

**Technical notes for using 3D Biotek's PCL and PS scaffolds**

Dear valued customer,

Our 3D Insert products are so new that you may be the first researcher to use them for culturing the cell types that interest you. We kindly appreciate any comments and experiences that you can share with us after using our scaffolds in your experiments.

Before you begin, we would like to share some experiences with you:

1. When seeding cells onto **3D Biotek's PCL and PS scaffolds**, take care to ensure that the cell suspension is pipetted directly onto the center of the scaffolds. Cell suspensions that contact the sides of tissue culture wells can significantly lower cell seeding efficiency.
2. A benefit of using **3D Biotek's PS scaffold** is the ability to monitor cell growth and morphology in real time with an inverted light microscope. Be sure to take advantage of this by simply changing the focus plane on a light microscope to view cells cultured on 4 different layers throughout the scaffold.
3. We have found that cells cultured on **3D Biotek's PCL and PS scaffolds** secrete significant extracellular matrix (ECM), similar to cells *in vivo*. Please keep this in mind when detaching cells for subsequent assays, since a longer trypsin incubation period or an alternative cell detachment protocol may be needed for some cell types.

For further inquiries, please contact technical support at:

E-mail: [techsupport@3dbiotek.com](mailto:techsupport@3dbiotek.com)  
Tel: 732-729-6270 (ext. 4108)