion Transcripts have greater sensitivity and use an easier, ready-to-go workflow. The standardized assay format and protocol with an optimized master mix results in less variability in assay setup and allows laboratories to generate more reproducible data.

The Philadelphia translocation (*BCR-ABL* fusion proteins)

BCR-ABL fusion proteins are

associated with the formation of the Philadelphia translocation (Ph) and are one of the most common genetic abnormalities studied in blood cancer research. At the molecular level, the Ph chromosome, or t(9;22) (q34;q11) translocation, results from the fusion of the *BCR* gene (chromosome 22), which forms the 5' end of the fusion transcript, to the *ABL* gene (chromosome 9), which forms the 3' end. In the vast majority of cases, the breakpoints in the *BCR* gene are found within three well-defined regions: the major breakpoint (*M-bcr*), minor

breakpoint (*m-bcr*), and micro breakpoint (μ -*bcr*). Depending on which breakpoints are used, three main chimeric proteins of different sizes are generated (Table 1, Figure 1). These BCR-ABL chimeric proteins (p190, p210, p230) show increased, deregulated tyrosine kinase activity, which appears to deregulate normal cytokine-dependent signal transduction leading to inhibition of apoptosis, independent of growth factors.

Real-time PCR detects translocations and quantifies expression

Current methods for identifying translocations include FISH and karyotyping, neither of which can be used to quantify the expression level of the fused gene as real-time PCR does. Real-time quantitative PCR can provide an appropriate monitoring strategy for analyzing *BCR-ABL* expression levels in the samples under study [5].

Improved *BCR-ABL* PCR assays

Real-time PCR is the gold standard for quantitative measurement of nucleic acid. In collaboration with Thermo Fisher Scientific, EAC researchers developed primers and probes to detect specific *BCR-ABL*

fusion transcripts [1,2]. Recently we have improved on these primer and probe designs, creating new TaqMan Gene Expression Assays for all of the *BCR-ABL* fusion transcripts (Table 3). Selected transcripts were annotated, and bases located at the fusion transcript breakpoint, known SNPs, and repetitive sequences were masked. TaqMan™ minor groove binder (MGB) Assays were then designed using the Applied Biosystems™ bioinformatics design pipeline. The assays were designed such that the primers and probes would bind on either side of the fusion transcript breakpoint (Figure 2), and each assay design was checked by *in silico* quality control.

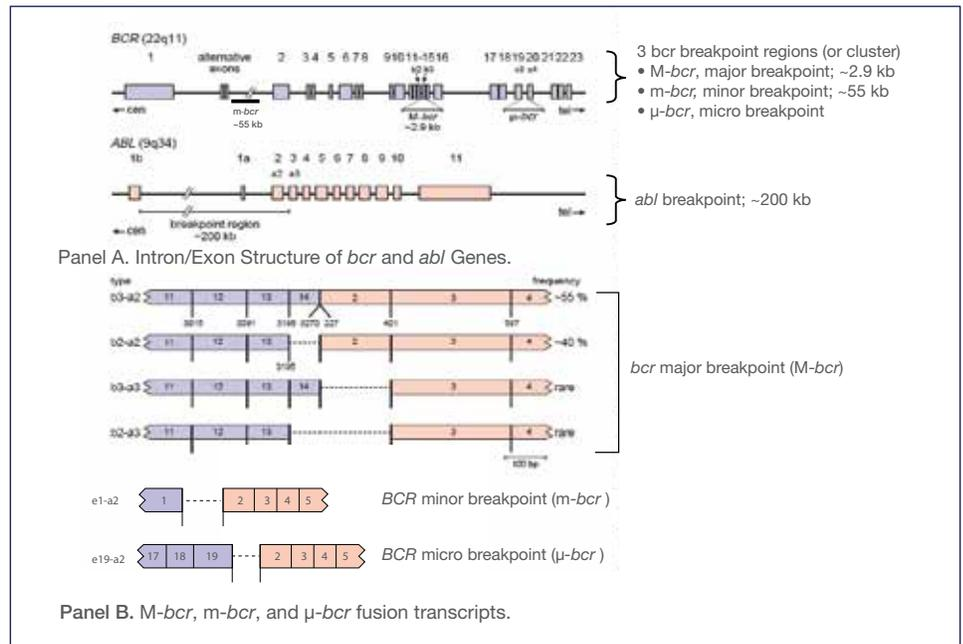


Figure 1. *BCR-ABL* chromosomal breakpoints and fusion gene transcripts. (A) Schematic diagram of the exon/intron structure of the *BCR* and *ABL* genes involved in t(9;22) (q34;q11). The centromere (cen) and telomere (tel) orientation, exon numbering, and relevant breakpoint regions are indicated, including the micro breakpoint cluster region (μ -*bcr*). (B) Schematic diagram of *BCR-ABL* major (M), minor (m), and micro (μ) transcripts. For M-*bcr*, the b3-a2 and b2-a2 transcripts are found most frequently, but sporadic cases with b3-a3 and b2-a3 transcripts have been reported. (Parts of this figure are used with permission from *Leukemia*.)

Table 1. *BCR-ABL* fusion transcripts and resulting fusion proteins.

Breakpoint designation	Chrm 22 (<i>BCR</i> gene) break location	Chrm 9 (<i>ABL</i> gene) break location	Variant transcript designation	Chimeric protein size (name)
M- <i>bcr</i> (exons 12-16)	Intron 13	Intron 1	b2-a2 (e13-a2)	210 kDa (p210)
	Intron 14	Intron 1	b3-a2 (e14-a2)	" "
	Intron 13	Intron 2	b2-a3 (rare) (e13-a3)	" "
	Intron 14	Intron 2	b3-a3 (rare) (e14-a3)	" "
m- <i>bcr</i>	Intron 1	Intron 1	e1-a2	190 kDa (p190)
μ - <i>bcr</i> (rare)	Intron 19	Intron 1	e19-a2	230 kDa (p230)

TaqMan Gene Expression Assays vs. EAC assays

As proof of principle, the TaqMan Gene Expression Assay designs were first tested using plasmids containing the translocation variant and human samples containing the translocation event. Amplification only occurred in samples containing the fusion transcript, confirming assay specificity (data not shown).

Subsequently, researchers in the Hematopathology Unit of the Hospital Clinic in Barcelona used human samples to compare the TaqMan Gene Expression Assays for

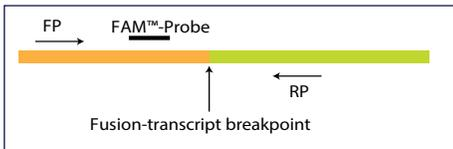


Figure 2. Design for TaqMan Gene Expression Assays for fusion transcripts. Design of primer and probe binding locations [2].

Approximately 10 bp surrounding the breakpoint were masked to avoid designing the probe across this region, since the precise sequence around the breakpoint can be ambiguous. Thus, probes did not span transcript breakpoints. FP = forward primer; RP = reverse primer.

BCR-ABL fusion transcripts and the EAC primer and probe designs [1,2]. (Note: The *BCR-ABL* TaqMan Gene Expression Assays include assays for several fusion transcripts for which there were no corresponding EAC designs.) The experimental procedure is provided in the sidebar, *Technical details for fusion assays*. Results follow.

Detection of *BCR-ABL* fusion transcript

Although it has been recommended to use a fixed threshold value using EAC designs (threshold 0.1), better results have been obtained using TaqMan Gene Expression Assays with Automatic Analysis provided by the Applied Biosystems™ 7900HT Fast Real-Time PCR System. Standard curves generated by amplification of dilutions of each of the fusion transcripts (cloned into Ipsogen and Invitrogen™ pCR2.1 TOPO™ TA plasmid vectors) using both TaqMan Gene Expression Assays and the *ABL* endogenous control TaqMan Assay were very reproducible. PCR efficiencies were close to 100% with $R^2 > 0.99$ (Figure 3).

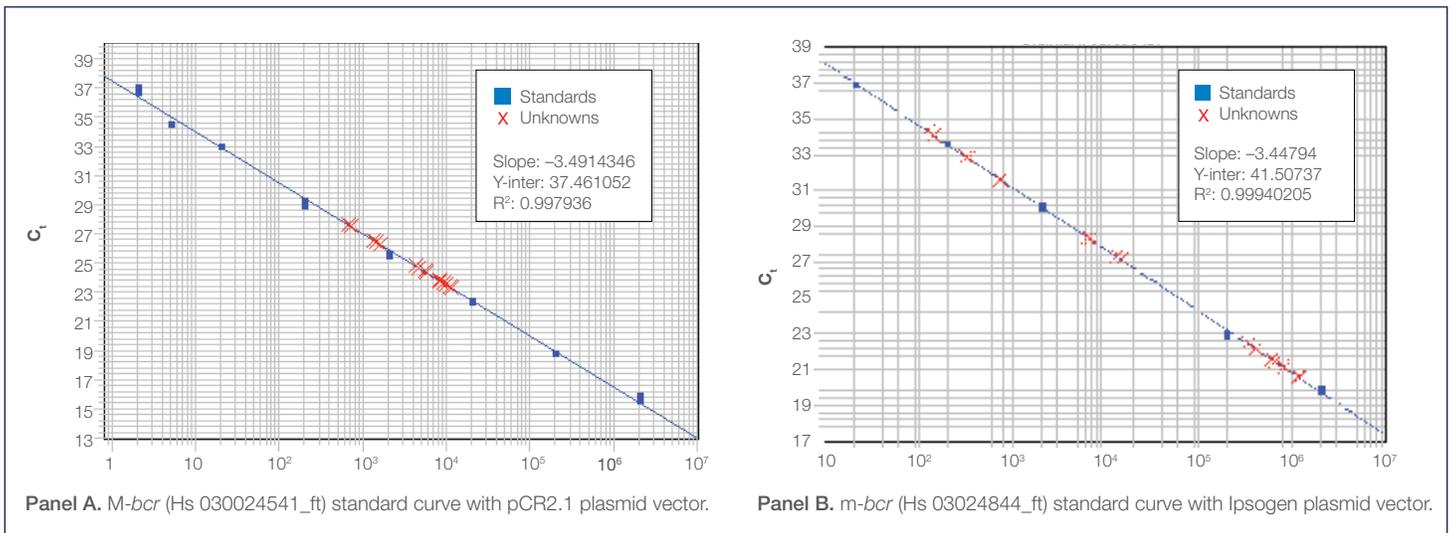


Figure 3. Standard curves obtained using TaqMan Gene Expression Assays. (A) Two replicates of each of eight dilutions (2×10^6 , 2×10^5 , 2×10^4 , 2×10^3 , 2×10^2 , 2×10 , 5, and 2 copies/ μL) of the *M-bcr* breakpoint cloned into the pCR2.1 plasmid vector were amplified, using 2 μL of each dilution. (B) Two replicates of each of five dilutions (2×10^6 , 2×10^5 , 2×10^3 , 2×10^2 , and 2×10 copies/ μL) of the *m-bcr* breakpoint cloned into the Ipsogen plasmid vector were amplified using 2 μL of each dilution.

M-*bcr* and m-*bcr* analysis

M-*bcr* and m-*bcr* fusion transcript quantification results were generated using the Ipsogen plasmid vector and the pCR2.1 plasmid vector (this second vector has only been used for M-*bcr*). A statistical study demonstrated that quantitative results obtained for M-*bcr* and m-*bcr* fusion transcripts with TaqMan Gene Expression Assays for Fusion Transcripts were identical to those obtained with EAC assays for all 20 samples analyzed, independent of the plasmid vector used (Ipsogen or pCR2.1 plasmid vector) (Figure 4 and Figure 5).

To check the sensitivity of TaqMan Gene Expression Assays compared with the EAC assays, samples amplified with TaqMan Gene Expression Assays were reanalyzed using a fixed threshold value of 0.1. For M-*bcr*, the C_t values obtained with TaqMan Gene Expression Assays were 0.49 to 1.97 lower (mean: 0.85) than C_t values obtained using EAC assays (Figure 6, Panel A). For m-*bcr*, the C_t values obtained with TaqMan Gene Expression Assays were 1.76 to 2.53 (mean: 2.22) lower than C_t values obtained using the EAC assays (Figure 6, Panel B). This analysis demonstrates the greater sensitivity of TaqMan Gene Expression Assays

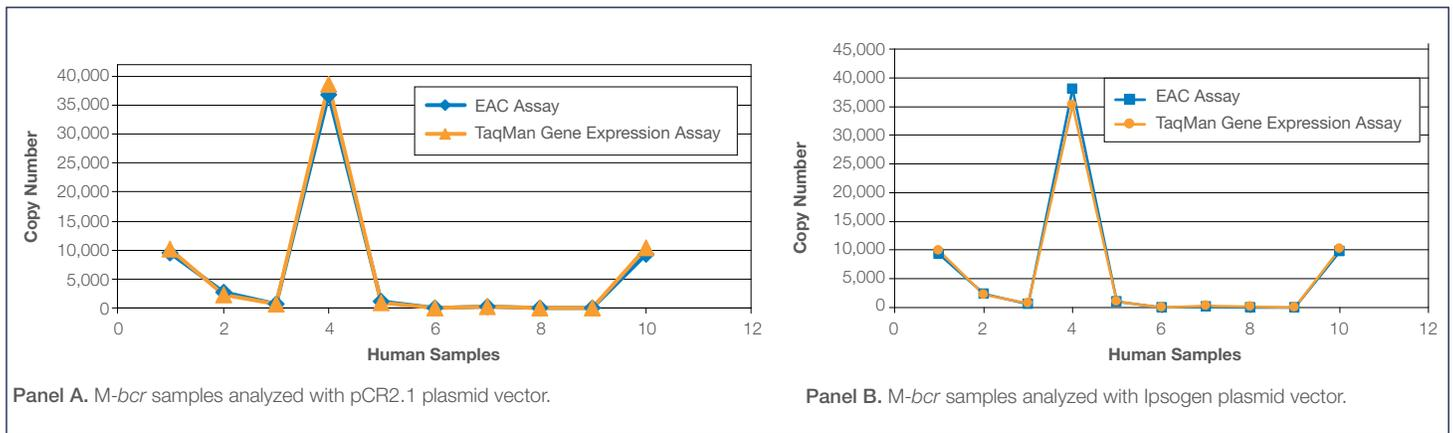


Figure 4. M-*bcr* copy number using TaqMan Gene Expression Assays vs. EAC assay. Copy number analysis was performed using TaqMan Gene Expression Assay (Hs03024541_ft) and the comparable EAC assay with either (A) pCR2.1 plasmid vector or (B) Ipsogen plasmid vector. The TaqMan Gene Expression Assay gave the same results as the EAC assay in both cases.

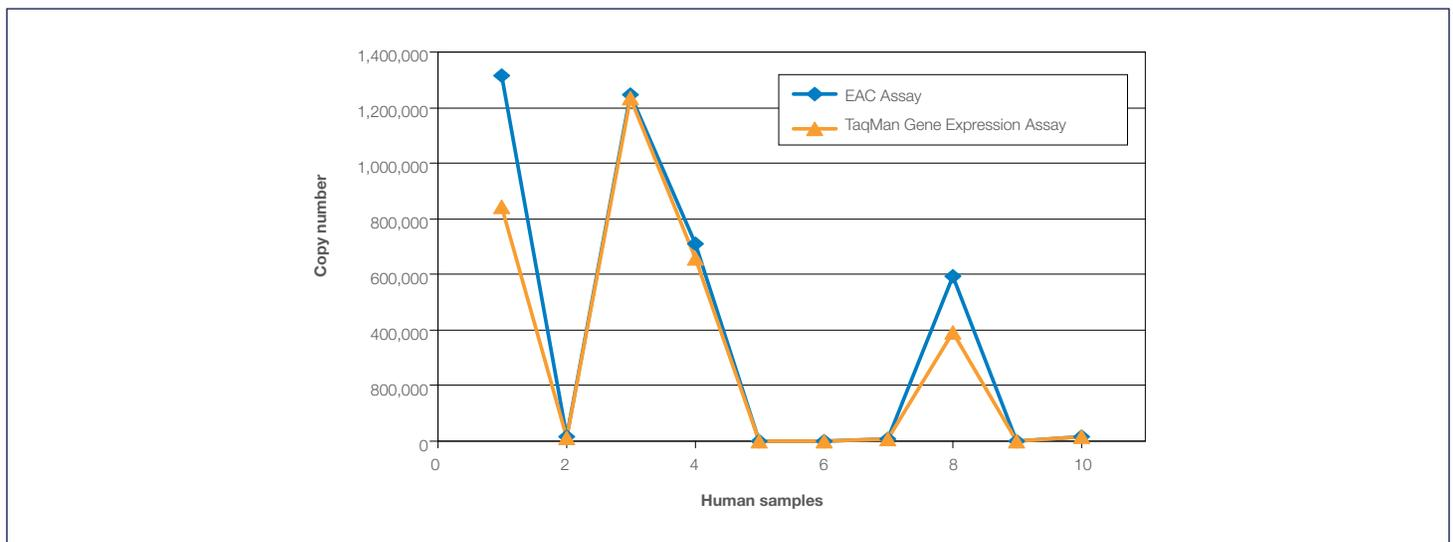
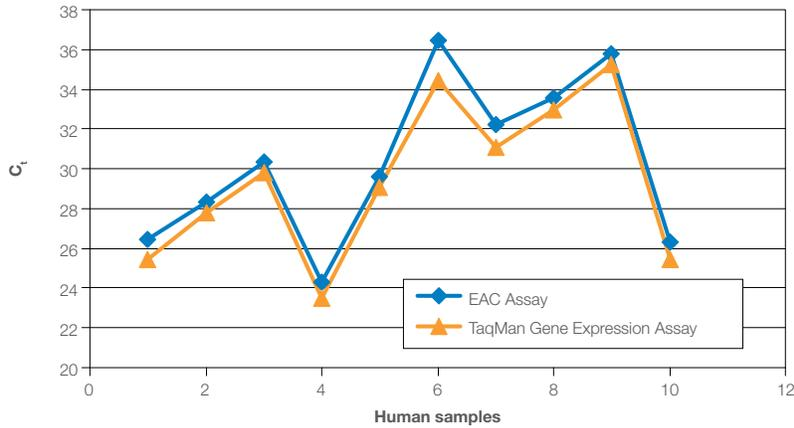


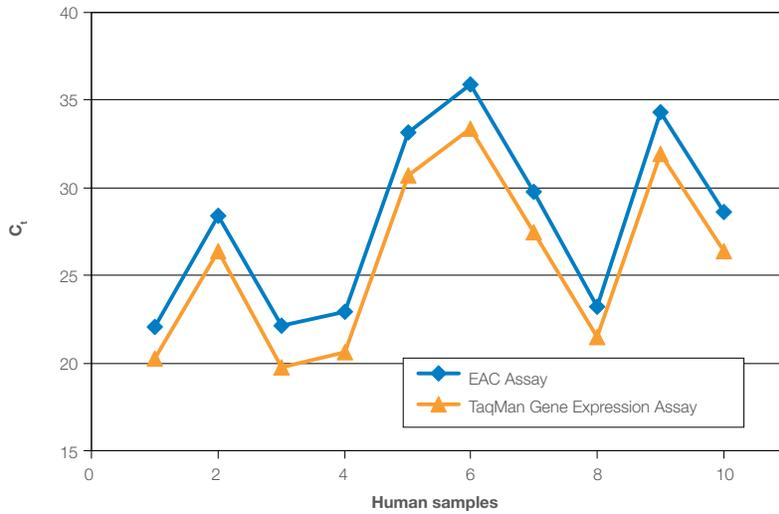
Figure 5. m-*bcr* copy number using TaqMan Gene Expression Assay vs. EAC assay. Copy number analysis was performed using TaqMan Gene Expression Assay Hs03024844_ft and the comparable EAC assay with the Ipsogen plasmid vector. The TaqMan Gene Expression Assay gave the same results as the EAC assay.

versus the EAC assays. Besides, quantitative results obtained with the fixed threshold value for both *M-bcr* and *m-bcr* translocations were almost identical to those obtained with the 7900HT Fast Real-Time PCR System's Automatic Analysis. The greater sensitivity

achieved with TaqMan Gene Expression Assays has also made possible the detection of the *m-bcr* fusion transcript in a sample that previously could not be amplified using the EAC probe and primers (data not shown).



Panel A. *M-bcr* samples.



Panel B. *m-bcr* samples.

Figure 6. C_t values for *M-bcr* using TaqMan Gene Expression Assays vs. EAC assays. Reanalysis of *M-bcr* samples (A) TaqMan Gene Expression Assay Hs03024541_ft and comparable EAC assay using a fixed 0.1 threshold value) and of *m-bcr* samples (B) TaqMan Gene Expression Assay Hs03024844_ft and comparable EAC assay). The TaqMan Gene Expression Assays were consistently more sensitive than the EAC assays.

Technical details for fusion assays

Samples. Peripheral blood samples were a subset of routine samples collected by the Hematopathology Unit of the hospital, and were kept anonymous. Informed consent was obtained in accordance with the Institutional Ethics Committee of the Hospital Clinic (Barcelona, Spain) and the Helsinki declaration.

RNA and cDNA preparation. Dr. Colomer and colleagues used 10 human samples each for analysis of *M-bcr* and *m-bcr* expression. Leukocytes from *BCR-ABL*-positive peripheral blood samples were isolated by 2% dextran sedimentation. Total RNA was extracted from the leukocytes using Invitrogen™ TRIzol™ Reagent following the manufacturer's instructions. Total RNA (1 µg; quantified by Thermo Scientific™ NanoDrop™ technology) was reverse transcribed into cDNA (50 µL reactions) using random primers and Invitrogen™ M-MLV Reverse Transcriptase (or Invitrogen™ SuperScript™ I or II) following the protocol published by the EAC Consortium [2] (Table 2).

Fusion transcript assays. *M-bcr* (b2-a2, b3-a2; TaqMan Gene Expression Assay Hs030024541_ft) and *m-bcr* (e1-a2; TaqMan Gene Expression Assay Hs03024844_ft) fusion transcript quantification was performed in two different reactions using both TaqMan Gene Expression Assays and the EAC fusion transcript primers and probes. The Abelson (*ABL*) TaqMan Gene Expression Assay (Hs99999002_mH) and EAC *ABL* primers and probes were used together to amplify the endogenous control.

Quantitative real-time PCR (qPCR) for *BCR-ABL* and *ABL* control transcripts was performed in duplicate on the 7900HT Fast Real-Time PCR System using standard run conditions. Reactions (25 µL) included Applied Biosystems™ TaqMan™ Gene Expression Master Mix and cDNA (2 µL). A known positive and negative control were amplified for each assay.

Table 2. EAC reverse transcription protocol.

Incubate 1 µg total RNA in 10 µL H ₂ O at 70°C for 10 min
Cool on ice and add other reagents to a final volume of 20 µL
100 U Reverse transcriptase (either M-MLV or SuperScript I or II)
RT buffer (according to the reverse transcriptase used)
1 mM dNTP
10 mM DTT
25 µM Random hexamers
20 U RNase inhibitor
Incubate subsequently at:
Room temperature for 10 min
42°C for 45 min
99°C for 3 min
Place the sample at 4°C
Dilute the final cDNA with 30 µL of H ₂ O

Data analysis. For analysis of both the *M-bcr* breakpoint and *ABL* endogenous control gene data, standard curves were created using the Ipsogen plasmid vector and/or the pCR2.1 + *BRC-ABL* plasmid vector [3]. For analysis of the *m-bcr* breakpoint data, standard curves were created using the Ipsogen plasmid vector. Detection of the rare *M-bcr* b3-a3 transcript was tested with a TaqMan Gene Expression Assay (Hs 03043652_ft) on *M-bcr* b3-a3-positive samples using the protocol described above. No plasmid was available for this last fusion transcript.

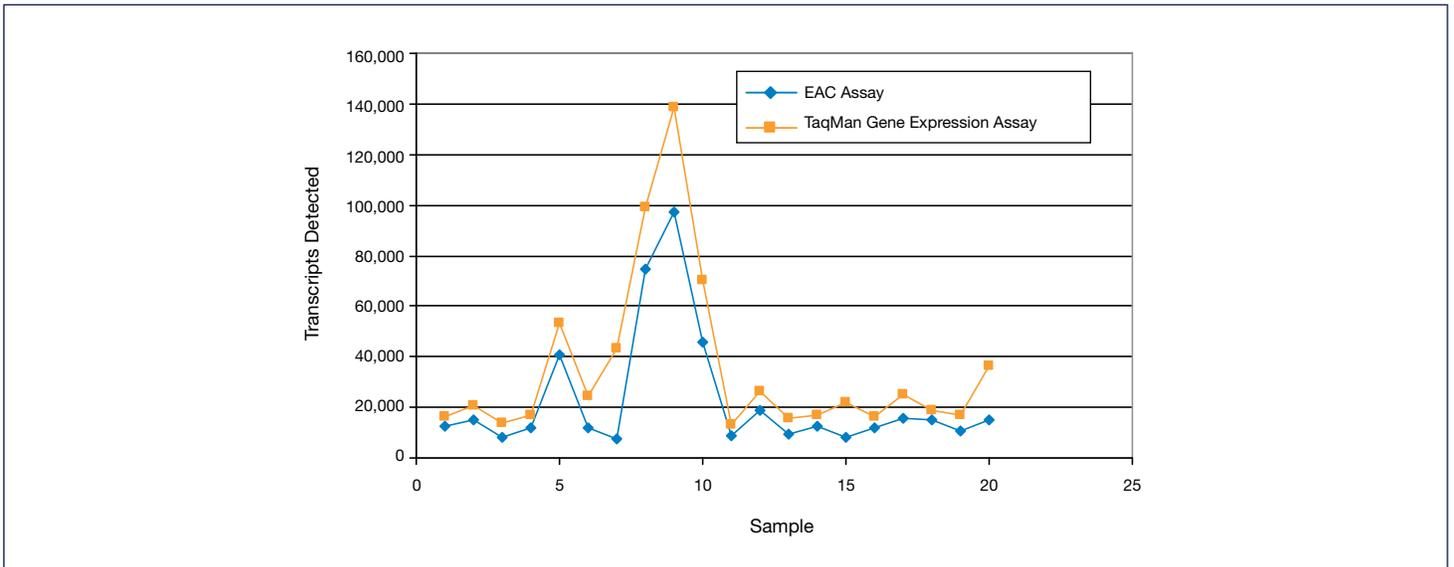


Figure 7. Better detection of *ABL* expression levels using TaqMan Gene Expression Assays. Samples (20) were amplified using probe and primer designs for the *ABL* endogenous control transcript. Higher *ABL* expression levels were detected with the TaqMan Gene Expression Assay Hs99999002_mH than with the EAC designs (pCR2.1 plasmid vector). The same data was obtained using the Ipsogen plasmid (data not shown).

Analysis of *ABL* expression

The expression levels of the *ABL* endogenous control obtained with TaqMan Assay Hs99999002_mH were higher than those obtained using the EAC designs in all 20 samples (on average 1.7 times higher). Samples were analyzed independently with the Ipsogen and the pCR2.1 plasmid vector (Figure 7).

Analysis of rare fusion transcript forms

A TaqMan Gene Expression Assay (Hs03024652_ft) was able to detect the rare *M-bcr* b3-a3 transcript in human *M-bcr* b3-a3-positive samples. Figure 8 shows a time course of *M-bcr* b3-a3 transcript expression in the positive samples taken at time points out to one year, when the rare transcript was not longer detected. EAC designs were not available for detection of this transcript.

Ready-to-use TaqMan Assays provide many advantages for detecting fusion transcripts

We have recently released a novel set of 165 TaqMan Gene Expression Assays for quantitation of human fusion transcripts. These assays were developed with our validated bioinformatics pipeline used to design the 1.3 million TaqMan Gene Expression Assays currently available. As with other TaqMan Gene Expression Assays, the fusion transcript assays undergo a synthesis quality control test using mass spectrometry to verify primer and probe sequence and concentration.

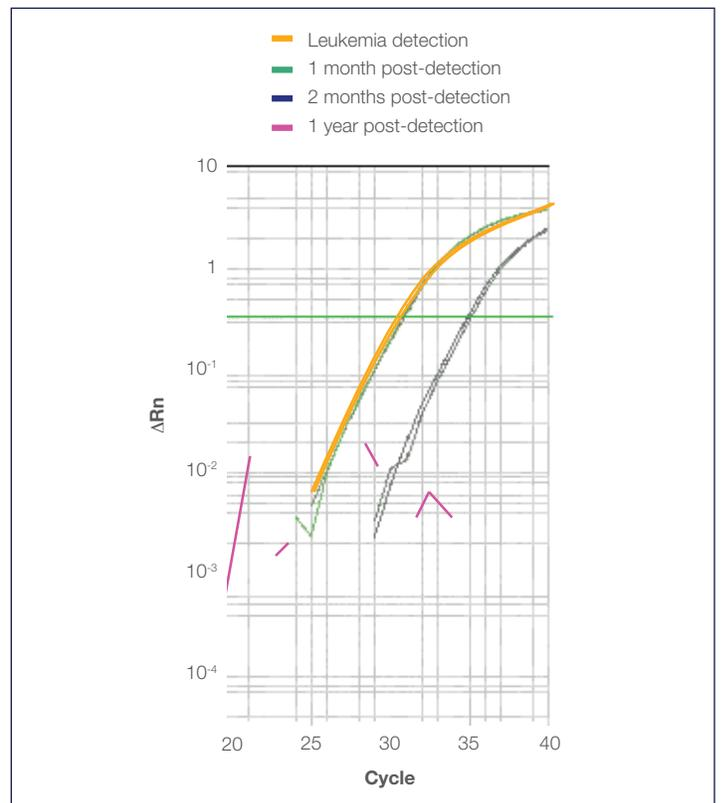


Figure 8. TaqMan Assays make possible detection of rare *M-bcr* transcript. Detection of the rare b3-a3 *M-bcr* translocation in a positive sample, over a time course out to 1 year. Amplifications were performed using the TaqMan Gene Expression Assay Hs03024652_ft.

Simplifying fusion transcript detection and quantitation with better tools

TaqMan Gene Expression Assay products enable researchers to conduct fusion studies quickly and easily by eliminating the time-consuming processes involved in assay development. As all of the assays are ready to use (primer and probes formulated in single-tube, 20X mix), it is easier to set up the reaction and to compare results with other researchers and labs directly and accurately. See Table 3 and page 8 for ordering information.

You can search online for the assay you need at thermofisher.com/taqmanfusion

Table 3. *BCR-ABL* and *ABL* endogenous control TaqMan Gene Expression Assays.

BCR-ABL TaqMan Gene Expression Assays		
Assay ID	Transcript	Assay Accession No.
Hs03024541_ft	b2-a2	AJ131467.1
Hs03024541_ft	b3-a2	AJ131466.1
Hs03024844_ft	e1-a2	AF113911.1
Hs03205538_ft	e19-a2	AM491363.1
Hs03043652_ft	b3-a3	AM491360.1
Hs03043652_ft	b3-a2	AJ131466.1
Hs03024646_ft	b2-a3	AY043457.1
Hs03024646_ft	b3-a3	AM491360.1
Hs03024646_ft	b2-a2	AJ131467.1
Endogenous Control		
Gene name	Assay ID	
Abelson (<i>ABL</i>)	Hs 99999002_mH	

In this study, Dr. Colomer and colleagues use TaqMan Gene Expression Assays for *M-bcr* (Hs030024541_ft) and *m-bcr* (Hs03024844_ft) to detect and quantify *BCR-ABL* fusion transcripts. The TaqMan Assays were able to detect *M-bcr* and *m-bcr* transcripts in the same samples as EAC probe and primers designed to these targets. However, the TaqMan Gene Expression Assays provided more sensitivity, yielding transcript

amplification in samples not amplifiable with the EAC designs. In addition, the single-tube 20X format of TaqMan Gene Expression Assays and associated Applied Biosystems reagents made the assays easier to process, and saved valuable time. “When used with the EAC program’s standardized protocol, these TaqMan Gene Expression Assays would help eliminate much of the variability seen across different laboratories due

to individual primer and probe preparation protocols. Having standardized reverse transcription quantitative PCR (RT-qPCR) assays would harmonize the current technology for detecting *BCR-ABL* transcripts, saving time and providing more reproducibility in results,” notes Dr. Colomer.

Slightly better results have been achieved using the automatic threshold software on the 7900HT Fast Real-Time PCR System to analyze results obtained with TaqMan Gene Expression Assays. However, as it is recommended to use a fixed threshold value with the EAC primer and probe designs (threshold = 0.1), the data for all samples was reanalyzed using this fixed threshold. Under these conditions, TaqMan Gene Expression Assays were shown to be more sensitive, allowing the amplification of a sample that was negative using EAC primer and probe designs.

The TaqMan Assays were also able to successfully detect the rare *M-bcr* transcript b3-a3 and the μ -*bcr* transcript e19-a2 in human samples (data not shown).

It is especially important to note that these particular assays provide new tools to researchers, since there are no comparable EAC probe and primer designs for these specific translocations.

We provide researchers a standardized, easy-to-use workflow to quantify the different *BCR-ABL* fusion transcripts. This workflow makes it possible to obtain rapid and reproducible results within and between laboratories.

TaqMan Gene Expression Master Mix

The TaqMan Gene Expression Assay workflow can be further simplified by incorporating TaqMan Gene Expression Master Mix. The Gene Expression Master Mix comes concentrated with all needed reagents premixed, decreasing hands-on time for dilution, mixing, and pipetting. TaqMan Gene Expression Master Mix delivers sensitive and specific detection across a broad range of template quantities, down to a single copy of target. For ease of use, TaqMan Gene Expression Master Mix uses universal thermal cycling conditions and users can set up the reaction at room temperature. See page 8 for ordering information.

LeukoLOCK™ Total RNA Isolation System

The Ambion™ LeukoLOCK Total RNA Isolation System is an innovative method for cellular fractionation of whole blood, and total RNA stabilization and extraction from the leukocyte population. It has been optimized for use with human blood. Blood is a storehouse of cellular information; however, the presence of globin mRNA in RNA prepared from whole blood can interfere with downstream expression profiling applications. The LeukoLOCK system employs filter-based leukocyte-depletion technology to isolate leukocytes from whole blood, and Ambion™ *RNAlater*™ to stabilize the cells on the filter. By excluding red blood cells, the RNA that is purified from captured leukocytes is inherently depleted of globin mRNA, which improves sample utility for expression profiling and other applications. See page 8 for ordering information.

High-Capacity cDNA Reverse Transcription Kit

The Applied Biosystems™ High-Capacity™ cDNA Reverse Transcription Kit delivers extremely high-quality, single-stranded cDNA from total RNA. It contains all components necessary for the quantitative conversion of 0.02 to 2 µg total RNA in a single 20 µL reaction to single-stranded cDNA. Downstream applications include real-time PCR, standard PCR, and microarrays. See page 8 for ordering information.

Custom Plating Service

The Applied Biosystems™ TaqMan™ Custom Plating Service offers the convenience of pre-plated TaqMan Gene Expression Assays, Custom Assays, and Custom Probe/Primer Sets in 96- or 384-well plates. Set up custom configurations using TaqMan Gene Expression Assays (Inventoried, Made-to-Order, and Custom) and Custom TaqMan™ Probes and Primers. You can select from a variety of reaction volumes and receive assays in dried or liquid formulation.

Acknowledgments

We acknowledge Beatriz Cabot, Applied Biosystems Senior Field Application Specialist, Molecular Biology, and Marisa Checa, Applied Biosystems Field Application Specialist, Molecular Biology, for their generous contribution in this collaboration.

References

1. p190 primers and probe: ENF401, ENR561, ENF541; p210 primers and probe: ENF501, ENR561, ENP541.
2. Gabert J, Beillard E, van der Velden VH et al. (2003) Standardization and quality control studies of 'real-time' quantitative reverse transcriptase polymerase chain reaction of fusion gene transcripts for residual disease detection in leukemia—a Europe Against Cancer program. *Leukemia* 17:2318–2357.
3. Beillard E, Pallisgaard N, van der Velden VH et al. (2003) Evaluation of candidate control genes for diagnosis and residual disease detection in leukemic patients using 'real-time' quantitative reverse-transcriptase polymerase chain reaction (RQ-PCR)—a Europe Against Cancer program. *Leukemia* 17:2474–2486.
4. Hughes T, Deininger M, Hochhaus A et al. (2006) Monitoring CML patients responding to treatment with tyrosine kinase inhibitors: review and recommendations for "harmonizing" current methodology for detecting BCR-ABL transcripts and kinase domain mutations and for expressing results. *Blood* 108:28–37.
5. Melo JV (1996) The diversity of BCR-ABL fusion proteins and their relationship to leukemia phenotype. *Blood* 88:2375–2384.

Ordering information

Description	Size	Cat. No.
LeukoLOCK Total RNA Isolation System	20 rxn*	AM1923
High-Capacity cDNA Reverse Transcription Kit	200 rxn*	4368814
TaqMan Gene Expression Assays for Fusion Transcripts		
Inventoried	250 rxn	4331182 [†]
Made-to-order	360 rxn	4351372
TaqMan Gene Expression Master Mix, 1 Mini-Pack (1 x 1 mL)	40 rxn*	4370048
TaqMan Universal PCR Master Mix, 1-Pack (1 x 5 mL)	200 rxn*	4304437
7900HT Fast Real-Time PCR System with Standard 96-well Block Module	1 instrument	4329003

*Available in other sizes or in bundles.

[†]See list below for specific assays.

TaqMan Gene Expression Assays for *BCR-ABL* fusion transcripts

Assay ID	Transcript	Assay Accession No.
Hs03024541_ft	b2-a2	AJ131467.1
Hs03024541_ft	b3-a2	AJ131466.1
Hs03024844_ft	e1-a2	AF113911.1
Hs03205538_ft	e19-a2	AM491363.1
Hs03043652_ft	b3-a3	AM491360.1
Hs03043652_ft	b3-a2	AJ131466.1

Endogenous control

Gene name	Assay ID
Abelson (<i>ABL</i>)	Hs99999002_mH

Find out more at thermofisher.com/taqmanfusion

A license to perform the patented 5' Nuclease Process for research is obtained by the purchase of (i) both Licensed Probe and Authorized 5' Nuclease Core Kit, (ii) a Licensed 5' Nuclease Kit, or (iii) license rights from Applied Biosystems. The TaqMan Gene Expression Assay contains Licensed Probe. Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,538,848, 5,723,591, 5,876,930, 6,030,787, 6,258,569, and 5,804,375 (claims 1–12 only). The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. Separate purchase of an Authorized 5' Nuclease Core Kit would convey rights under the applicable claims of US Patents Nos. 5,210,015 and 5,487,972, and corresponding patent claims outside the United States, which claim 5' nuclease methods. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel.

For Research Use Only. Not for use in diagnostic procedures. © 2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license. **CO016435 0715**

ThermoFisher
SCIENTIFIC