

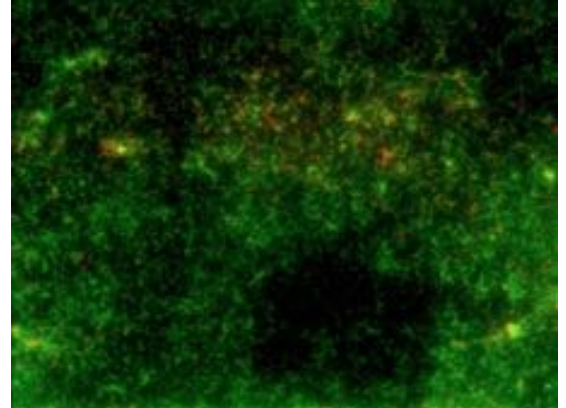
Biofilm Research *accelerated*

Unique Assays and Dramatic Productivity Gains

BioFlux™ Systems

BioFlux is a unique microfluidic platform delivers, going beyond what is possible with traditional systems. Key advantages:

- Simultaneously grow up to 96 biofilms under controlled shear conditions
- Provide controlled media flow for up to 21 days
- Drastically reduce media consumption:
 - Reduce use of expensive reagents like stains and drug compounds
 - Grow oral biofilms using saliva as media
- Use 2-inlet designs to perform:
 - Multiple parallel bacterial chemotaxis experiments
 - Multiple parallel competitive biofilm development assays
- Control temperature, gas content, nutrient delivery, even grow anaerobic biofilms
- Accurately control shear forces to quantify bacterial adhesion
- Quick and easy setup with no messy pumps or tubing



Biofilm Research *accelerated*

Unique Assays and Dramatic Productivity Gains

Over a hundred labs and biotech companies are using the BioFlux to perform world-class research. It offers the biological relevance of a laminar flow cell, with the throughput and convenience of standard microplates. The system comes complete with all the tools necessary to generate your data quickly and reliably.

Learn more about BioFlux

- **Product Information**
fluxionbio.com/bioflux
- **Applications**
fluxionbio.com/bioflux-applications
- **Live Demos**
fluxionbio.com/livedemos
- **Webinars**
fluxionbio.com/bioflux-webinars
- **Application Notes & Publications**
support.fluxionbio.com



© 2016 Fluxion Biosciences, Inc. All rights reserved.



Toll Free (USA): +1 (866) 266-8380
Phone: +1 (650) 241-4777
Email: sales@fluxionbio.com