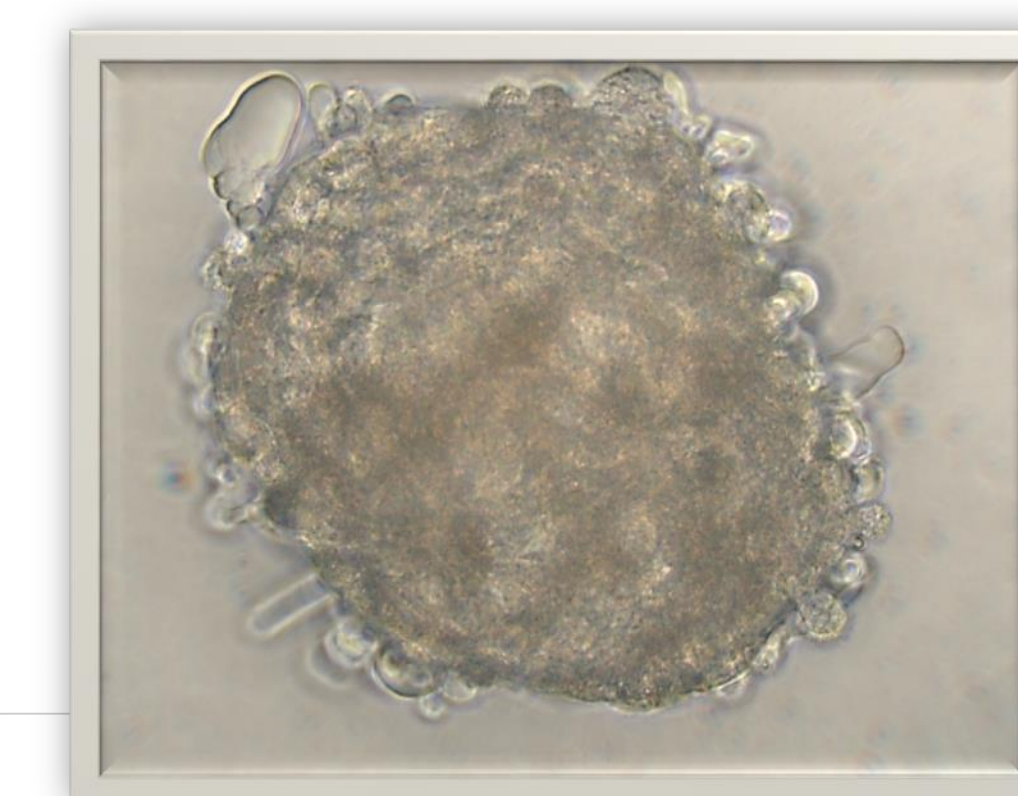


In-vitro evaluation of primary animal hepatic function using a 3D-spheroid culture system

Julia Tritapoe, Sujoy Lahiri, Michael Millet, Theresa Nguyen, Mark J. Powers and David Kuninger

Thermo Fisher Scientific, 7300 Governors Way, Frederick, Maryland, 21704



ABSTRACT

Recent evidence suggests that culturing primary hepatocytes in 3-dimensional (3D) hepatic spheroids offers major advantages in providing longevity and increased physiological relevance in modeling liver metabolism. Dog (Beagle), Mouse (CD-1), and Rat (Sprague Dawley) hepatocytes from Thermo Fisher Scientific were plated for spheroid qualification. 3D hepatic spheroids were characterized in weeks 1-4 in which the following were assessed: cell viability (as indicated by ATP synthesis), albumin production, phase I metabolic activity, and gene expression. Our data indicate that primary Dog and Rodent hepatocytes formed 3D spheroids by Day 5, and moreover were viable and functional for up to 4 weeks. In conclusion, our data indicate that animal 3D spheroid cultures are an accurate and sustainable *in vitro* model of hepatocyte function that maintain hepatic functions for a longer period of time in comparison with traditional 2D cultures.

INTRODUCTION

Conventional methods of culturing primary hepatocytes in 2-dimensional monolayer (2D) present limitations for the study of hepatic biology, liver function, and drug induced hepatotoxicity. Traditional 2D hepatocyte cultures rapidly de-differentiate resulting in the loss of hepatic specific function in approximately 5 days. While there are data to support that culturing primary human hepatic spheroids is a sustainable and robust *in vitro* model, limited information is available on the use of 3D animal spheroids in drug discovery. It is our hypothesis that growing animal hepatocytes in 3D spheroid cultures will more accurately reflect *in vivo* liver biology and maintain the liver functions for a longer period of time in comparison with the traditional 2D culture. Here we describe the incorporation of primary Dog and Rodent hepatocytes into 3D-culture systems that can be readily scaled for screening applications and studies where long term viability is important.

MATERIALS AND METHODS

Media and Reagents:

Plateable Dog (Beagle) (cat# DGCP10), Plateable Mouse (CD-1) Cryopreserved Hepatocytes (cat# MSCP10); Plateable Rat (Sprague-Dawley) Cryopreserved Hepatocytes (cat# RTCP10); Hepatocyte Plating medium – Supplemented Williams Media E, WEM (Gibco™ cat# A1217601) Hepatocyte Maintenance Medium – For Dog: Supplemented WEM (Gibco™ cat# A1217601) For Rodent: Supplemented DMEM Hepatocyte Thaw Medium - HTM (Gibco™ cat# CM7500) 3D Plates - Nunclon™ Sphera™ low attachment U-bottom 96-well Microplates (cat# 174925) 2D Collagen Plates (controls) – Gibco™ Collagen Coated 24-well plates (cat# A1142802)

Method

1. Cryopreserved Rat and Mouse Primary Hepatocyte (Gibco™) vials were thawed quickly in a 37°C water bath. Upon thawing the cells were promptly transferred into a 50 mL centrifuge tube containing HTM.
2. The tubes were centrifuged to pellet the cells. After centrifugation the supernatant was discarded.
3. The cell pellet was gently re-suspended in 1 mL hepatocyte plating media, followed by additional 3 mL. Cell counting was performed with Trypan blue using hemocytometer.
4. Appropriate volume of cell suspension was prepared to contain desired number of cells. Using a multichannel pipette 200uL of the cell suspension was added into individual well of Nunclon™ Sphera™ low attachment U-bottom 96-well Microplates. Cell numbers between 375 and 6000 cells/well were used.
5. The Sphera™ microplates were centrifuged at 200xg for 2 minutes to allow cells to group at the bottom of the wells. **NOTE:** Dog cells were centrifuged at 500g for 4 minutes.
6. The Sphera™ Microplates were transferred to an incubator (37° C, 5% CO₂, humidified) and allowed to sit undisturbed in Plating Media for 5 days. **NOTE:** Dog cells were centrifuged at 125g for 2 minutes on Days 1 and 2. (Figure 1)
7. By 5-days the hepatocytes formed spheroid with compact spherical mass of cells (Figure 1 and 2). On day 5, upon confirmation of spheroid formation, 50% the plating media was exchanged with hepatocyte maintenance media.

RESULTS

Figure 1. Work Flow of assembly and characterization of primary hepatocyte into 3D-spheroid

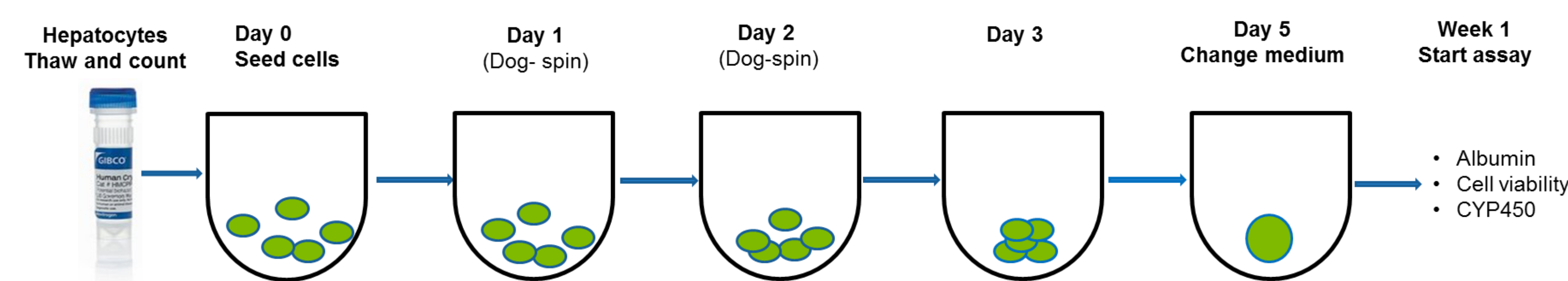
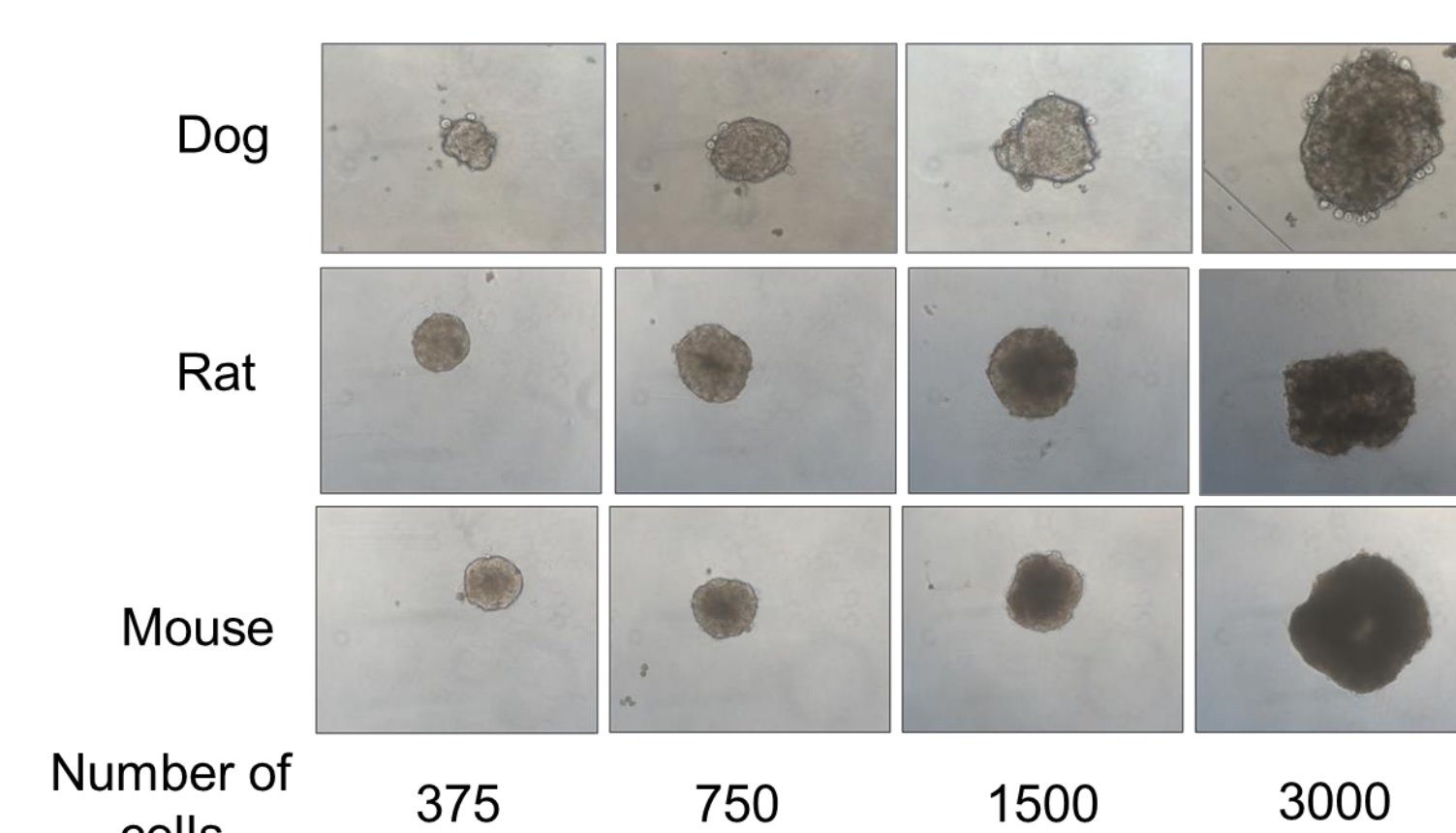
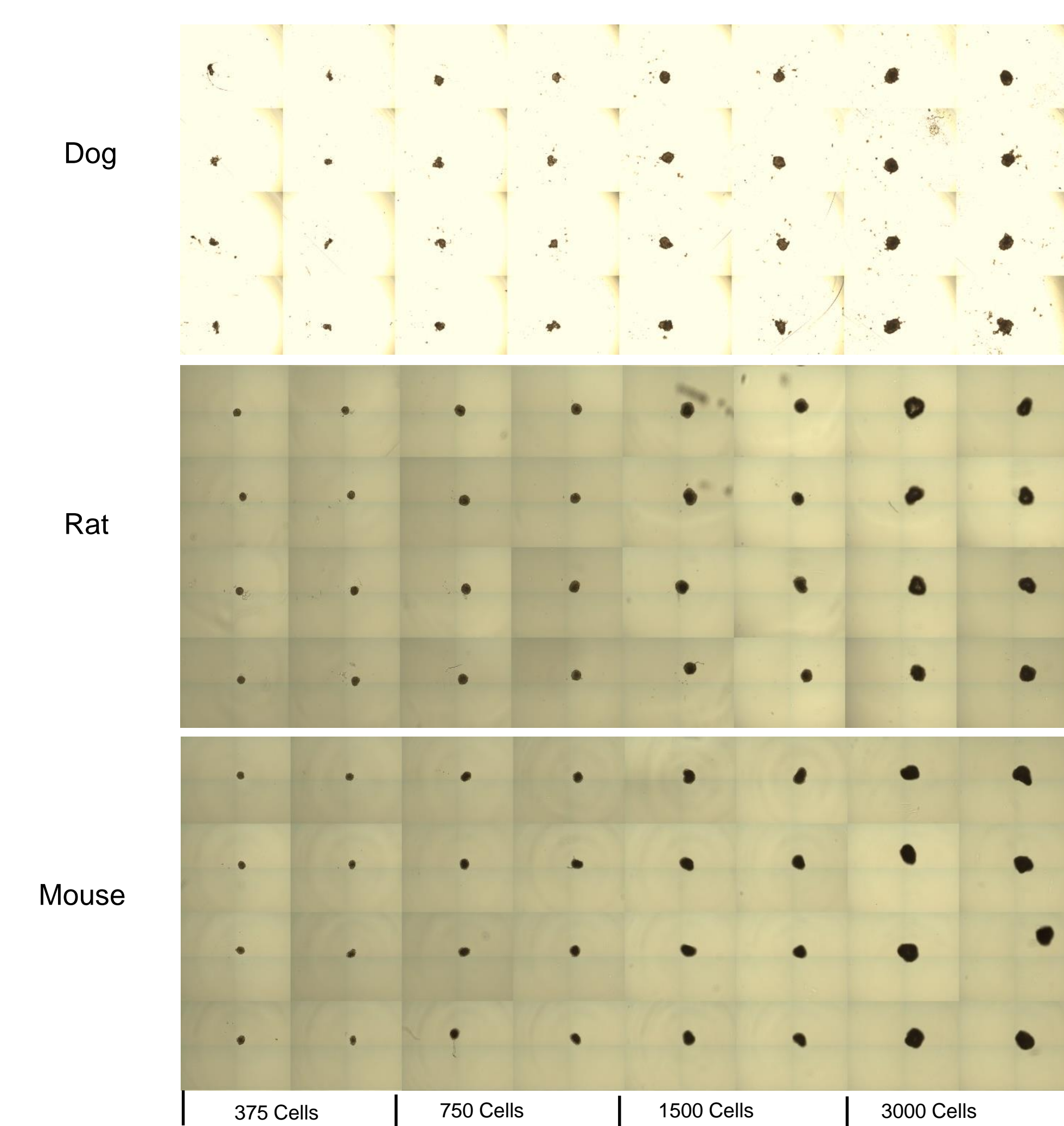


Figure 2. Dog and Rodent Hepatic Spheroids using Variable Cell numbers



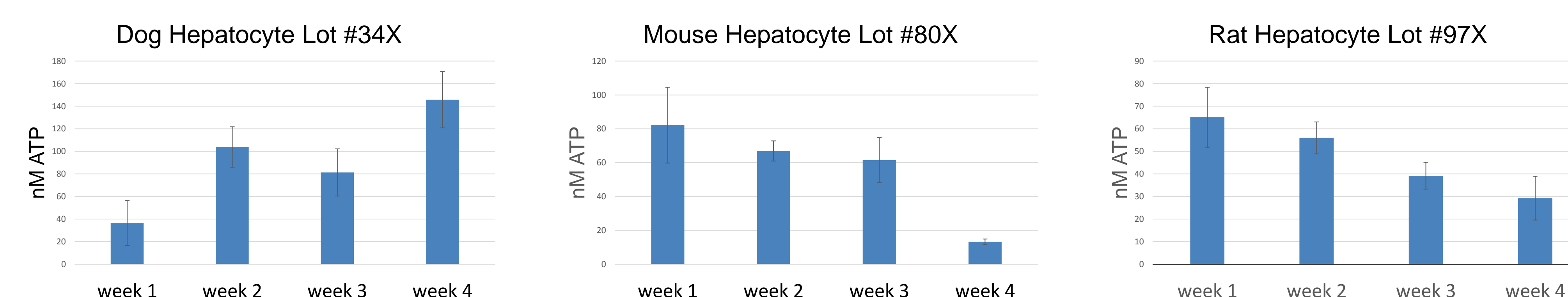
Dog and Rodent primary hepatocytes were seeded at different cell densities. During week 1 the spheroids were imaged using bright field microscopy at 10x optical zoom. Representative images from one lot per species are presented. **Data not shown:** Plating media supplemented with 20% serum was tested in the formation of dog spheroids, however this change did not impact the workflow outlined in Figure 1. or the results shown above.

Figure 3. Optimization of Hepatic Spheroids: Varying Cells/Well



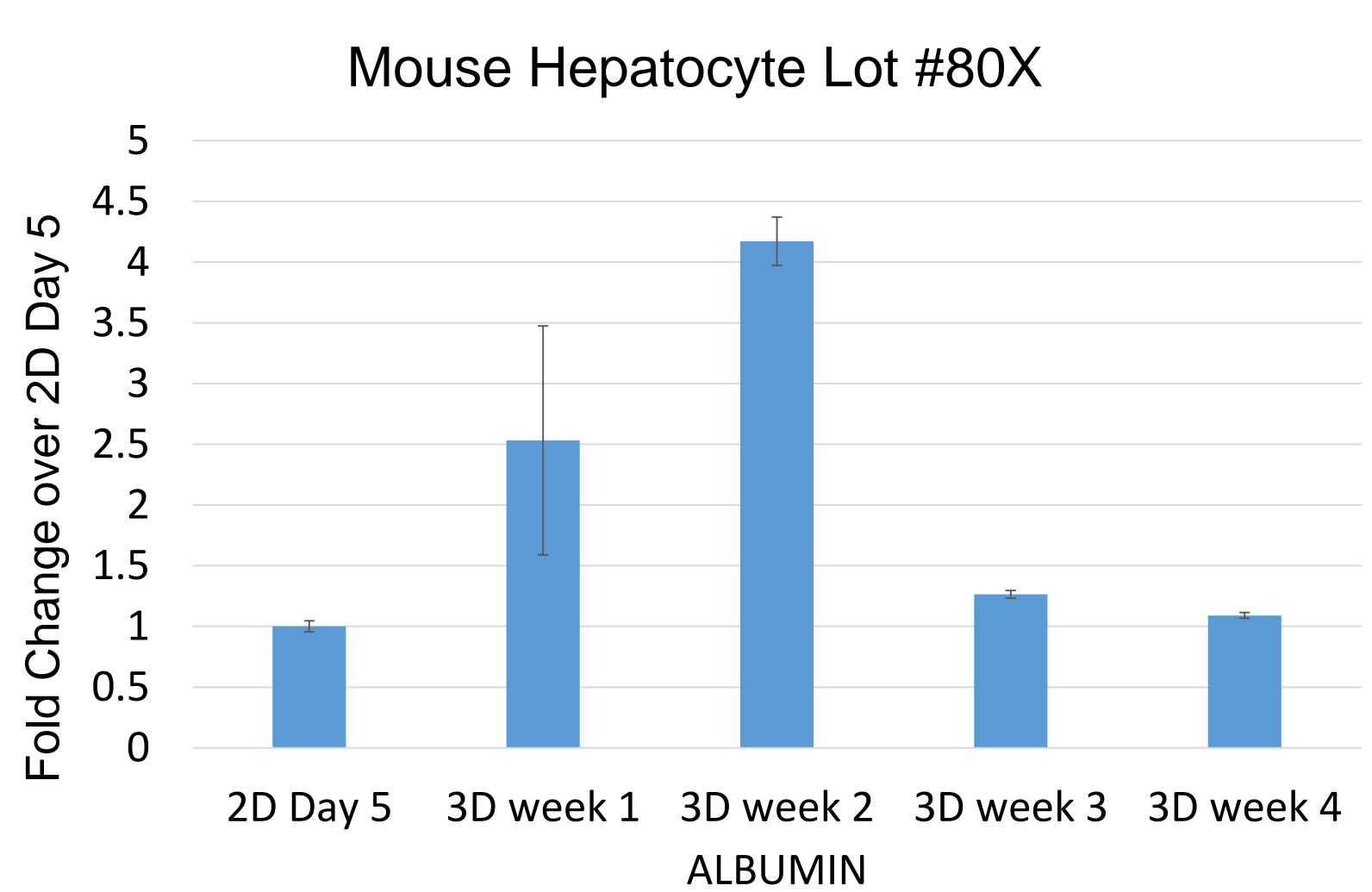
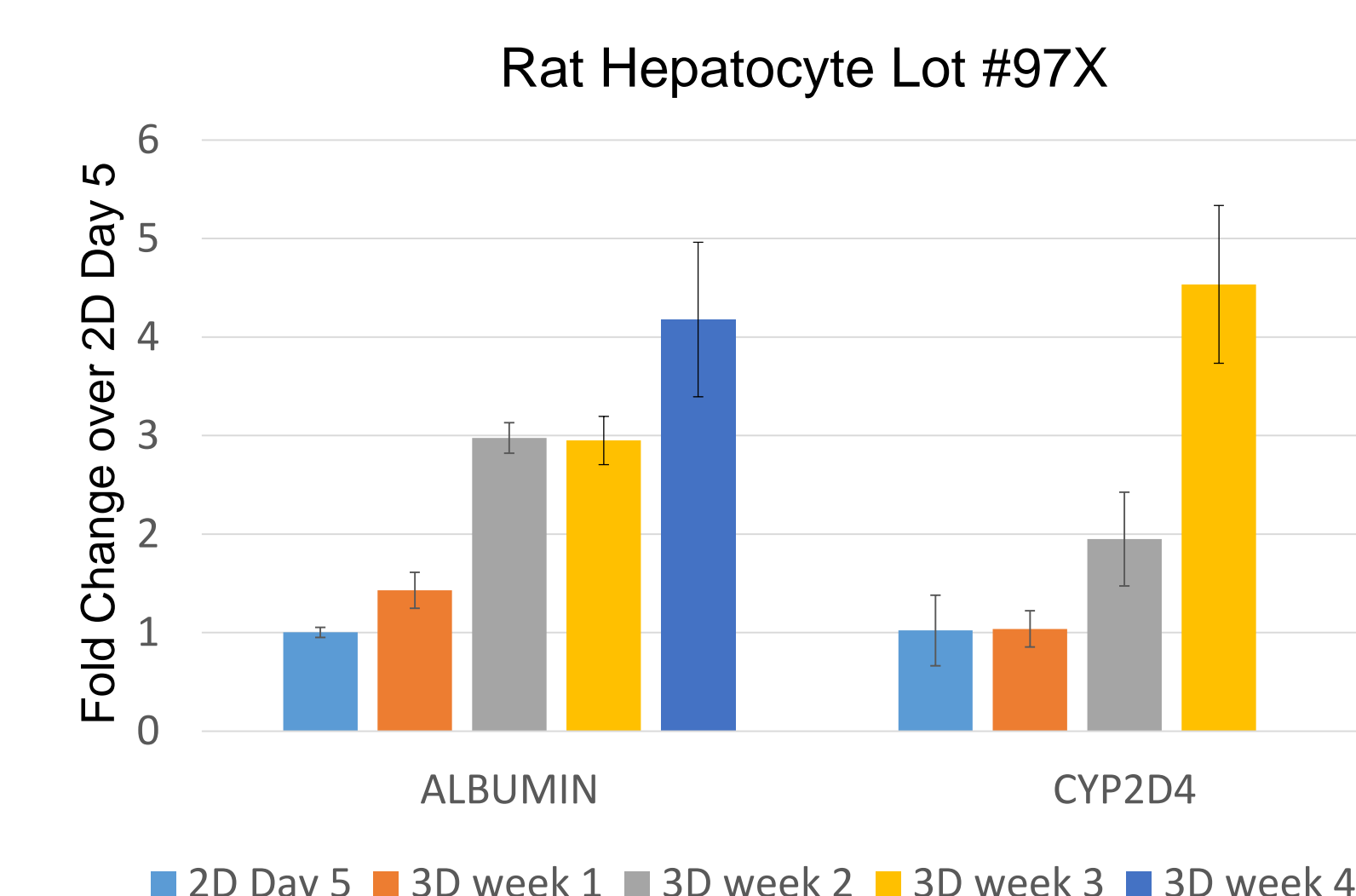
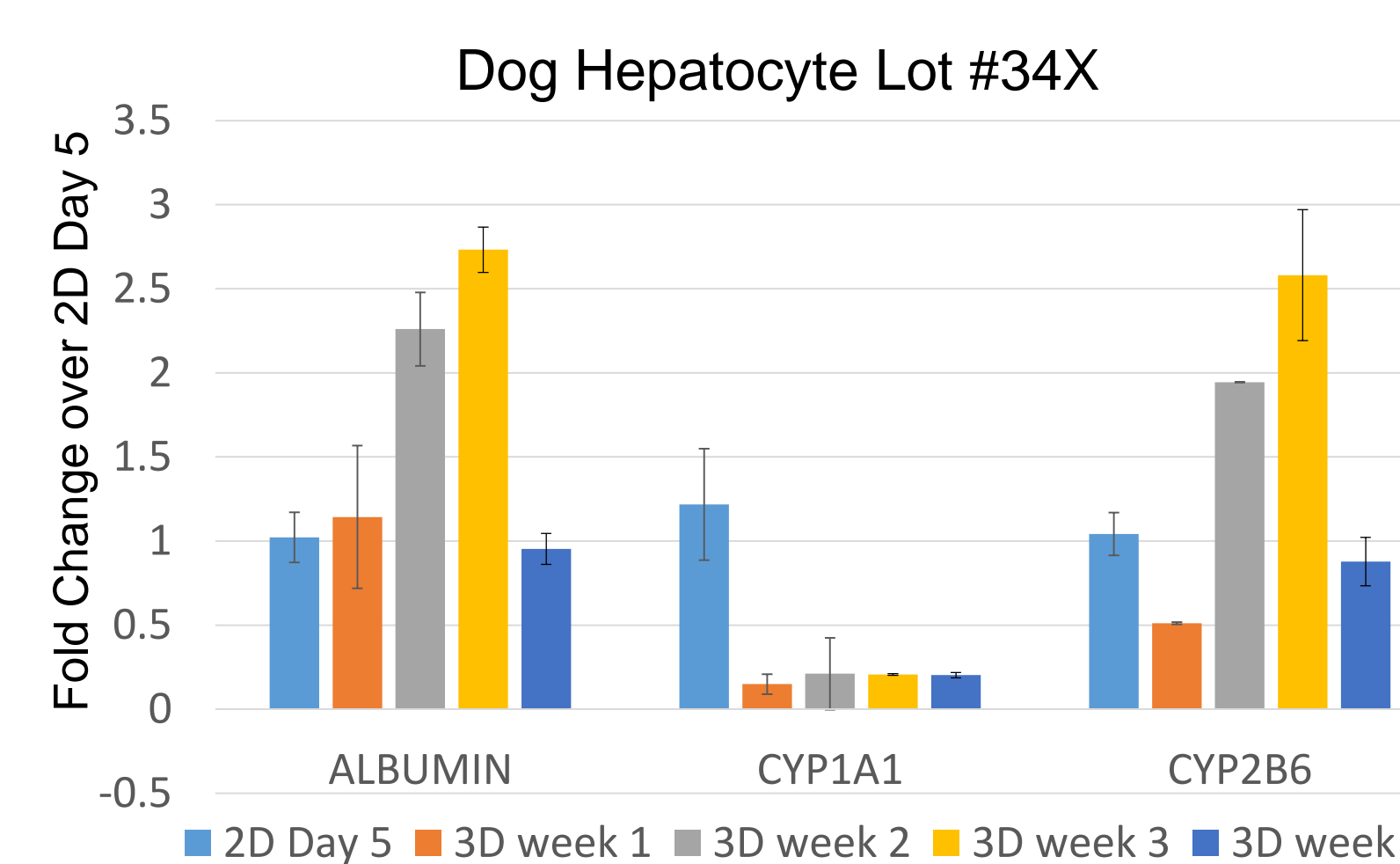
Primary Hepatocytes were seeded in Nunclon™ Sphera™ low attachment U-bottom 96-well Microplates in variable cell numbers.

Figure 5. ATP synthesis by 3D spheroids



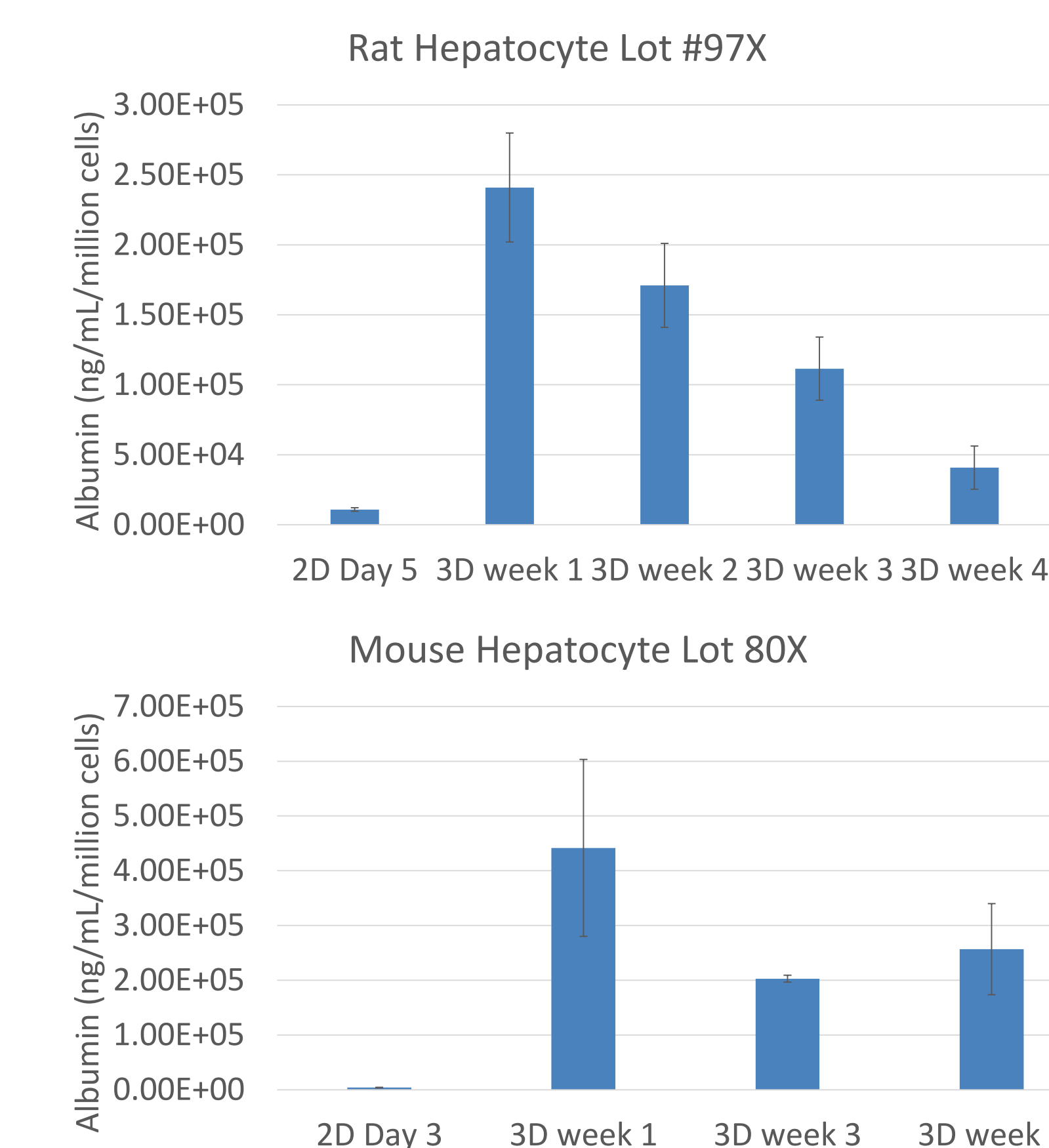
In weeks 1-4 of the spheroid culture the viability of the spheroids was assessed by measuring the ATP levels of 3 individual spheroids using a CellTiter-Glo® 3D cell viability assay (Promega). Results are mean +/- SEM, n = 3. **Data not shown:** modified parameter in ATP assay for Mouse and Rat Hepatocyte lots #80X and #97X with additional agitation, however this modification did not impact results shown above.

Figure 4. Comparison of Gene Expression between 2D and 3D cultures



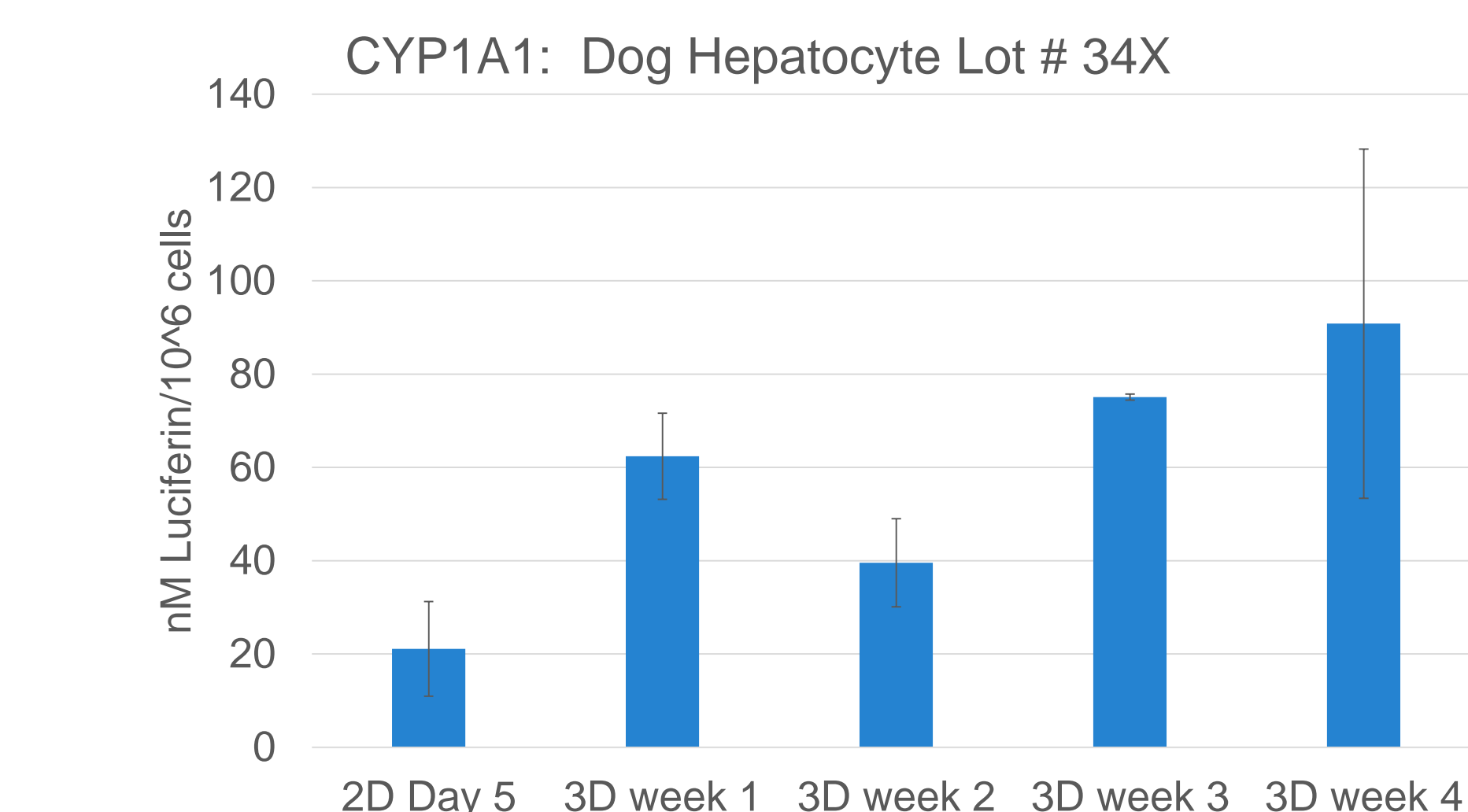
Gene expression levels were measured by qRT-PCR to compare the levels of gene expression between day 5 of 2D culture and various days of 3D culture.

Figure 6. Albumin Production



The concentration of albumin secreted was measured by an ELISA, (Abcam) and normalized to the total number of cells per well. Results are mean +/- SEM, n = 3.

Figure 7. CYP450 Activity



In weeks 1-4, 8 spheroids were pooled, CYP1A1 activity was measured using the P450 ProGlo assay (Promega). Results are mean +/- SEM, n = 2.

CONCLUSIONS

- Gibco™ Primary Dog, Mouse and Rat Hepatocytes can easily be assembled into a 3D culture in 5 days using Nunclon™ Sphera™ low attachment U-bottom 96-well Microplates, Gibco™ plating media & plating supplements and our recommended protocol.
- All animal hepatocyte lots tested for spheroid formation have formed spheroids. (2 lots of dog primary hepatocytes, 4 lots of mouse primary hepatocytes, and 3 lots of rat primary hepatocytes formed 3D-spheroids.)
- Dog spheroid formation requires additional spin steps on Days 1 and 2 for optimal results- these may be useful additional steps for lots refractory to spheroid generation.
- From data not shown- we've found multiple assays (CYP450 assays, ATP, others) or sample prep workflows designed for 2D that appear to require modification if they are to be useful for 3D applications.

FUTURE DIRECTIONS

- Explore alternative methods to assess viability and to measure CYP450 activity in animal spheroids.