

Harvesting Organoids for Biochemical Analysis



Trevigen®

Better Products. Better Results.

Sol Degese¹, Gabe Benton¹.

¹Organoid Resource Lab (ORL), Trevigen, Inc., 8405 Helgerman Court, Gaithersburg, MD 20877

Introduction

3D Cell culture models that resemble the architecture of their tissues of origin are the next generation of medical research models. The need for a more accurate prediction of a patient's response to a certain treatment has become key in the development of organoids. The ability to keep stem cells alive outside of the body allowed these organoids or mini-organs to be passaged, mainly in a three-dimensional matrix and never through plastic, almost indefinitely in some cases. Commonly, in order to passage or harvest these embedded organoids, proteases are employed to degrade these extracellular proteins. However, proteases also degrade proteins on the cell surface, and protease activity may carry over into subsequent cultures or lysate preparations.

Cultrex® Organoid Harvesting Solution provides a non-enzymatic method for depolymerizing extracellular



Organoid Harvesting Solution, 100 ml presentation. Catalog number 3700-100-01

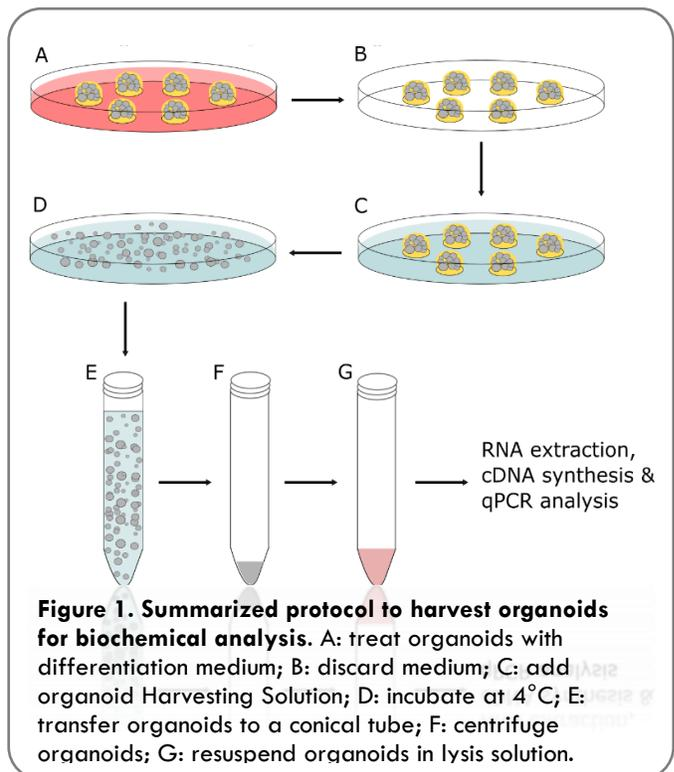
matrix proteins to allow for harvesting of intact organoids for passaging, cryopreservation, or biochemical analysis. Cultrex® Organoid Harvesting Solution is a non-enzymatic solution and is compatible with the main biochemical analysis techniques, such as quantitative PCR (qPCR) and western blotting. In this report, the Organoid

Resource Lab (ORL) provides you with a protocol to harvest organoids and analyze changes in gene expression by qRT-PCR.

Protocol

Briefly, Mouse Small Intestine organoids (mSI) were differentiated by removing Wnt3A in the culture medium and adding DAPT, a notch inhibitor. After several days,

the organoids were harvested, total RNA was isolated, and after cDNA synthesis, the expression of several markers of enteric differentiation was measured by qPCR.



Mouse Small Intestine (mSI) Organoid culture. Cultrex® Organoid Progenitor Cells: Mouse Small Intestine (catalog number 3750-001-01) were thawed, and DMSO from the freezing solution was removed by centrifugation. The cell pellet was resuspended in Cultrex® BME-R1 (catalog number 3433-005-R1) and cultured in mSI organoid culture medium for at least 3 passages. The composition of mSI organoid culture medium was the following: 50% of the final volume was completed with L Wnt3A conditioned medium, 1X B27 supplement, 1X N2 supplement, 2 mM GlutaMAX-I, 10 mM HEPES, 10 mM Nicotinamide, 1 mM N-Acetylcysteine, 10 nM [Leu15]-Gastrin I Human, 1 mg/ml HA-R-Spondin1-Fc (produced using Cultrex® Rspo1 cells, catalog number 3710-001-K), 100 ng/ml Recombinant Mouse Noggin, 500 nM A83-01, 7.5 µg/ml Human

Insulin, 10 μM SB 202190, 10 $\mu\text{g}/\text{ml}$ Human Transferrin, 50 ng/ml Recombinant Mouse EGF, Advanced DMEM/F-12 Cell Culture Medium was added to complete to final volume.

Differentiation of mSI organoids. In Fig.1, A, organoids were differentiated. The composition of the differentiation medium was identical to the organoid culture medium with the exception that it did not contain L Wnt3A and DAPT (a Notch inhibitor) was added fresh every medium change at a final concentration of 5 μM . The difference in volume was adjusted by adding Advanced DMEM/F12 Cell Culture Medium. For each qPCR data point, 6 domes of BME-R1 containing mSI organoids were arranged in a well of a 6-well tissue culture plate (each dome was composed by 50 μl of matrix). To detect significant morphological changes mSI organoids had to be maintained and passaged for several weeks in differentiation medium, however, to perform qPCR of enteric markers mSI organoids were maintained in differentiation medium for only 2 or 7 days.

Organoid harvesting, RNA extraction, and cDNA synthesis. In Fig.1, B - G the organoids are harvested and processed. In order to extract RNA and analyze gene expression, organoids were harvested using Cultrex[®] Organoid Harvesting Solution (catalog number 3700-100-01). Culture or differentiation medium was discarded, and the wells were washed with 5 ml of cold (4°C) PBS and incubated for 30 to 60 minutes with 5 ml of cold (4°C) Organoid Harvesting Solution. During this time, the plates were placed inside a container with ice or in a cold room with gentle shaking in order to achieve matrix depolymerization. Once the matrix was dissolved and the organoids were released, the solution was transferred to a conical centrifuge tube and centrifuged for 5 minutes at 500 x g at 4°C. The supernatant was discarded. The organoid pellet was resuspended in 1 ml of TriZol[™] (Thermo Fisher). Total RNA from organoids was extracted following manufacturer's instructions, and approximately 500 ng were used to synthesize cDNA from each sample using iScript[™] cDNA synthesis kit (Bio-Rad). Alternatively, the organoid pellet can be resuspended in a different lysis buffer depending on the application.

Quantitative PCR. Each cDNA sample was diluted 10 times with nuclease-free water and used as template for a quantitative PCR reaction. Each reaction had a final volume of 25 μl and contained 12.5 μl of 2X iQ[™] SYBR[®] Green Supermix (Bio-Rad), 0.5 μl of forward and reverse qPCR primer mix 500 nM, 9.5 μl of

nuclease-free water and 2.5 μl of diluted cDNA. Each sample was measured in triplicate using AB StepOnePlus thermocycler. All primer pairs were ordered from IDT DNA technologies. Beta-2 microglobulin was used as internal control.

Results

In Figure 2 qPCR results are shown. Lgr5, a G-protein coupled receptor, expressed mainly in stem cells is reduced after only 2 days of differentiation. This result is expected as the stem cell population decreases, and new crypts and villi start to form. On the other hand, the expression of Defa5, Muc2, and Tff3, all markers of intestinal differentiated cell types^{1,2} was increased over time.

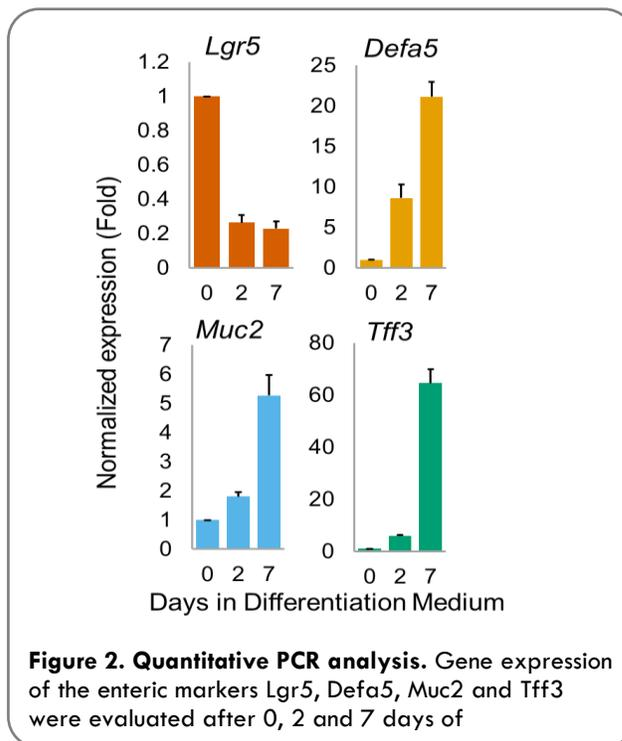


Figure 2. Quantitative PCR analysis. Gene expression of the enteric markers Lgr5, Defa5, Muc2 and Tff3 were evaluated after 0, 2 and 7 days of

References

1. VanDussen, K. L. *et al.* Development of an enhanced human gastrointestinal epithelial culture system to facilitate patient-based assays. *Gut* **64**, 911-920, doi:10.1136/gutjnl-2013-306651 (2015).
2. Yin, X. *et al.* Niche-independent high-purity cultures of Lgr5⁺ intestinal stem cells and their progeny. *Nat Methods* **11**, 106-112, doi:10.1038/nmeth.2737 (2014).

Contact Us

Trevigen, Inc.
8405 Helgerman Court
Gaithersburg, MD 20877
301-216-2800
info@trevigen.com
www.trevigen.com