

Reverse Transfection of Mammalian Cells for Functional Screening: Results from Testing the BioOdyssey™ Calligrapher™ MiniArrayer

Saman Honarnejad,^{1,3} Rainer Pepperkok,³ and Holger Erfle^{2,3}

¹ Department of Biotechnology, Mannheim University of Applied Sciences, Windeckstrasse 110, D-68163, Mannheim, Germany; ² MitoCheck Project Group, ³ Cell Biology/Biophysics, European Molecular Biology Laboratory, Meyerhofstrasse 1, D-69117, Heidelberg, Germany

Introduction

RNA interference (RNAi) is a recent advancement that can reduce the expression of specific target genes, with potential applications on a genome-wide scale. On the other hand, overexpression of organelle-specific markers (GFP-tagged cDNAs) using reverse transfection can be the method of choice for functional screening-based experiments. By using 8-well Lab-Tek Chamber Slides (Nunc), it is possible to have eight replicates of a subgrid, consisting of different spotted small interfering RNAs (siRNAs) or cDNAs that are ready to transfect and use for different applications. One such use in drug discovery is to increase the efficiency of transfection or treatments by influencing downregulation or overexpression by adding certain biological or chemical compounds to each well, and simultaneously monitoring the effects by using time-lapse high-throughput screening. However, to achieve this, robust methodologies that allow automated and efficient delivery of siRNAs or cDNAs into living cultured cells, and reliable quality control of siRNA or cDNA function, must be in place.

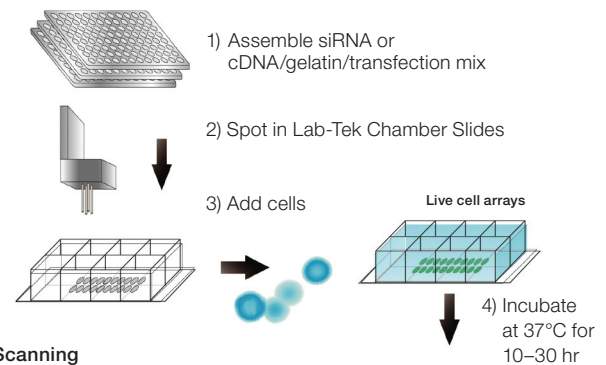
Here, we describe optimization of reverse transfection of cultured cells with siRNA or cDNA for subsequent high-content screening microscopy applications (Figure 1). All the necessary transfection components are mixed prior to robotic spotting on noncoated 8-chambered cover glass tissue culture dishes, which are ideally suited for time-lapse microscopy applications in living cells. The addition of fibronectin to the spotting solution improves cell adherence and transfection efficiency. After cell seeding, no further cell culture manipulation, such as medium changes, is needed. Calculation of the cell density improves autofocus performance for high-quality data acquisition and cell recognition.

In this report, we confirm known mitotic phenotypes after knockdown with specific siRNAs, such as INCENP (multinucleated cells, segregation problems), PLK1 (prometaphase arrest), and TPX2 (prometaphase arrest),

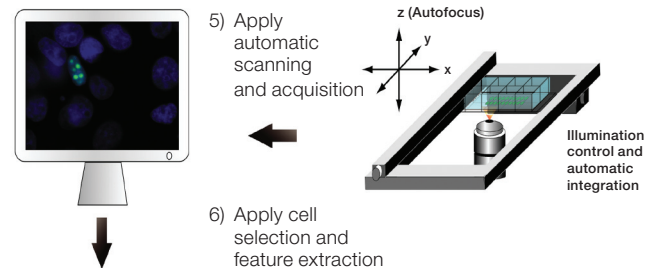
with a transfection efficiency of up to 80%. We also studied localization of different GFP-tagged cDNAs in different organelles and constructs in mammalian cells, with a transfection efficiency of up to 43% (data not shown).

Methods

Arraying



Scanning



Classification



Fig. 1. Reverse transfection and automatic image acquisition workflow. 1, different solutions of transfection mix, cDNA or siRNA, and gelatin are prepared in 384-well plates; 2, solutions are spotted in arrays into chambered cover glass tissue culture dishes (Lab-Tek Chamber Slides, Nunc), using the BioOdyssey Calligrapher miniarrayer; 3, cultured cells are added to the Chamber Slides and transfected with cDNA or siRNA on the arrayed spots; 4, after 10–30 hr of incubation, the fluorescent signal of the expressed GFP-labeled protein can be visualized; 5, on the automatic scanning platform, a motorized stage and motorized z-stepper perform scanning of live cells; 6, known phenotypes are selected and further extracted with different software algorithms; 7, phenotypes are identified by trained classifier (Conrad et al. 2004).

Preparation of Transfection Mix

siRNA — In a 384-well plate, 3 μl of 0.4 M sucrose in Opti-MEM I with GlutaMAX (Invitrogen Corporation) was mixed with 3.5 μl of Lipofectamine 2000 (Invitrogen) and 5 μl of siRNA (390 ng/ μl). The mixture was incubated at room temperature for 20 min, so that the transfection complexes or siRNA (Integrated DNA Technologies, IDT) could form. Then 7.25 μl of gelatin/fibronectin mixture with a ratio of 100/1 (v/v), was pipetted into each well, bringing the total volume to 18.75 μl . The mixture was incubated again for 15 min followed by short centrifugation (Erfle et al. 2007).

cDNA — The same protocol that was used for preparing the siRNA mix was used to prepare the cDNA mix, with the exception that 2 μl of cDNA (500 ng/ μl) was used and the total volume of the mixture was 16.75 μl .

Spotting

The BioOdyssey Calligrapher miniarrayer was used to print eight replicate subgrids with 25 (5 \times 5) spots in each well of a Lab-Tek Chamber Slide, for a total of 200 spots printed on a single slide (Figure 2). One solid pin used for spotting was washed with distilled water. Humidity was set at 57% to avoid sample evaporation.

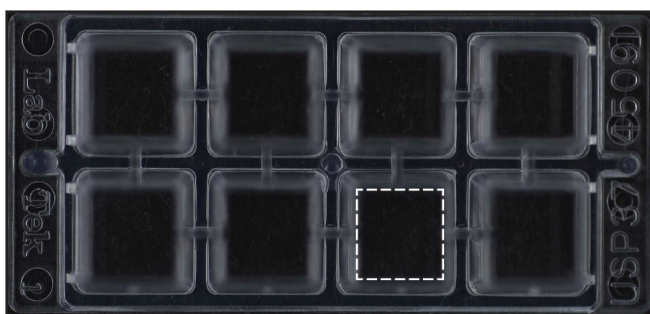


Fig. 2. 8-well Lab-Tek Chamber Slide. Each slide is 74 \times 22 mm with a thickness of 0.13–0.17 mm and culture area of 0.8 cm^2 . The dashed square presents a well of the Chamber Slide (see Figure 3 for a zoomed spotted version).

The middle spot/point of each well was marked with a pen before seeding of cells to provide the position coordinates for the microscope software (Figure 3). Each well contained several different siRNAs (Figure 4).

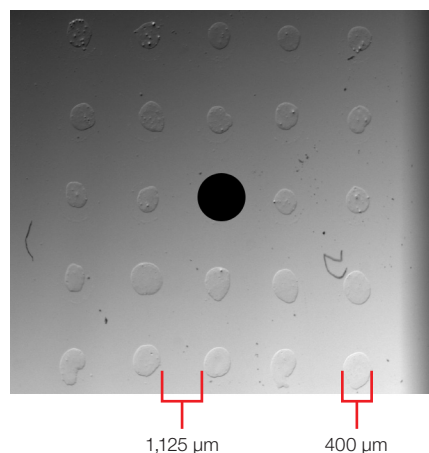


Fig. 3. Dried spots in one well of a Lab-Tek Chamber Slide. A total of 25 spots with a distance between each spot of $\sim 1,125 \mu\text{m}$ and diameters of $\sim 400 \mu\text{m}$. The black dot in the center is a mark made by a pen on the external surface of the slide to provide the position for the microscope software.

INCENP-A1	INCENP-A2	INCENP-A3	TPX2-A1	TPX2-A2
TPX2-A3	TPX2-B1	TPX2-B2	TPX2-B3	PLK1-A1
PLK1-A2	PLK1-A3	●	PLK1-B1	PLK1-B2
PLK1-B3	INCENP-B1	INCENP-B2	INCENP-B3	TPX2-C1
TPX2-C2	TPX2-C3	INCENP-C1	INCENP-C2	INCENP-C3

Fig. 4. Different siRNAs spotted in a single well of a Lab-Tek Chamber Slide.

Seeding of HeLa “Kyoto” (H2B-GFP) Cells

Cells were passaged the day before they were seeded. For each well of a Lab-Tek Chamber Slide, approximately 3 drops (100 μl) of a cell suspension was determined to be optimal for ease of pipetting and to prevent desiccation. HeLa “Kyoto” cells (Neumann et al. 2006) were seeded at a density of 0.4×10^{-5} cell/ μl using a 10 ml pipet to plate drops in the middle of the well while shaking slightly (in a figure-eight movement) to obtain a homogenous distribution of the cells on the surface of each well. Because the seeding procedure stresses the cells, they were immediately returned to the 5% CO_2 incubator and incubated for ~ 24 –30 hr at 37°C.

Acquisition of Images

Images were taken with a high-throughput automated microscope (scan[^]R screening station, Olympus Europa GmbH). By fixing the midpoint of each well, and calculating the distance between spots and number of subgrids, the microscope captures images from the position of interest using a laser autofocus system (Figure 5).

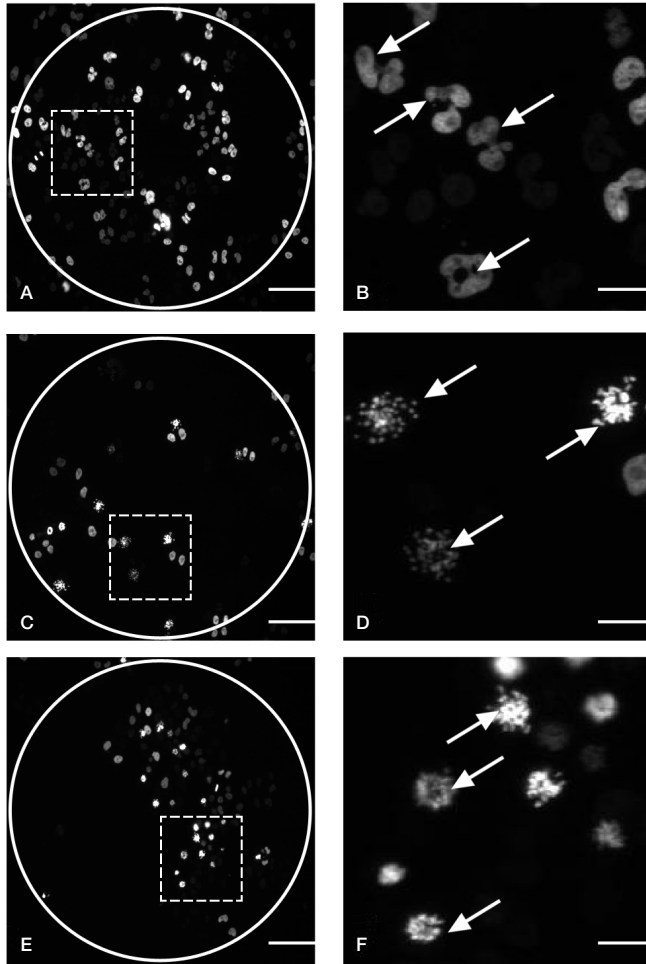


Fig. 5. Knockdown of three targets in HeLa "Kyoto" (H2B-GFP) cell line. Confirmation of known mitotic phenotypes, such as INCENP (multinucleated cells, segregation problems) (A and B); PLK1 (prometaphase arrest) (C and D), and TPX2 (prometaphase arrest) (E and F), with an efficiency of up to 80%. Panels B, D, and F are high-magnification versions of panels A, C, and E, respectively. Arrows point to phenotypes with mitotic defects. Bars, 20 μ m (A, C, and E) and \sim 5 μ m (B, D, and F).

Conclusions

We have demonstrated that effective gene silencing can be obtained using reverse transfection in plates with multiple wells. Each well of the Lab-Tek Chamber Slide possesses a number of different siRNAs, allowing us to assay multiple conditions simultaneously. In order to use these plates, an instrument that is capable of accurate spotting is required. The BioOdyssey Calligrapher miniarrayer fulfilled this requirement, allowing accurate placement of spots within the grids that were then imaged after cells were transfected. The BioOdyssey Calligrapher miniarrayer can be easily integrated into a reverse transfection protocol to successfully perform functional screening experiments.

Acknowledgements

We thank IDT for providing us with the siRNA oligonucleotides, and Bio-Rad Laboratories for the collaboration and giving us the chance to test the BioOdyssey Calligrapher miniarrayer. In addition, we want to thank the members of the MitoCheck Project Group and the Rainer Pepperkok team at the European Molecular Biology Laboratory (EMBL).

References

Conrad C et al., Automatic identification of subcellular phenotypes on human cell arrays, *Genome Res* 14, 1130–1136 (2004)

Erfle H et al., Reverse transfection on cell arrays for high content screening microscopy, *Nat Protoc* 2, 392–399 (2007)

Neumann B et al., High-throughput RNAi screening by time-lapse imaging of live human cells, *Nat Methods* 3, 385–390 (2006)

Bibliography

Erfle H and Pepperkok R, Arrays of transfected mammalian cells for high content screening microscopy, *Methods Enzymol* 404, 1–8 (2005)

Erfle H and Pepperkok R, Production of siRNA- and cDNA-transfected cell arrays on noncoated chambered coverglass for high-content screening microscopy in living cells, *Methods Mol Biol* 360, 155–161 (2007)

Erfle H et al., siRNA cell arrays for high-content screening microscopy, *Biotechniques* 37, 454–458, 460, 462 (2004)

Liebel U et al., A microscope-based screening platform for large-scale functional protein analysis in intact cells, *FEBS Lett* 554, 394–398 (2003)

Mishina YM et al., Multiplex GPCR assay in reverse transfection cell microarrays, *J Biomol Screen* 9, 196–207 (2004)

Pepperkok R and Ellenberg J, High-throughput fluorescence microscopy for systems biology, *Nat Rev Mol Cell Biol* 7, 690–696 (2006)

Simpson JC et al., Systematic subcellular localization of novel proteins identified by large-scale cDNA sequencing, *EMBO Rep* 1, 287–292 (2000)

Starkuviene V et al., High-content screening microscopy identifies novel proteins with a putative role in secretory membrane traffic, *Genome Res* 14, 1948–1956 (2004)

Wheeler DB et al., Cell microarrays and RNA interference chip away at gene function, *Nat Genet* 37, S25–30 (2005)

Wiemann S et al., The German cDNA network: cDNAs, functional genomics and proteomics, *J Struct Funct Genomics* 4, 87–96 (2003)

Ziauddin J and Sabatini DM, Microarrays of cells expressing defined cDNAs, *Nature* 411, 107–110 (2001)

Certain arrays and/or the method of preparation, analysis, or use may be covered by patents or other intellectual property rights owned by third parties.

Chamber Slides and Lab-Tek are trademarks of Thermo Fisher Scientific, Inc. GlutaMAX, Lipofectamine, and Opti-MEM are trademarks of Invitrogen Corporation. scan[^]R is a trademark of Olympus Europa GmbH.

Information in this tech note was current as of the date of writing (2007) and not necessarily the date this version (rev A, 2007) was published.



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