

MYCN
TP53

Real-Time qPCR as a Tool for Evaluating RNAi-Mediated Gene Silencing

Authors: Tom Van Maerken, Pieter Mestdagh,
Sarah De Clercq, Filip Pattyn, Nurten Yigit,
Anne De Paepe, Jean-Christophe Marine,
Frank Speleman, and Jo Vandesompele

Real-time quantitative PCR (rt-qPCR) is the method of choice for accurate, sensitive, and specific quantitation of nucleic acid sequences. Applications of this technology are numerous, both in molecular diagnostics and in virtually all fields of life sciences, including gene expression profiling, measurement of DNA copy number alterations, genotyping, mutation detection, pathogen detection, measurement of viral load, disease monitoring, and assessment of drug response. Several ingredients are essential to the successful and reliable completion of an rt-qPCR assay, such as careful primer design and evaluation, template preparation, the use of a robust normalization strategy, and accurate data analysis. This article describes how rt-qPCR can be implemented as a tool to monitor silencing efficiency and functional effects of RNA interference (RNAi)-mediated gene knockdown, using examples from our research on neuroblastoma. For detailed information on the experiments that contributed to this research, including instruments, reagents, and procedures, request bulletin 5692.

BIO-RAD

Neuroblastoma and the *MYCN* and *TP53* Cancer Genes

Neuroblastoma is a childhood cancer derived from precursor cells of the adrenergic system, arising in the adrenal medulla or in sympathetic ganglia. Although a relatively rare form of cancer, neuroblastoma is among the most fatal of childhood diseases. Indicators of mortality include age at diagnosis (the outcome for children with neuroblastoma is most favorable when diagnosed before the age of one year, even when the disease has metastasized), tumor stage, and level of *MYCN* protein activity (the most fatal clinicogenetic subtype of neuroblastoma is characterized by amplification of the *MYCN* oncogenic transcription factor) (Vandesompele et al. 2005). The mechanisms by which this transcription factor exerts its oncogenic activity and confers an unfavorable prognosis are poorly understood.

Another intriguing feature of neuroblastoma is the remarkably low frequency of *TP53* mutations at diagnosis (Tweddle et al. 2003). Previous studies have shown that reactivation of the p53 pathway by the selective small-molecule MDM2 antagonist nutlin-3 constitutes a promising novel therapeutic approach for neuroblastoma (Van Maerken et al. 2006). To gain insight into the mechanism of action of these two pivotal genes in neuroblastoma pathogenesis and to create model systems for future exploration of targeted therapeutics in relationship to *MYCN* and *TP53* status, RNAi was used as an experimental tool for suppressing the expression of these genes. Because neuroblastomas are notoriously difficult to transfect, we introduced an siRNA model with accurate detection of silencing efficiency and the resulting effects. In particular, for study of *MYCN* function, this model is believed to be more relevant, because traditional systems with forced overexpression of this gene in single-copy cells seem to lack the proper cellular context to mimic endogenous amplification and hyperactivity. Our final goal is to disentangle *MYCN*'s transcriptional web, in order to interfere with its oncogenic signaling pathways, while leaving the beneficial pathways unaltered.

From Experimental Design to Analysis of an rt-qPCR Assay

Purity and integrity of the template are critical factors to the success of an rt-qPCR assay. Several commercial kits are available for producing clean RNA samples. Contaminants should be avoided or removed, as they can greatly influence the reverse-transcription step or the actual PCR. The presence of PCR inhibitors can be determined by a variety of methods, including the simple and fast PCR-based SPUD assay (Nolan et al. 2006). An oligonucleotide target sequence with no homology to human DNA is spiked into human RNA samples and a water control at a known concentration. rt-qPCR quantitation of the oligonucleotide template in both the RNA samples and the (negative) water control is indicative of possible enzymatic inhibitors present in the RNA extract. For assessment of RNA

The Many Faces of Disease

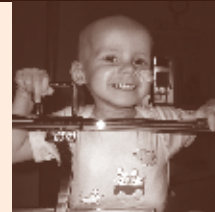
With the focus of his research in neuroblastoma, a very deadly form of childhood cancer, Professor Jo Vandesompele often gets asked if he meets the children behind the research. His answer: "No, we see a tube." In fact, most researchers spend countless hours with analytical tools, but little time, if any, interacting with people affected by disease. That's why the scientists that comprise Vandesompele's lab at Ghent University in Belgium are introducing a pilot program, where parents of children who have died or are suffering from neuroblastoma will be invited to speak to researchers about their experiences.

"Most of us don't maintain a sense of what we're doing research for," says Vandesompele. "A sample is brought from a hospital lab. We begin extracting molecules and conducting procedures that have nothing to do with the child the sample came from, a child who might be dying. There's a disconnect there that shouldn't be." The parent program is meant to bridge this disconnect.

The idea sprang from travels to international conferences, where parents involved in disease-related support groups occasionally give talks. Vandesompele's colleagues realized that in terms of motivating progress toward a cure, even the world's best scientists can't match the words of a parent whose child has died. And it's not just that parents have heartbreaking tales to tell. It's also that they have a passion for raising money to support research, and that they're truly interested in what's happening in the field.

"Yes, we're doing science," says Vandesompele, "but being connected to the human aspects of research can motivate scientists to be much more precise, closer to perfect in what we are doing. Passion brings us to a level unattainable based on intellectual skills alone."

Soon, at least in Belgium, researchers will begin to be able to match a name and a face to a test tube.



integrity, electrophoresis and PCR-based methods are available (Fleige and Pfaffl 2006, Nolan et al. 2006). Figure 1 shows an electropherogram of high-quality RNA assessed using the Experion™ automated electrophoresis system. Sharp peaks at 18S and 28S and no nonspecific peaks are desired results when determining whether or not RNA samples are intact.

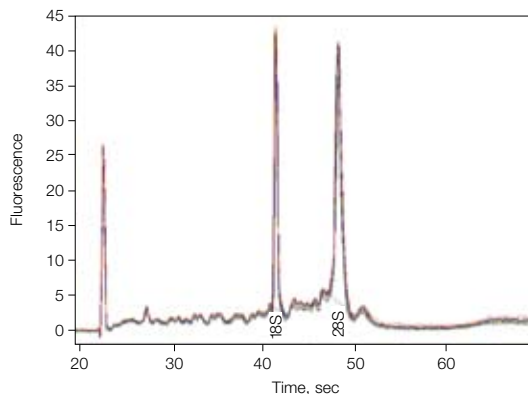


Fig. 1. Electropherogram generated using the Experion system depicting high-quality RNA sample.

To control for inevitable experimental variation due to factors such as the amount and quality of starting material, enzymatic efficiencies, and overall cellular transcriptional activity, use of a reliable normalization strategy in which these factors are taken into account is necessary. In principle, internal reference genes offer the best way to deal with the multiple sources of variables that might exist between different samples. A truly accurate normalization can only be achieved when multiple reference genes are utilized, as use of a single reference gene results in relatively large errors in a considerable proportion of the sample set (Vandesompele et al. 2002). Care should be taken when selecting the genes to be used for normalizing the expression levels since no universal set of always-applicable reference genes exists. Different sample origins and experimental manipulations might require another set of genes to be used as reference genes. The selection and validation of reference genes can be done using the geNorm algorithm (see sidebar), which determines the most stable genes from a set of tested candidate reference genes in a given sample panel and calculates a normalization factor (Vandesompele et al. 2002).

Bioinformatics-based quality assessment of newly designed rt-qPCR primers can considerably improve the likelihood of obtaining specific and efficient primers. A number of quality control parameters have been integrated in Ghent University's RTPPrimerDB in silico assay evaluation pipeline (Pattyn et al. 2006). This pipeline allows a streamlined evaluation of candidate primer pairs, with automated BLAST specificity search, prediction of putative secondary structures of the amplicon, indication of which transcript variants of the gene of interest will be amplified, and search for known SNPs in the primer annealing regions. This in silico evaluation, however, does not preclude the need for experimental validation after synthesis of the primers. Ideally, experimental evaluation addresses specificity, efficiency, and dynamic range of the assay using a broad dilution series of template (Figure 2).

Processing and analysis of the raw rt-qPCR data represent a multistep computational process of averaging replicate C_T values, normalization, and proper error propagation along the entire calculation track. This process might prove cumbersome and deserves equal attention as the previous steps in order to get accurate and reliable results. This final procedure has been automated and streamlined in Biogazelle's qBasePlus software (www.biogazelle.com, see sidebar), a dedicated program for the management and analysis of rt-qPCR data (Hellemans et al. 2007).

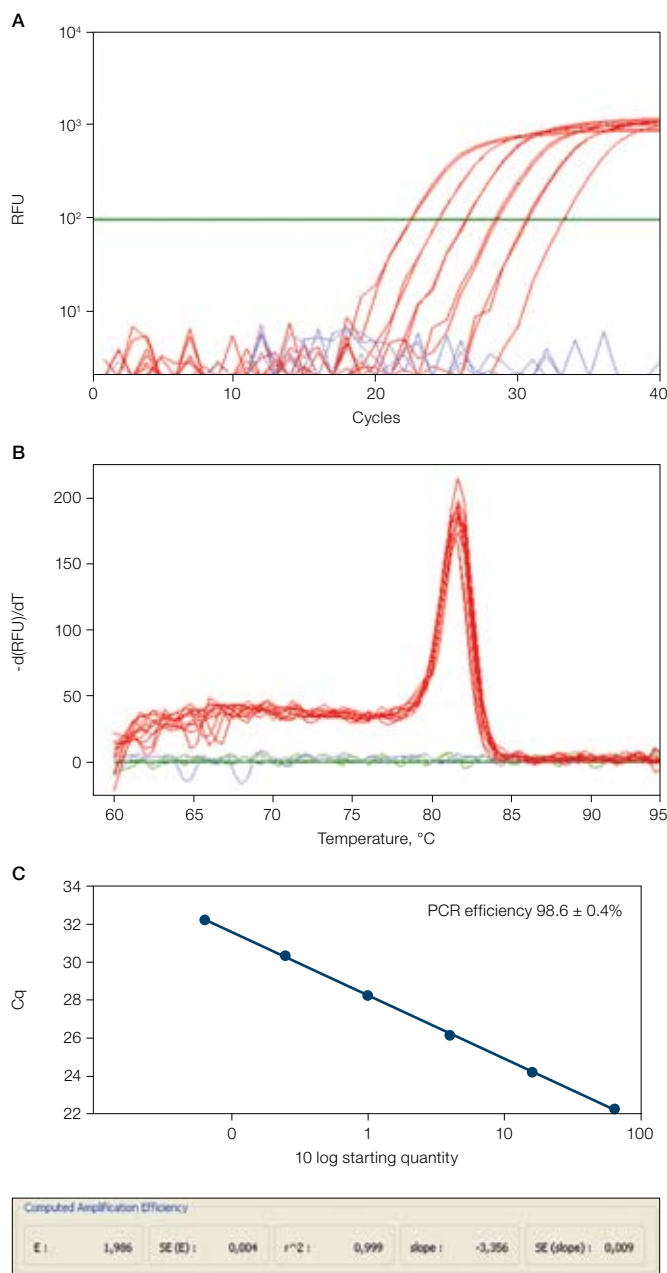


Fig. 2. Experimental validation of newly designed rt-qPCR primers. **A**, PCR efficiency and dynamic range of the rt-qPCR assay was tested using a 4-fold serial dilution of six points of reverse transcribed human qPCR reference total RNA (64 ng down to 62.5 pg) and TP53_P2 primers; **B**, specificity of the TP53_P2 primers was assessed by generating a melting curve of the PCR product; **C**, standard curve and PCR efficiency estimation (including the error) according to the qBasePlus software. Cq, quantitation cycle value generated in RDML software (see sidebar).

Programming Progress



The year 2000 is a milestone that symbolizes movement toward the height of progress, particularly in science and technology. But back in 2000, Professor Jo Vandesompele (then a doctorate student beginning what would become a career devoted to the study of the genetics of neuroblastoma at Ghent University in Belgium) was attempting to conduct sophisticated analysis of genetic research results with rudimentary tools.

"In 2000," says Vandesompele, "evaluating candidate reference genes with respect to their expression stability was impossible." Moreover, the concept of accurate normalization using multiple reference genes did not exist. "The problem of housekeeping gene variability was significantly underestimated at that time," he explains. In addition, he remembers calculating qPCR analyses by hand with Excel software, "a slow and error-prone process that required insight into mathematics and various quantitation models."

With no other solutions available, Vandesompele and colleagues set out to develop the first of many software, web, and database tools that continue to help drive progress in genetic research — not just in their lab, but in labs across the world. Launched in 2002, geNorm software is a tool used for the identification of stably expressed reference genes (<http://medgen.ugent.be/genorm/>). This launch was quickly followed by development of RTPrimerDB in 2003, a real-time PCR primer and probe database containing published and validated assays, as well as an integrated *in silico* PCR assay evaluation pipeline (<http://medgen.ugent.be/rtpriermdb/>).

In 2004, Jan Hellemans, a PhD student in the University's Center for Medical Genetics laboratory, began automating the arduous mathematical computations associated with qPCR analysis by programming a few simple macros in Excel. These initial macros evolved into the qBase 1.0 qPCR data analysis software package (<http://medgen.ugent.be/qbase/>). Since then, several thousand copies have been downloaded and used worldwide. In 2007, the Excel version began being phased out by qBasePlus, a professional Java-based application that runs 20 times faster and is more intuitive than the original platform. All current versions of these programs are available at no charge, and even this latest tool developed by Biogazelle, a Ghent University spin-off company, will offer both free and reasonably priced licensing packages.

That these programs have revolutionized the synthesis of real-time PCR data is unquestionable. What is surprising, at least to Vandesompele, is that "what were once just tools to measure gene expression levels in scarce tumor biopsies from children with neuroblastoma in our laboratory, have now grown in scope to form an independent research line."

And while researchers in this lab continue to try to find new ways to combat neuroblastoma, so will they continue to discover tools to aid achievement of reliable and meaningful results through bioinformatics. Future plans include establishment of an international consortium to finalize a standard exchange format for real-time PCR data (coined RDML, previews of this effort can be seen at www.rdml.org). In addition, they are developing a web-based primer design portal that will enable researchers to design high-quality assays in a high-throughput environment.

rt-qPCR for Assessment of siRNA Silencing Efficiency: Anti-MYCN siLentMer™ siRNA Duplexes

Human IMR-32 neuroblastoma cells were transfected with different anti-MYCN siLentMer siRNA duplexes or a nonspecific control siRNA, and the MYCN transcript level was determined 48 hours posttransfection by rt-qPCR. Our results indicate the importance of primer location for evaluation of siRNA silencing efficiency, in agreement with a previous independent report (Shepard et al. 2005). The target mRNA sequence is cleaved by the RNA-induced silencing complex (RISC) near the center of the region complementary to the guiding siRNA (Elbashir et al. 2001). Complete nucleolytic degradation of the resulting fragments is not always guaranteed, which might result in underestimation of siRNA silencing efficiency if primers are used that do not span the siRNA target sequence, as observed for this gene (Figure 3).

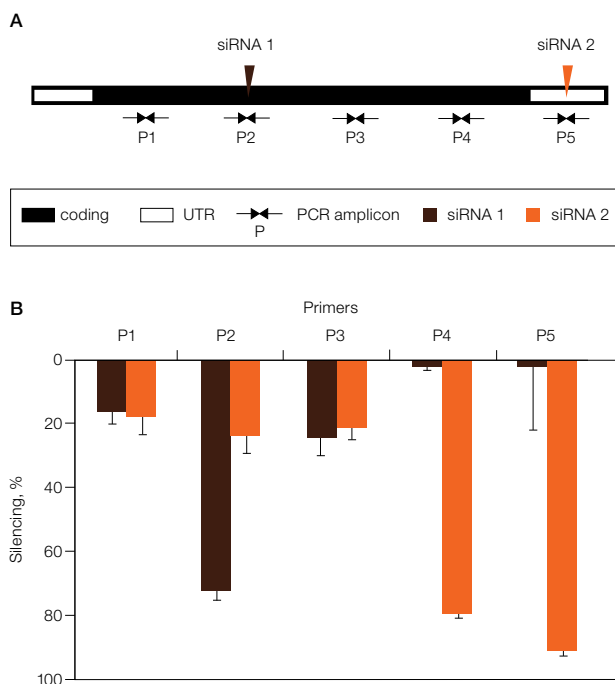


Fig. 3. Importance of primer location for rt-qPCR assessment of siRNA silencing efficiency. **A**, schematic representation of the MYCN mRNA structure with location of siRNA targeted sequences and primer binding sites; **B**, percentage silencing of MYCN gene expression measured by five different primer pairs (P1–P5) in IMR-32 cells 48 hr posttransfection with anti-MYCN siLentMer siRNA duplexes (siRNA 1 or siRNA 2), compared to cells transfected with a nonspecific control siRNA.

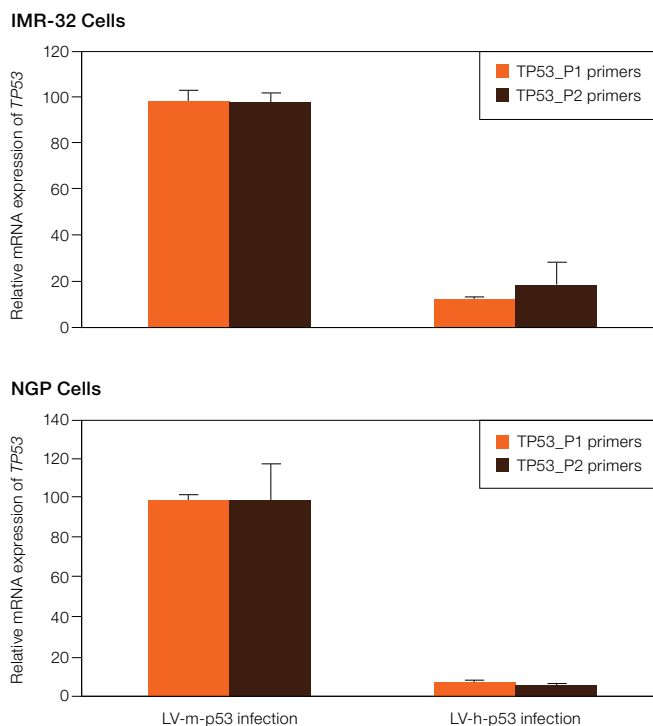


Fig. 4. Assessment of shRNA-mediated *TP53* knockdown efficiency by rt-qPCR. IMR-32 and NGP cells were infected with a lentivirus carrying an shRNA construct specific for either the human *TP53* gene (LV-h-p53) or the murine *Trp53* gene (LV-m-p53) as a control. Efficiency of *TP53* gene silencing was evaluated by rt-qPCR using two different primer pairs (TP53_P1 and TP53_P2). Bars indicate mRNA expression levels of *TP53* relative to the respective LV-m-p53 cells; error bars depict standard error of the mean (duplicated PCR reactions for *TP53* and three reference genes).

rt-qPCR for Monitoring of shRNA Silencing Efficiency and Functional Effects: Lentiviral-Mediated shRNA Knockdown of *TP53*

For generation of stable *TP53* knockdown variants of neuroblastoma cell lines with wild-type p53, we infected IMR-32 and NGP cells with a lentiviral vector encoding an shRNA directed specifically against human *TP53* (LV-h-p53) or against the murine *Trp53* gene (LV-m-p53, negative control). rt-qPCR analysis with two different primer pairs demonstrated that expression of *TP53* was reduced by 81–87% in IMR-32-LV-h-p53 cells and by 92–94% in NGP-LV-h-p53 cells compared to the respective LV-m-p53 controls (Figure 4). Functionality of the *TP53* knockdown variants was validated by rt-qPCR and cell viability analysis after treatment of the cells with nutlin-3, a small-molecule compound that selectively disrupts the interaction between p53 and its negative regulator *MDM2*, resulting in stabilization and accumulation of the p53 protein and activation of the p53 pathway (Vassilev et al. 2004). Transactivation of p53 target genes such as *BBC3* (*PUMA*) and *MDM2* by nutlin-3 and nutlin-3 induced downregulation

Designed for the Way You Want to Work

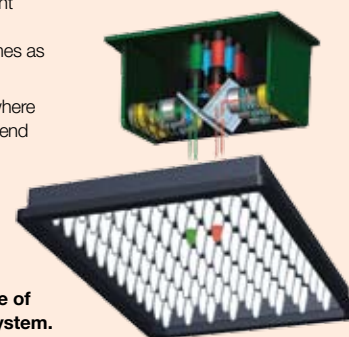
Use of a high-performance real-time qPCR system is important to accurately measure the effectivity of your siRNA knockdown. The CFX96™ real-time PCR detection system (used in the experiments discussed in this article) builds on the power and flexibility of the C1000™ thermal cycler, adding an easy-to-install interchangeable reaction module to create an exceptional real-time PCR system.

The system's thermal performance combined with an innovative optical design ensure accurate, reliable data. The powerful yet intuitive software accelerates every step of your real-time PCR research, shortening the time between getting started and getting great results.

The CFX96 system's solid-state optical technology (six filtered LEDs and six filtered photodiodes) provides sensitive detection for precise quantitation and target discrimination. Scanning just above the sample plate, the optics shuttle individually illuminates and reads fluorescence from each well with high sensitivity and no crosstalk. The optical system always collects data from all wells during data acquisition, so you can enter or edit well information on your own schedule.

With the CFX96 system, you can:

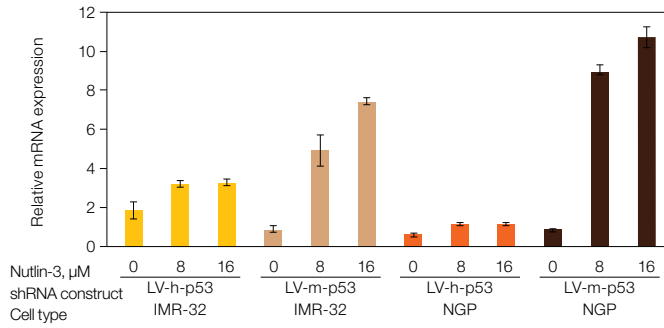
- Be up and running fast — quick installation and factory-calibrated optics let you set up the system in seconds
- Perform more experiments — fast thermal cycling produces results in <30 minutes
- Save research time — thermal gradient feature lets you optimize reactions in a single experiment
- Minimize sample and reagent usage — reliable results are obtained with sample volumes as low as 10 μ l
- Analyze results when and where you want — software can send e-mail notification with attached data file when the run is finished
- Trust your results — Security Edition software integrates the CFX96 system with good laboratory practice (GLP) standards for data collection and analysis
- Expand your throughput when you need to — up to 4 instruments can be controlled by a single computer



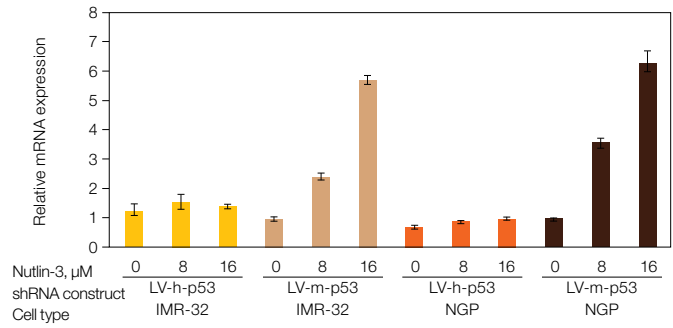
Six-channel optics shuttle of the CFX96 system.

of *TP53* mRNA level, a consequence of the ability of the p53 protein to negatively regulate its own transcriptional expression after accumulation (Hudson et al. 1995), was largely prevented by lentiviral-mediated expression of shRNA against human *TP53* (Figure 5). At the cellular level, silencing of human *TP53* severely attenuated the nutlin-3 induced reduction in cell viability observed in nontransduced parental cells, in contrast to control infection with LV-m-p53 (Figure 6). These results firmly demonstrate potent and selective impairment of p53 function in IMR-32-LV-h-p53 and NGP-LV-h-p53 cells.

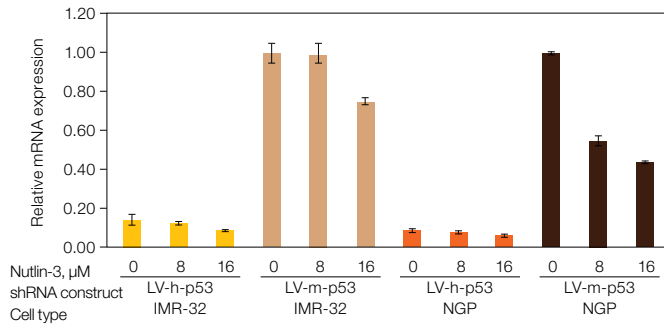
A. *BBC3 (PUMA)* expression



B. *MDM2* expression



C. *TP53* expression (using TP53_P1 primers)



D. *TP53* expression (using TP53_P2 primers)

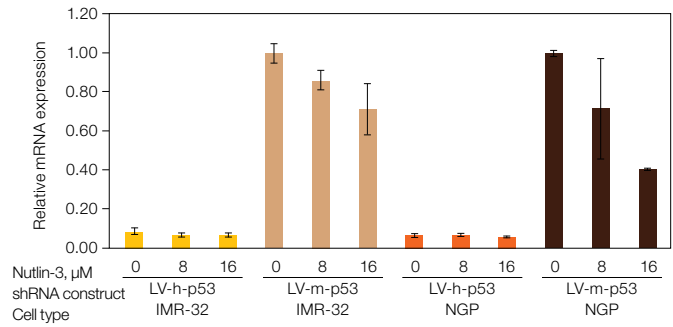
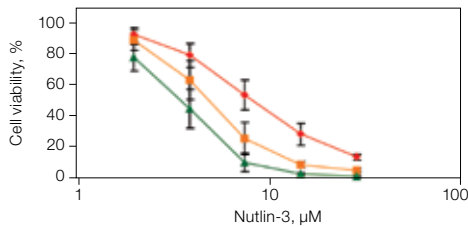
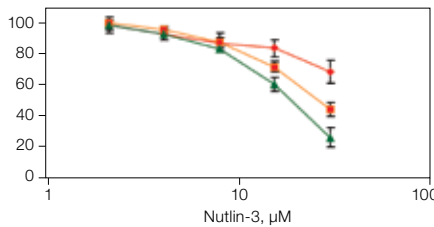


Fig. 5. Functional validation of shRNA-mediated *TP53* knockdown through rt-qPCR analysis of transcript levels of p53-regulated genes after nutlin-3 treatment. IMR-32 and NGP cells were infected with a lentiviral vector encoding an shRNA directed specifically against either the human *TP53* gene (LV-h-p53) or the murine *Trp53* gene (LV-m-p53). Cells were treated with 0, 8, or 16 μM nutlin-3 for 24 hr, and expression of *BBC3 (PUMA)* (A), and *MDM2* (B), p53-regulated genes, and *TP53* was determined by rt-qPCR. Two different primer pairs (TP53_P1 and TP53_P2) were used for quantitation of *TP53* transcript levels (C, D). Bars indicate mRNA expression levels relative to the respective vehicle-treated (0 μM nutlin-3) LV-m-p53 infected cells, mean of two different rt-qPCR measurements; error bars show standard error of the mean.

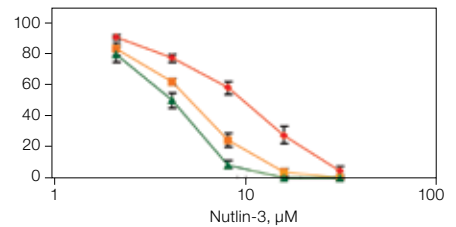
A. Uninfected IMR-32 cells



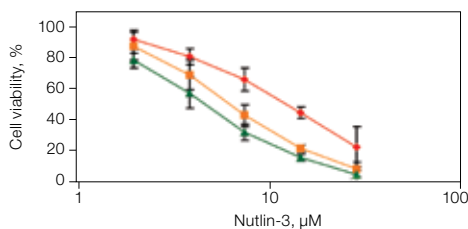
B. IMR-32 cells infected with LV-h-p53



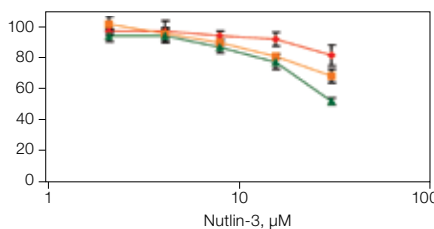
C. IMR-32 cells infected with LV-m-p53



D. Uninfected NGP cells



E. NGP cells infected with LV-h-p53



F. NGP cells infected with LV-m-p53

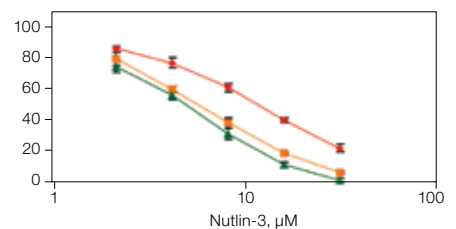


Fig. 6. Functional validation of shRNA-mediated *TP53* knockdown through cell viability analysis after treatment of IMR-32 and NGP cells with nutlin-3. Effect of nutlin-3 on viability of uninfected cells (A, D), LV-h-p53 infected cells (B, E) and LV-m-p53 infected cells (C, F). Exponentially growing cells were exposed to 0–32 μM of nutlin-3 for 24 (→), 48 (→), and 72 (→) hr, and the percentage cell viability with respect to vehicle-treated cells was determined. Error bars indicate standard deviation of mean cell viability values of three independent experiments.

Conclusions

rt-qPCR analysis provides a convenient and reliable method for evaluation of knockdown efficiency and functional consequences of RNAi-mediated gene silencing. Successful application of this monitoring tool requires careful attention to be given to all different steps in the rt-qPCR workflow, including primer design and evaluation, template preparation, normalization strategy, and data analysis, as discussed in this article.

Similar studies will be conducted in the future to evaluate results achieved using additional cell lines and varying combinations of multiple siLentMer duplexes, durations of effect, and concentrations of active siLentMer duplexes.

References

- Elbashir SM et al., RNA interference is mediated by 21- and 22-nucleotide RNAs, *Genes Dev* 15, 188–200 (2001)
- Fleige S and Pfaffl MW, RNA integrity and the effect on the real-time qRT-PCR performance, *Mol Aspects Med* 27, 126–139 (2006)
- Hellemans J et al., qBase relative quantification framework and software for management and automated analysis of real-time quantitative PCR data, *Genome Biol* 8, R19 (2007)
- Hudson JM et al., Wild-type p53 regulates its own transcription in a cell-type specific manner, *DNA Cell Biol* 14, 759–766 (1995)
- Nolan T et al., Quantification of mRNA using real-time RT-PCR, *Nat Protoc* 1, 1559–1582 (2006)
- Nolan T et al., SPUD: a quantitative PCR assay for the detection of inhibitors in nucleic acid preparations, *Anal Biochem* 351, 308–310 (2006)
- Pattyn F et al., RTPrimerDB: the real-time PCR primer and probe database, major update 2006, *Nucleic Acids Res* 34, D684–D688 (2006)
- Shepard AR et al., Importance of quantitative PCR primer location for short interfering RNA efficacy determination, *Anal Biochem* 344, 287–288 (2005)
- Tweddle DA et al., The p53 pathway and its inactivation in neuroblastoma, *Cancer Lett* 197, 93–98 (2003)
- Vandesompele J et al., Unequivocal delineation of clinicogenetic subgroups and development of a new model for improved outcome prediction in neuroblastoma, *J Clin Oncol* 23, 2280–2299 (2005)
- Vandesompele J et al., Accurate normalization of real-time quantitative RT-PCR data by geometric averaging of multiple internal control genes, *Genome Biol* 3, RESEARCH0034 (2002)
- Van Maerken T et al., Small-molecule MDM2 antagonists as a new therapy concept for neuroblastoma, *Cancer Res* 66, 9646–9655 (2006)
- Vassilev LT et al., In vivo activation of the p53 pathway by small-molecule antagonists of MDM2, *Science* 303, 844–848 (2004)
- Grant support: the Fund for Scientific Research – Flanders (FWO) grants 011F4004 (T. Van Maerken, Research Assistant), G.1.5.243.05 (J. Vandesompele) and G.0185.04; GOA grant 12051203; grants from the Belgian Foundation against Cancer (S. De Clercq), the Ghent Childhood Cancer Fund, and the Ghent University Research grant (B.O.F.) 01D31406 (P. Mestdagh).
- Contact information: Jo Vandesompele, Center for Medical Genetics, Ghent University Hospital, MRB, De Pintelaan 185, 9000 Ghent, Belgium. Phone: 32-9-332-5187; Fax: 32-9-332-6549; E-mail: Joke.Vandesompele@UGent.be

BLAST is a trademark of the National Library of Medicine. Excel is a trademark of Microsoft Corporation.



LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only.

The dye(s) used in Expiration kits are manufactured by Molecular Probes, Inc. and are licensed for research use only. The siLentMer products are manufactured by Integrated DNA Technologies, Inc. (IDT) and are for research use only. For custom siRNA synthesis, contact IDT. Notice regarding Bio-Rad thermal cyclers and real-time systems — Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents Nos. 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160–163 and 167 only), and 6,703,236 (claims 1–7 only), or corresponding claims in their non-U.S. counterparts, owned by Applera Corporation. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA. Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics.



**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000