

# Analysis of IL-4 Dependent Gene Expression in Namalwa Cells by siRNA Transfection: An Example of Pathway Analysis Using the Gene Pulser MXcell™ Electroporation System

Claudia M. Litterst,<sup>1</sup> Tobias Münz,<sup>2</sup> Edith Pfützner,<sup>2</sup> and Luis Ugozzoli<sup>1</sup>  
<sup>1</sup>Bio-Rad Laboratories, Inc., Hercules, CA 94547 USA; <sup>2</sup>Department of Biochemistry and Biophysics, Friedrich-Schiller-Universität Jena, Center for Molecular Biomedicine, Hans-Knöll-Straße 2, D-07745 Jena, Germany

## Introduction

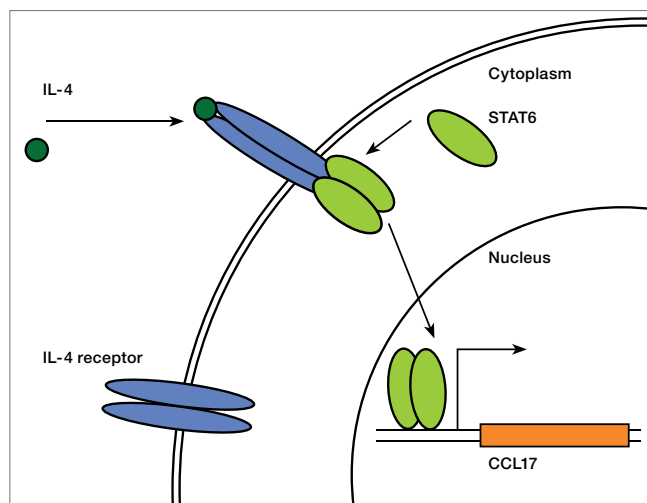
siRNA-mediated knockdown of gene expression is a powerful functional genomics tool to interrogate signaling pathways in basic research and drug discovery. However, such studies can be technically challenging since they require efficient siRNA delivery to achieve significant reduction of protein levels. In addition, the siRNA transfection procedure must not perturb the cell physiology and alter the signaling pathway under investigation.

Figure 1 shows a simplified overview of the IL-4 signaling pathway. Binding of IL-4 to the IL-4 receptor leads to dimerization of the signal transducer and activator of transcription 6 (STAT6). The activated STAT6 dimer translocates to the nucleus and drives the expression of its target genes, for example, thymus and activation-regulated chemokine (TARC/CCL17, hereafter referred to as CCL17) (Wirnsberger et al. 2006). CCL17 is a high-affinity ligand for the CC chemokine receptor 4 (CCR4 receptor) and acts as a selective chemoattractant for T cells expressing this receptor (Imai et al. 1997, Imai et al. 1999). It has been suggested that CCL17 produced by B cells plays an important role in the pathogenesis of allergic disease (Lin et al. 2003).

The difficult-to-transfect Burkitt lymphoma cell line Namalwa, is a recognized model system to study the IL-4 signaling pathway (Rousset et al. 1988). In this article we describe the development of an optimized electroporation protocol for siRNA transfection of Namalwa cells using the Gene Pulser MXcell electroporation system along with Gene Pulser® electroporation buffer for siRNA transfection, and the application of siRNA-mediated gene silencing in a functional study on IL-4 dependent gene expression.

## Methods

Namalwa cells (American Type Culture Collection, ATCC #CRL-1432) were grown in GIBCO RPMI Media 1640 (Invitrogen Corporation) with 7.5% FBS and 1% pyruvate.

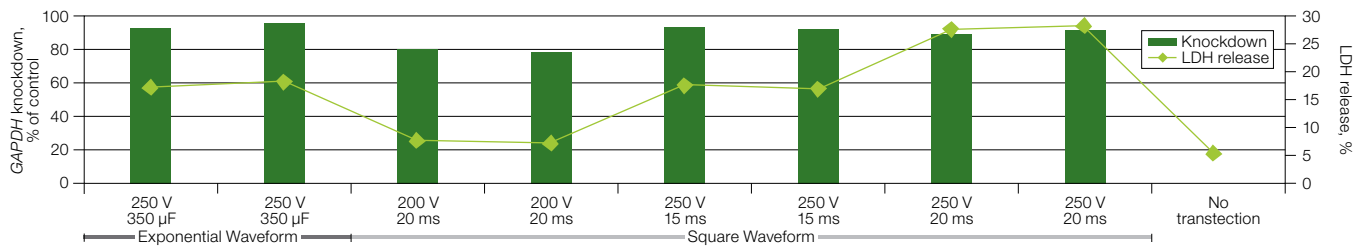


**Fig. 1. Overview of the IL-4 signaling pathway.** Binding of IL-4 to its receptor leads to dimerization and activation of STAT6. Activated STAT6 translocates to the nucleus where it binds to the promoter region of its target genes, for example, CCL17, and activates transcription.

Recombinant human IL-4 was purchased from R&D Systems, Inc. and was used at 10 ng/ml final concentration. STAT6 siRNAs were purchased from Integrated DNA Technologies, Inc. (IDT); all other siRNAs were from Bio-Rad.

For electroporation, cells were washed in PBS and resuspended in Gene Pulser electroporation buffer (Bio-Rad Laboratories, Inc.) at a density of  $5 \times 10^6$  cells/ml. siRNAs were added at 100 nM final concentration. The cell suspension was mixed and 150  $\mu$ l suspension/well was transferred to the appropriate wells of a 96-well electroporation plate (Bio-Rad). Electroporations were carried out on the Gene Pulser MXcell electroporation system (Bio-Rad). After electroporation, cell suspensions were immediately transferred to growth media ( $7.5 \times 10^5$  cells/ml final concentration) and incubated for 24 hr.

LDH levels in cell culture supernatants were determined using the CytoScan LDH Cytotoxicity Assay (G Biosciences). Relative LDH levels were calculated in reference to untreated lysed cells according to the manufacturer's recommendations. RNA was



**Fig. 2. Optimization of electroporation conditions.** Namalwa cells were electroporated with nonsilencing control siRNA and *GAPDH* siRNA using the Gene Pulser MXcell system with the settings at 1000  $\Omega$ . Capacitance was set to 1000  $\mu$ F for square-wave conditions. Twenty-four hours after transfection, cell supernatants were collected and subjected to LDH release assay. Total RNA was prepared from cells and relative levels of *GAPDH* mRNA were determined by RT-qPCR using the CFX384 real-time PCR detection system. Relative *GAPDH* mRNA levels were determined with respect to the levels measured in cells transfected with the nonsilencing control siRNA for each condition.

purified using the Aurum™ total RNA 96 kit, then cDNA was prepared with the iScript™ cDNA synthesis kit (both kits from Bio-Rad). Real-time PCR reactions were performed using iQ™ supermix on the CFX384™ real-time PCR detection system, and data was analyzed using CFX Manager™ software (all from Bio-Rad). Primers for amplification of the actin gene, *GAPDH*, and the *CCL17* gene were purchased from IDT. Primer sequences for *CCL17* were described previously (Radstake et al. 2005). Primer sequences for *GAPDH* and the actin gene can be obtained upon request. Actin was used as a reference gene for the analysis shown in Figure 2, and *GAPDH* was the reference gene for the experiment shown in Figure 3A.

For western blotting, cells were lysed in lysis buffer (1% Triton, 20 mM Tris-HCl, pH 8, 0.1% SDS, 150 mM NaCl) containing Protease Inhibitor Cocktail Set III (Calbiochem). Protein concentrations were determined using the DC protein assay (Bio-Rad). Total protein (50  $\mu$ g/sample) was separated on 10% Ready Gel® Tris-HCl gels (Bio-Rad) and transferred to PVDF membranes using the Mini-PROTEAN® Tetra system (Bio-Rad). Membranes were probed with BD Transduction Laboratories STAT6 (BD Biosciences) and alpha-tubulin (Sigma-Aldrich, Inc.) antibodies. Western blots were developed using HRP-conjugated secondary antibodies according to standard procedures.

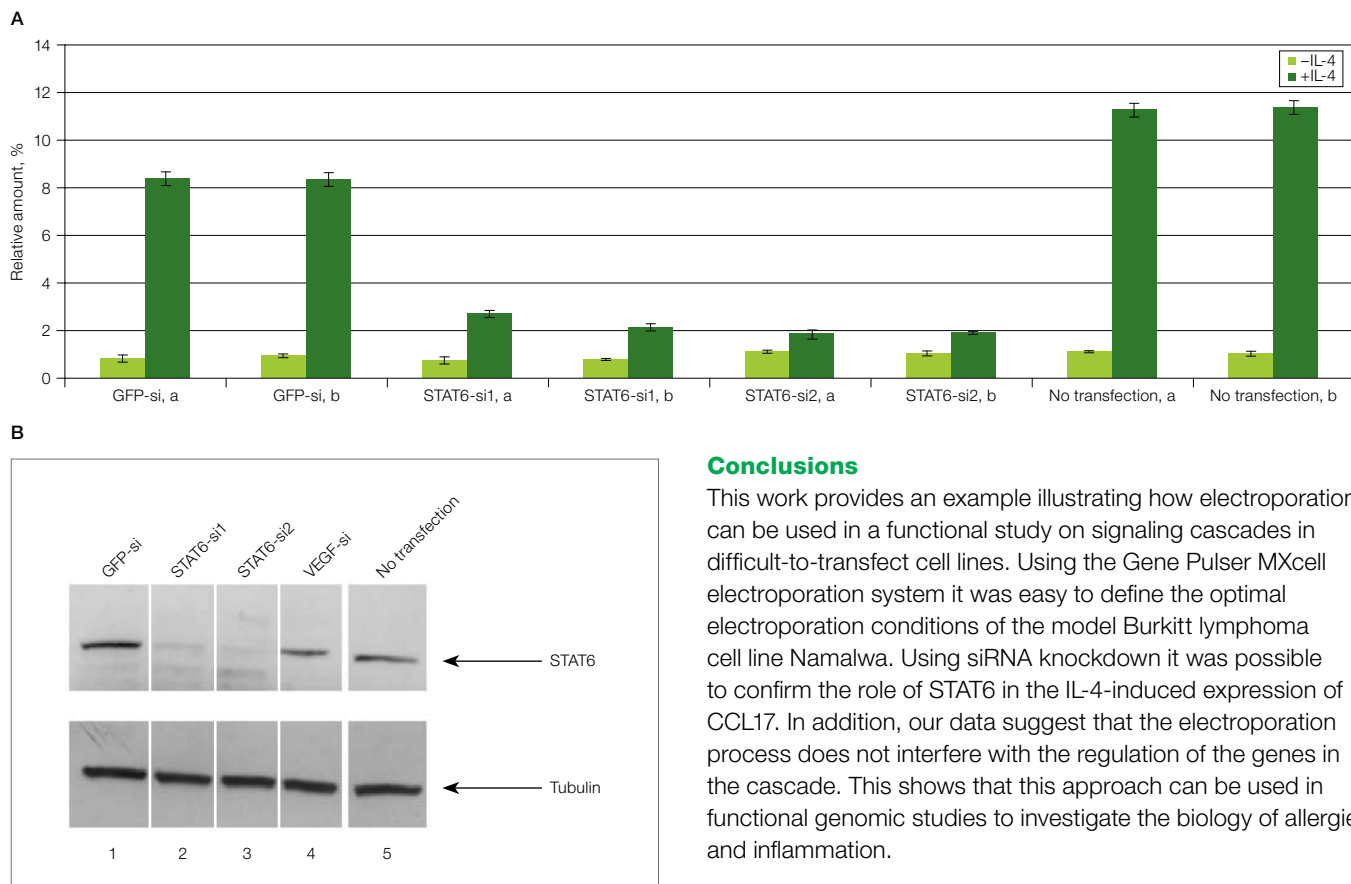
## Results and Discussion

Preliminary experiments were first performed to identify optimal electroporation conditions for siRNA delivery into Namalwa cells. A range of settings (shown in Figure 2), that had been previously optimized for electroporation of plasmid DNA in Namalwa cells (Litterst and Ugozzoli 2009), was screened. Namalwa cells were electroporated with *GAPDH* and a nonsilencing control siRNA for each condition. Twenty-four hours after transfection, cytotoxicity and knockdown of *GAPDH* expression were analyzed.

As shown in Figure 2, strong *GAPDH* knockdown (>90%) was obtained with both exponential-decay and square-wave conditions (250 V/350  $\mu$ F and 250 V/15 ms, respectively). LDH levels (a measure of cell lysis) were comparable for both

of these conditions (approximately 17% relative LDH release), indicating that both settings result in efficient siRNA transfection of Namalwa cells with minimal effect on cell viability. Other square-wave conditions were either less efficient with respect to knockdown (200 V/20 ms) or resulted in higher cytotoxicity (250 V/20 ms) and are therefore not recommended for siRNA transfection of Namalwa cells. The exponential-decay condition (250 V/350  $\mu$ F) was also the optimal setting for DNA transfection (Litterst and Ugozzoli 2009), indicating that once conditions are established for DNA delivery, they can be used for efficient siRNA transfection as well. This condition was subsequently used for our IL-4 dependent gene expression analysis.

Namalwa cells were transfected with GFP siRNA and two STAT6 siRNAs as indicated in Figure 3A. Twenty-four hours after transfection, one set of replicates was treated with recombinant IL-4 and levels of *CCL17* mRNA were determined after 24 hr by reverse transcription quantitative PCR (RT-qPCR). As shown in Figure 3A, control samples transfected with GFP siRNA showed similar induction of *CCL17* expression (8-fold induction) upon IL-4 treatment as nontransfected cells (11-fold induction), indicating that the electroporation procedure does not significantly interfere with the IL-4 pathway and the cells retain their ability to upregulate *CCL17* after siRNA transfection. IL-4 treatment of samples transfected with STAT6 siRNAs induced expression of *CCL17* mRNA only 2-fold. To confirm that the reduced response of transfected cells to IL-4 was due to knockdown of STAT6, STAT6 protein levels were analyzed in transfected cells by western blotting. siRNA transfection of STAT6 resulted in almost complete depletion of the STAT6 protein (Figure 3B, lanes 2 and 3) compared to the cells transfected with control siRNA (Figure 3B, lanes 1 and 4) and nontransfected cells (Figure 3B, lane 5), while tubulin protein levels were similar in all samples. These results are in agreement with previous studies showing that IL-4 induced expression of *CCL17* is STAT6 dependent (Wirnsberger et al. 2006). Control samples transfected with *GAPDH* siRNA showed an up to 86% reduction of *GAPDH* mRNA after 48 hr (data not shown), indicating that the knockdown was still effective at the time expression of *CCL17* was measured.



**Fig. 3. STAT6 is required for IL-4-induced CCL17 expression.** **A**, cells were transfected with control siRNAs (GFP-si), STAT6 siRNA (STAT6-si1, STAT6-si2), or not transfected as indicated. Each transfection was performed in four replicates. Twenty-four hours after transfection, two samples per set of replicates were treated with IL-4 (■) and two samples were left untreated (□). Forty-eight hours after transfection, RNA was prepared from cells and CCL17 mRNA levels were determined by RT-qPCR. Relative levels of CCL17 expression were calculated with respect to the GFP siRNA-transfected sample without IL-4; **B**, cells transfected with STAT6 (STAT6-si1, STAT6-si2) and control siRNAs (GFP-si, VEGF-si) as indicated in 3A were lysed 48 hr after transfection. VEGF-siRNA was used as an additional unrelated control to monitor variations in protein levels. Total protein extract (50 µg) was resolved by SDS-PAGE. Protein levels of STAT6 and tubulin were analyzed by western blotting using specific antibodies.

## Conclusions

This work provides an example illustrating how electroporation can be used in a functional study on signaling cascades in difficult-to-transfect cell lines. Using the Gene Pulser MXcell electroporation system it was easy to define the optimal electroporation conditions of the model Burkitt lymphoma cell line Namalwa. Using siRNA knockdown it was possible to confirm the role of STAT6 in the IL-4-induced expression of CCL17. In addition, our data suggest that the electroporation process does not interfere with the regulation of the genes in the cascade. This shows that this approach can be used in functional genomic studies to investigate the biology of allergies and inflammation.

## References

- Imai T et al. (1997). The T cell-directed CC chemokine TARC is a highly specific biological ligand for CC chemokine receptor 4. *J Biol Chem* 272, 15036-15042.
- Imai T et al. (1999). Selective recruitment of CCR4-bearing Th2 cells toward antigen-presenting cells by the CC chemokines thymus and activation-regulated chemokine and macrophage-derived chemokine. *Int Immunol* 11, 81-88.
- Lin L et al. (2003). TARC and MDC are produced by CD40 activated human B cells and are elevated in the sera of infantile atopic dermatitis patients. *J Med Dent Sci* 50, 27-33.
- Litterst CM and Ugozzoli L (2009). Optimization of electroporation conditions for two different Burkitt lymphoma cell lines using the Gene Pulser MXcell™ system. *BioRadiations* 128, 22-23.
- Radstake TR et al. (2005). Increased expression of CCL18, CCL19, and CCL17 by dendritic cells from patients with rheumatoid arthritis, and regulation by Fc gamma receptors. *Ann Rheum Dis* 64, 359-367.
- Rousset F et al. (1988). Regulation of Fc receptor for IgE (CD23) and class II MHC antigen expression on Burkitt's lymphoma cell lines by human IL-4 and IFN-gamma. *J Immunol* 140, 2625-2632.
- Wirnsberger G et al. (2006). IL-4 induces expression of TARC/CCL17 via two STAT6 binding sites. *Eur J Immunol* 36, 1882-1891.

The ATCC trademark and ATCC catalog number CRL-1432 are trademarks of the American Type Culture Collection. BD Transduction Laboratories is a trademark of Becton, Dickinson and Company. CytoScan is a trademark of Geno Technology Inc. GIBCO is a trademark of Invitrogen Corporation.

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of iQ supermix includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from 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 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.

Information in this tech note was current as of the date of writing (2009) and not necessarily the date this version (rev A, 2009) was published.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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