

The Application of 3D Alginate Scaffolds for Liver Tissue Modeling

Alexandria Sams, Zhensheng Li, Mark Powers

Invitrogen Corporation • 1600 Faraday Avenue • Carlsbad, CA 92008 • USA

INTRODUCTION

Mammalian cell culture platforms have traditionally focused on optimizing specific targets of interest such as growth, productivity, or protein expression. While such approaches have enabled fundamental biological discoveries over the past several decades, they are typically incompatible with replicating *in vivo* cellular function. It is thought that this is in large part due to the inability of such systems to provide cultured cells with the proper physiological context. Recapitulating key aspects of the hierarchical environments in which cells reside *in vivo* -- mechanical, chemical, spatial, architectural, etc. -- has been shown to allow cells to function in a more functionally relevant manner. Approaches include cellular co-culture, perfusion bioreactors, and architecturally and chemically relevant biomaterials. Within our group, the design and development of extracellular matrices and scaffolds for the growth of cells in 3D is a particular area of focus.

One such material is AlgMatrix™, a lyophilized alginate sponge scaffold that drives intercellular interactions through the architectural properties of the interstitial pores. This novel scaffold is a platform technology to facilitate the creation of multicellular spheroidal structures: because cells do not significantly interact with alginate, they are able to interact more readily with one another. The technology is of particular interest in the field of liver biology, as hepatocellular spheroids have been shown to maintain expression of key hepatic markers, but are traditionally somewhat cumbersome to culture and maintain. For the purposes of these studies, primary rat hepatocytes were cultured in 24-well AlgMatrix™ plates and assessed for cell morphology, viability and functional parameters relative to standard culture formats.

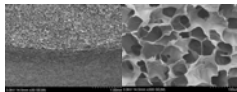
MATERIALS & METHODS

Cell culture: Briefly, hepatocytes (primary Sprague-Dawley rat hepatocytes or C3A human hepatoblastoma cells) were seeded by pipetting a concentrated cell suspension on top of the dry AlgMatrix™ scaffold. Following the complete absorption of media into the scaffold an additional volume of culture medium was added to each well. C3A cells were cultured in serum-containing media. Primary rat hepatocytes were cultured in serum-free media. Seeding concentrations and culture periods are detailed in each figure caption. All cultures were incubated in 5% CO₂ and 37°C.

For Collagen-Geltrex sandwich cultures, 24-well plates were coated with 30µg/mL collagen (Invitrogen, Carlsbad, CA). Primary rat hepatocytes were seeded (0.375×10⁶ cells/well). At ~5h post seeding serum-free media containing Geltrex™ (0.35mg/mL) (Invitrogen, Carlsbad, CA) was added to the cells. Cultures were incubated in 5% CO₂ and 37°C with daily medium changes.

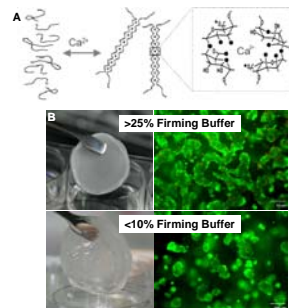
Caspase-3 Activity Determination: The quantification of caspase-3 proteolytic activity was assessed using a commercial kit (ApoTarget™ Caspase-3 Fluorometric Protease Assay, Invitrogen, Carlsbad, CA). The substrate is composed of a caspase-3 cleavage site and a fluorophore (ex: 400nm, em: 505nm).

FIGURE 1. Scaffold Properties



Alginate is a polysaccharide extracted from seaweed. AlgMatrix™ scaffolds are made from alginate cast in multiwell plates and lyophilized to optimize pore formation. These SEM images of sponges removed 96 well plates reveal interconnected pores with diameters ranging from 50 to 150µm. Cells residing in the inert alginate pores are able to interact with one another (rather than with the alginate biomaterial) to form 3D cellular structures.

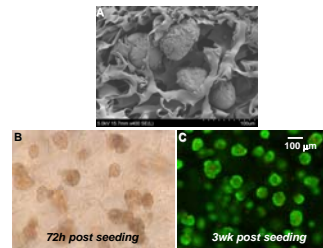
FIGURE 2. Tunable & Dissolvable



A) Alginate is a linear unbranched polymer composed of 1,4-linked β-D-mannuronic and α-L-guluronic acid residues. Ca²⁺ ions cause polyguluronic regions to dimerize to form hydrogels. B) AlgMatrix™ Firming Buffer is a calcium based solution that is applied prior to or concurrent with cell seeding. The ratio of firming buffer to medium in the solution dictates the firmness of the sponge and allows for tunable control of compliance/stiffness (an important factor in dictating cellular phenotype). AlgMatrix™ Dissolving Buffer is a sodium citrate solution that chelates calcium to liberate cells from sponges.

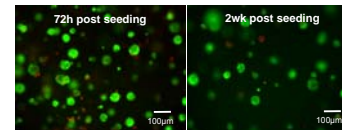
RESULTS

FIGURE 3. Spheroid Formation



C3A (human hepatoblastoma) cells were seeded (2×10⁶ cells/well) into a 6-well AlgMatrix™ plate and maintained for 3 wks with medium changes every 72h. A) SEM image illustrates coalescence of cells to form compact multicellular spheroids. B) Phase-contrast image shows C3A spheroid formation within 72h post cell seeding. C) Live/dead staining (2µM Calcein AM, 4µM Ethidium homodimer-1) indicates a significant # of viable hepatocytes present at 3 wks in culture.

FIGURE 4. Long-term Viability of Primary Hepatocytes



Freshly-isolated primary rat hepatocytes (Sprague-Dawley) were seeded (1×10⁶ cells/well) in a 24-well AlgMatrix™ plate and maintained with daily medium changes. Cell morphology was optimized at 10% Firming Buffer for use in primary hepatocyte applications. Live/dead staining (8µM Calcein AM, 4µM Ethidium homodimer-1) indicates that the spheroids contain a significant fraction of viable cells, through at least 2 wks of culture.

SUMMARY

- AlgMatrix™ is a chemically defined, highly porous (>90%) 3D scaffold; the hydrophilic nature of the material allows for rapid and efficient cell seeding and distribution.
- AlgMatrix™ stiffness can be controlled by the use of Firming Buffer (a calcium based solution). Cell recovery from AlgMatrix™ is achieved with the use of Dissolving Buffer, a non-enzymatic solution which dissolves the scaffold within a few minutes but leaves the cellular aggregates intact for further processing and/or analysis.
- Using various cell types we have demonstrated that cell aggregation within AlgMatrix™ leads to spheroid formation that is restricted by the fixed pore sizes (50-150µm) within the scaffold thereby preventing the development of hypoxic cores.
- We have demonstrated long-term viability using primary rat hepatocytes and C3A hepatoma cells. Cellular toxicity measurements demonstrate that the AlgMatrix™ culture environment promotes a healthier phenotype than the conventional sandwich culture.
- Our functional studies pertaining to CYP450 1A and 2B activity illustrate the superior capability of AlgMatrix™ to preserve liver-specific function and response in comparison to the conventional sandwich culture.

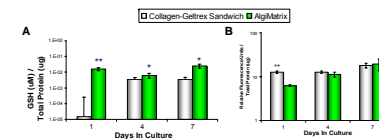
CONCLUSIONS

We have created a well-defined porous alginate matrix (AlgMatrix™) comprising interconnected pores that act as localized compartments in which cells form reproducible spheroidal aggregates. Our liver biology studies demonstrate that this scaffold offers an efficient platform for the creation of functionally relevant hepatocellular spheroids relative to conventional culture methods. Cellular toxicity measurements (intracellular glutathione levels, caspase-3 activity) demonstrate that the AlgMatrix™ culture environment is less stressful for hepatocytes in comparison to conventional sandwich culture. Functional performance analysis reveals that primary hepatocytes in AlgMatrix™ can respond specifically and quantitatively to model CYP450 inducers. Collectively, these results suggest that this platform is a potentially useful predictive tool in xenobiotic metabolism and toxicology research. We conclude that 3D cell culture scaffolds such as AlgMatrix™ serve as a foundation for the creation of more physiologically relevant culture systems.

ACKNOWLEDGEMENTS

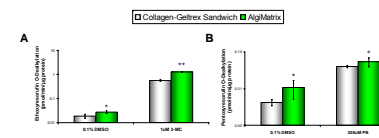
We wish to thank Ed LeCluyse, Stephen Ferguson, Cornelia Smith (Durham, NC), Jeanette Hill, and Jonathan Jackson (Austin, TX) for helpful discussions regarding primary hepatocyte work in AlgMatrix™.

FIGURE 5. Metrics of Cellular Stress: Primary Hepatocytes in AlgMatrix™ vs. Collagen-Geltrex™ Sandwich Culture



A) Intracellular levels of glutathione (GSH) are significantly higher for primary rat hepatocytes maintained in AlgMatrix™ throughout the culture period. Depletion of glutathione (GSH), a critical antioxidant, can lead to cell death. Therefore, GSH levels are used to assess toxicological responses. B) At the earliest time point (1 day post seeding) caspase-3 activity is significantly greater in hepatocytes seeded in the collagen-Geltrex™ sandwich culture. Caspase-3 is a key mediator of apoptosis. n=6. Two sample t-test, * denotes p<0.05, ** denotes p<0.001.

FIGURE 6. Functional Performance: CYP450 Activity



Among the CYP450 enzyme family the 1A and 2B subfamilies are particularly important with respect to drug & xenobiotic metabolism. A) EROD activity on day 5 post-seeding demonstrate that primary rat hepatocytes in AlgMatrix™ maintain significantly higher 1A activity (basal & induced expression). B) PROD activity on day 5 post-seeding demonstrates that primary rat hepatocytes in AlgMatrix™ maintain significantly higher 2B activity (basal & induced expression). n=6. Two sample t-test, * denotes p < 0.05, ** denotes p<0.001.