

Technical Bulletin #420

Effects of ECM on Cell Proliferation and Differentiation in Hepatocytes

Paula Flaherty, NPD Scientist

BD Biosciences, Two Oak Park, Bedford, MA 01730

Abstract

The cellular microenvironment *in vivo* is characterized by specific interactions between cells, extracellular matrix (ECM), and soluble factors, such as cytokines and hormones. To create physiologically relevant *in vitro* models that support normal cell structure and function, the *in vitro* microenvironment has to incorporate the major components of the *in vivo* microenvironment. By optimizing the medium and ECM, we have constructed a cell culture system that supports the maintenance of differentiated hepatocytes for up to 3 weeks *in vitro*. Hepatocytes were isolated from adult male Wistar rats and plated in Hepato-STIM™ medium onto amorphous collagen I, collagen I gel, or BD Matrigel™ Basement Membrane Matrix. Cells were assessed for morphology (by light and electron microscopy), growth rates, and differentiated function (cytochrome P450 1A1 activity). Cells plated onto amorphous collagen I demonstrated a flattened, squamous morphology and rapidly lost function within 3-5 days. Hepatocytes on collagen I gels exhibited a mixed morphology of squamous and spherical cells and lost function within 7-10 days. Hepatocytes on BD Matrigel Matrix demonstrated a spherical, clustered morphology and maintained well-differentiated function for up to 22 days. Electron microscopy at 1 and 4 weeks demonstrated healthy, functioning cells with ultrastructural characteristics typical of differentiated hepatocytes. The cells contain a complex cytoplasm and are arranged in clusters of 6-12 cells, in a single layer enclosing a lumen. Adjacent cells have interdigitation of cytoplasmic extensions, tight and gap junctions as well as intercellular spaces suggestive of bile canaliculi. In conclusion, we have demonstrated that the combination of an optimized culture medium (Hepato-STIM medium) and ECM (BD Matrigel Matrix) provides an optimized microenvironment (BD BioCoat™ Hepatocyte Differentiation Environment) that supports long-term maintenance of differentiated function in hepatocytes *in vitro*.

Materials and Methods

Cell Isolation: Primary rat hepatocytes were isolated from adult, male Wistar rats (180-200 gms) following a two-step collagenase perfusion modified from published procedures^{1,2,3}. Cellular viability was assessed by trypan blue exclusion. Cells were seeded at 8×10^5 cells/well into 6-well plates containing BD Matrigel Matrix, collagen I, or a collagen I gel (BD Biosciences). Culture medium was Hepato-STIM medium supplemented with 50 u/ml penicillin, 50 µg/ml streptomycin, 100mM L-glutamine, and 10 ng/ml EGF (BD Biosciences). Cultures were incubated at 37°C, 5% CO₂-95% air, and medium was changed every 2-3 days.

Cell Proliferation: Cell growth was assessed at various time points using hemacytometer cell counts after a 60 minute Dispase treatment of cells grown on BD Matrigel Matrix or collagen I, and a 30 minute 0.05% Collagenase treatment of cells grown in a collagen gel.

Cytochrome Expression: Cytochrome P450 1A1 activity was measured using a modified metabolic fluorescent assay⁴. Briefly, cultured cells were incubated with a noncytotoxic concentration (8µM) of 7-ethoxyresorufin for 30 minutes. Dicumarol (10µM) was added to the assay media to prevent further metabolism of the resorufin formed by the cytosolic enzyme diaphorase. Culture supernatants were incubated with β-glucuronidase for 2 hours at 37°C, and dealkylated resorufin released was quantified by spectrofluorimetry using 530 nm excitation and 590 nm emission filters.

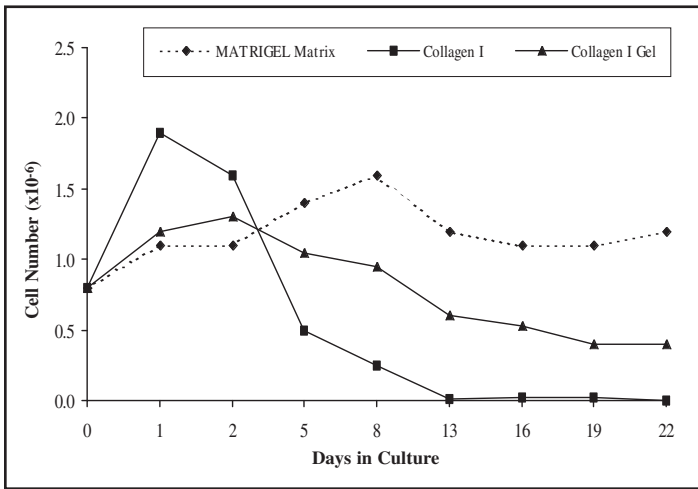
Light Microscopy: Samples were examined using an Olympus phase contrast microscope at a 10x magnification and photographed with Polaroid 667 film.

Scanning Electron Microscopy (SEM): Cells were fixed with 5% glutaraldehyde in a sucrose/cacodylate (0.1 M each) buffer, followed by 1% OsO₄ in buffer. Samples were dehydrated in graded ethanol, dried with HMDS, sputter-coated with gold palladium, and examined using an Amray 1800I microscope.

Transmission Electron Microscopy (TEM): Cells were fixed for 1.5 hours in K-II at 4°C, rinsed in 0.1 M cacodylate buffer, centrifuged, and suspended in 2% Bacto agar; 1 mm cubes were post-fixed in 2% OsO₄ in 0.1 M cacodylate for 2 hours. Samples were dehydrated in graded ethanol, infiltrated with propylene oxide and epon, and embedded in epon. Thick sections (1 µm) were stained with 0.5% toluidine blue (figure 5). Thin sections (70-90 nm) were stained with saturated uranyl acetate and Sato's lead stain (figures 6-9). Sections were examined and photographed using a Philips CM-10 microscope.

Results

Effect of ECM on hepatocyte proliferation



Effect of ECM on cytochrome P450 expression

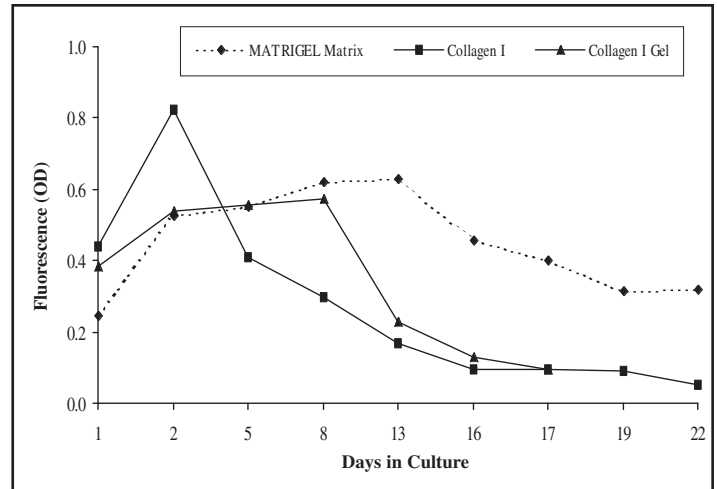


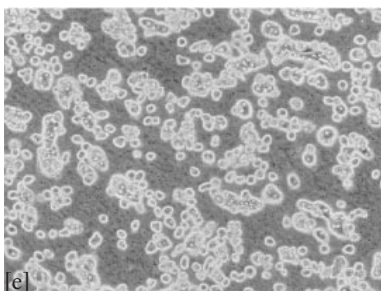
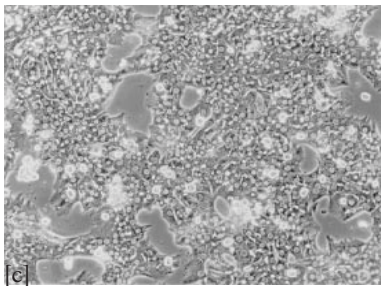
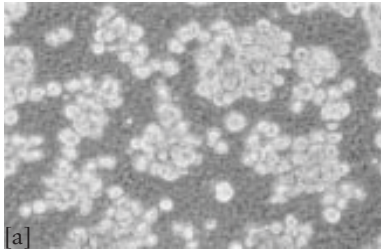
Figure 1: Effects of ECM on cell numbers. Note the decrease in the number of cells cultured on amorphous collagen I between days 2 and 13, and the gradual loss in the number of cells cultured in collagen I gels. In contrast, hepatocytes cultured on BD Matrigel Matrix maintain a constant cell number through day 22.

Figure 2: Effects of ECM on cytochrome P450 1A1 activity. Note the decrease in cytochrome P450 activity within 8-13 days for cells cultured on amorphous collagen I or in collagen I gels. In contrast, hepatocytes cultured on BD Matrigel™ Matrix maintain cytochrome P450 expression through day 22.

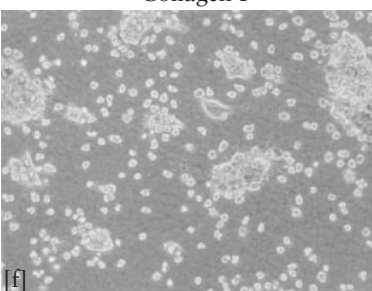
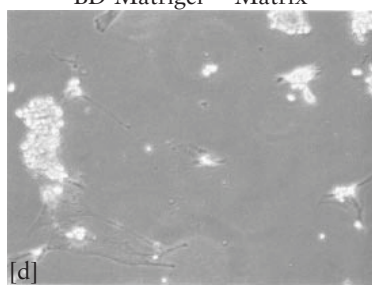
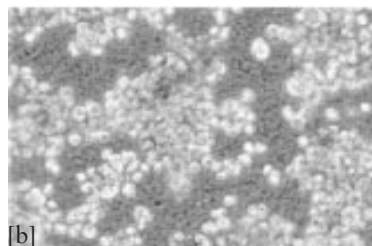
Effects of ECM on cell morphology: micrographs of hepatocytes cultured on various culture substrata

Light Micrographs

Day 1

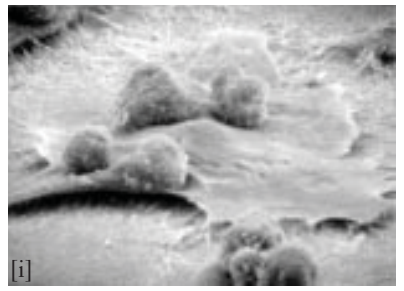
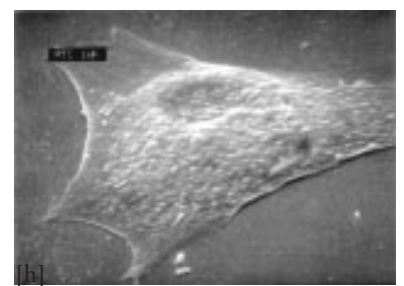


Day 21



Scanning Electron Micrographs

Day 2



BD Matrigel™ Matrix

Collagen I

Collagen I Gel

Figure 3: Light micrographs of primary rat liver epithelial cells cultured in a serum-free, supplemented medium for one day (a, c, e) or three weeks (b, d, f) on polystyrene plates containing BD Matrigel™ Matrix (BD BioCoat Hepatocyte Differentiation Environment) (a and b), collagen I (c and d), or collagen I gels (e and f). Note the lack of tight clusters of cells grown in collagen I gels, despite their spherical cell shape (magnification = 100x).

Figure 4: SEMs of hepatocytes cultured for two days on BD Matrigel™ Matrix (g), collagen I (h), or collagen I gels (i). Note the clusters of spherical cells for hepatocytes cultured on BD Matrigel™ Matrix, typical of differentiated cells (g).

Analyses of cell morphology by light and scanning electron microscopy demonstrate that a culture system, in which key components of the microenvironments (ECM and culture medium) have been optimized, can support a differentiated phenotype of primary rat hepatocytes. Furthermore, the results indicate that the combination of BD Matrigel™ Matrix and Hepato-STIM medium (forming the BD BioCoat Hepatocyte Differentiation Environment) outperforms other ECM and medium combinations examined. Using this system, differentiated cell morphology can be maintained for at least three weeks.

Transmission Electron Micrographs of Primary Rat Hepatocytes Cultured on BD Matrigel™ Matrix and on Optimized Serum-Free Medium

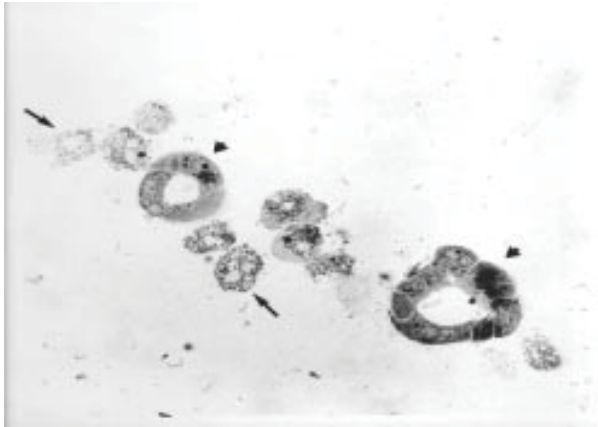


Figure 5: Thick sections of hepatocytes cultured on BD Matrigel Matrix in a serum-free medium (BD BioCoat Hepatocyte Differentiation Environment). Healthy cells are arranged in clusters (arrowheads) of 6-12 cells in a single layer (compare to figure 4a), enclosing a lumen, giving the appearance of a duct. In contrast, single cells (arrows) are dead or dying.

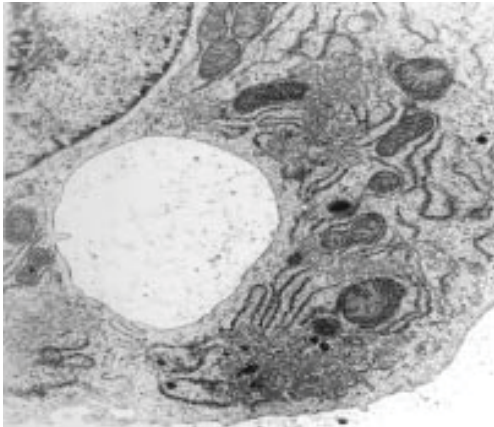
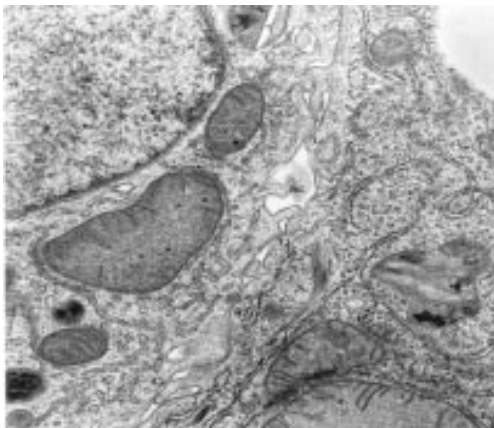


Figure 8 and 9: Thin sections of hepatocytes cultured in the BD BioCoat Hepatocyte Differentiation Environment for four weeks show healthy, functioning cells with frequent interdigitation of cytoplasmic extensions. Apposing cell membranes show tight junction as well as spaces or lumens between cells with microvilli present, suggestive of bile canaliculi.

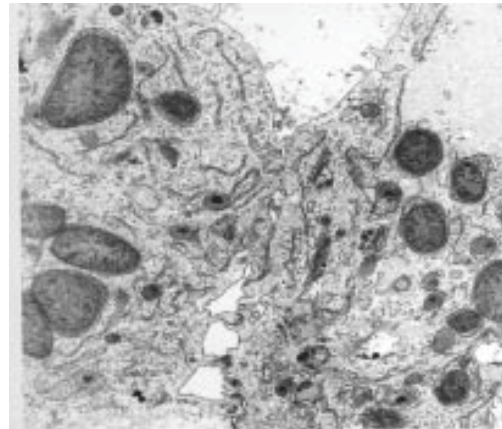
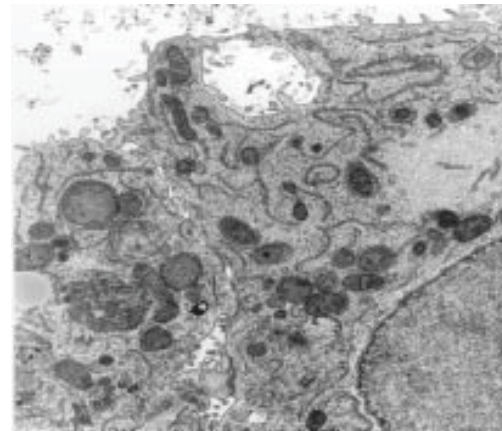


Figure 6 and 7: Thin sections of hepatocytes cultured for one week in the BD BioCoat Hepatocyte Differentiation Environment show healthy, functioning cells containing a complex cytoplasm, typical of differentiated cells. Nuclei are active with a single large nucleolus and the cytoplasm contains many mitochondria with few cristae and occasional calcium deposits as in normal hepatocytes. Rough ER is arranged in regular stacks of single strands, occasionally arranged near the lumen, outer cell surfaces, and surrounding mitochondria. The cytoplasm also contains golgi complexes, lysosomes, ribosomal rosettes, vesicles, lipid droplets, microfilaments, and microtubules. Hepatocytes contain large pools of flocculent material representing lost glycogen rinsed from tissue during dehydration.

Conclusions

- During a three-week culture period in the BD BioCoat Hepatocyte Differentiation Environment, the cell number of the hepatocytes changes only slightly indicating absence of cell proliferation (Figure 1), characteristic of differentiated cells^{5,6,7,10}.
- Cellular functions typical of differentiated hepatocytes are maintained for at least three weeks as shown by continued expression of cytochrome P450 (Figure 2), a hemoprotein involved in the oxidative metabolism of various compounds, an important function of the liver.
- Tight clusters of spherical cells with a lumen appear within the first day of culture in the BD BioCoat Hepatocyte Differentiation Environment. Cells will not spread, but remain in clusters throughout the entire culture period of three weeks (Figures 3a, 3b, and 4a).
- TEMs at one and four weeks of hepatocytes cultured in the BD BioCoat Hepatocyte Differentiation Environment show healthy functioning clusters of 6-12 cells in a single layer surrounding a lumen. Junctional complexes between apposing cell membranes are suggestive of bile canaliculi.

Discussion

The study of hepatocytes in culture is of great importance to research investigating mechanisms that control liver-specific functions. However, hepatocytes often dedifferentiate and die in response to inadequate culture environments, precluding such systems from being used as relevant *in vitro* models.

BD Matrigel™ Matrix, a reconstituted basement membrane, has been shown to suppress cell growth and prevent growth-associated dedifferentiation, as evidenced by the disappearance of growth-related transcription factors and appearance of liver-specific transcription factors in cells cultured on BD Matrigel Matrix⁸. Hepatocytes cultured in BD Matrigel Matrix-based systems have been shown to maintain liver-specific functions *in vitro* for longer periods of time than do hepatocytes cultured in most collagen-based systems⁹⁻¹². Hepatocytes cultured in BD Matrigel Matrix-based systems have been shown to express high levels of plasma proteins, such as albumin⁹ and the multidrug resistance-associated plasma membrane glycoprotein gp95¹⁰, and retain the ability to respond to a number of xenobiotics^{9,11} and hormones¹² with increased expression of members of the cytochrome P450 supergene family or metabolic enzymes, respectively.

Changes in cell shape as modulated by ECM have been

shown to regulate expression of growth and differentiation in rat hepatocytes^{8,13,14}. Culture of hepatocytes on collagen I induces expression of Immediate-Early Growth Response genes (Cjun, jun-B, C-myc) expressed in proliferating cells, and downregulates expression of transcription factors (e.g., C/EBPα) shown to be expressed in differentiated cells. In a collagen I sandwich, albumin mRNA expression is maintained, but C/EBPα expression is downregulated. In contrast, hepatocytes cultured on a reconstituted basement membrane (BD Matrigel Matrix) show restored C/EBPα expression and low-level expression of C-jun, jun-B, and C-myc, indicating that the cells are growth-arrested in GO phase^{8,14}.

The BD BioCoat Hepatocyte Differentiation Environment is an integrated culture system of BD BioCoat BD Matrigel Matrix plates (6-well) and a fully defined, serum-free optimized medium that is supplemented with hormones (dexamethasone and insulin), growth factors (EGF), and other defined metabolites. We have shown that this system has been optimized to support cultures of differentiated hepatocytes in a serum-free environment for at least 3 weeks. Due to its ability to support a liver-specific phenotype, this Environment can be used to create *in vitro* models to study various liver functions, such as liver-specific gene expression, metabolism and transport of drugs or natural compounds, drug toxicity, and altered or impaired liver function due to disease.

References

1. Berry, M.N. and Friend, D.S., *J. Cell Biol.* 43:506 (1969).
2. Rana, B., et al., *Mol. Cell. Biol.* 14:5858 (1994).
3. Seglan, P.O., *Metho. Cell Biol.* 13:29 (1976).
4. Schuetz, E.G., et al., *J. Cell. Physiol.* 134:309 (1988).
5. Aiken, J., et al., *J. Ped. Surg.* 25:140 (1990).
6. Schuetz, J.D. and Schuetz, E.G., *Cell Growth & Diff.* 4:31 (1993).
7. Donato, M.T., et al., *Anal. Biochem.* 213:29 (1993).
8. Kane, R.E., et al., *In Vitro Cell Dev. Biol.* 27A:953 (1991).
9. Bissell, D.M., et al., *J. Clin. Invest.* 79:801 (1987).
10. Mann, D.J., et al., *J. Mol. Endocrin.* 8:235 (1992).
11. Ben-Ze'ev, A., et al., *Proc. Natl. Acad. Sci. USA* 85:2161 (1988).
12. Hansen, L.K., et al., *Mol. Biol. Cell* 5:967 (1994).
13. DiPersio, C.M., et al., *Mol. Cell. Biol.* 11:4405 (1991).
14. Bucher, N.L.R., et al., *Mol. Biol. Cell* 5:312a (1994).

BD Biosciences

Two Oak Park
Bedford, MA 01730 USA
tel: 877.232.8995
fax: 800.325.9637

BD

2280 Argentia Road
Mississauga, Ontario
Canada L5N 6H8
tel: 866.979.9408
fax: 800.565.0897

BD

Akasaka Garden City,
Akasaka 4-15-1, Minato-ku,
Tokyo, 107-0052 Japan
tel: (81) 24 593 5405
fax: (81) 24 593 5761

BD Biosciences

Singapore Branch
30 Tuas Avenue 2
Singapore 639461
tel: (65) 6861 0633
fax: (65) 6860 1590

BD Biosciences

Erembodegem-Dorp 86
9320 Erembodegem, Belgium
tel: (32) 53 720 211
fax: (32) 53 720 450
contact_bdb@europe.bd.com

BD Biosciences

4 Research Park Drive
Macquarie University Research Park
North Ryde NSW 2113 Australia
tel: (800) 656 100
fax: (612) 8875 7200
aus_customerservice@bd.com