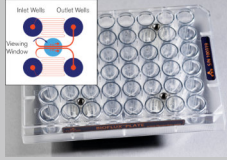
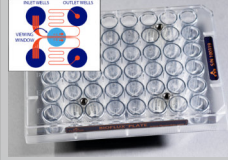
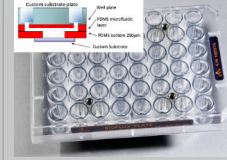
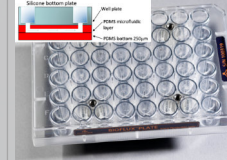

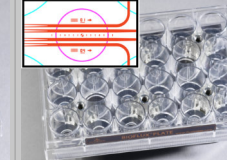



BioFlux Plate Selection Guide

	48-well 0-20 dyn/cm ²	48-well 0-200 dyn/cm ²	48-well 0-20 dyn/cm ² custom bottom	48-well 0-20 dyn/cm ² PDMS bottom	24-well 0-20 dyn/cm ²	24-well 0-20 dyn/cm ² PDMS bottom	6-well 0-20 dyn/cm ²
							
Throughput	24 experimental channels	24 experimental channels	24 experimental channels	24 experimental channels	8 experimental channels	8 experimental channels	3 experimental channels
Plate bottom	170µm cover slip glass	170µm cover slip glass	User-defined	250µm PDMS	170µm cover slip glass	250µm PDMS	170µm cover slip glass
Inlet / outlet well design	One inlet / one outlet	One inlet / one outlet	One inlet / one outlet	One inlet / one outlet	Two inlets / one outlet	Two inlets / one outlet	One inlet / one outlet
Well volume	1 mL	1 mL	1 mL	1 mL	3.5 mL	3.5 mL	17.7 mL
Shear range	Up to 20 dyn/cm ²	Up to 200 dyn/cm ²	Up to 5 dyn/cm ²	Up to 20 dyn/cm ²	Up to 20 dyn/cm ²	Up to 20 dyn/cm ²	Up to 20 dyn/cm ²
Maximum run time at 0.5 dyn	24.2 hours	24.2 hours	24.2 hours	24.2 hours	62.7 hours	62.7 hours	216.8 hours
Applications	Biofilms, standard cell adhesion, lower shear platelet/thrombosis assays	Vasculature modeling, thrombosis, vWF mediated platelet tethering	User-defined	Medical device infections, medical implants, environmental and industrial biofilms	Invasion, chemotaxis, competitive biofilms, drug dosing, wound healing	Invasion, chemotaxis, wound healing, drug screening for medical device infections	Long-term culturing of biofilms, leukocytes, platelets
BioFlux system compatibility	One / 200 / DCIS / 1000z	One / 200 / DCIS / 1000z	One / 200 / DCIS / 1000z	One / 200 / DCIS / 1000z	200 / DCIS / 1000z	200 / DCIS / 1000z	DCIS / 1000z
Part number	One / 200: 910-0004-5Pack DCIS / 1000z: 910-0047-5Pack	One / 200: 910-0033-5Pack DCIS / 1000z: 910-0048-5Pack	One / 200: 910-0119-5Pack DCIS / 1000z: 910-0120-5Pack	One / 200: 910-0115-5Pack DCIS / 1000z: 910-0121-5Pack	200: 910-0050-5Pack DCIS / 1000z: 910-0051-5Pack	200: 910-0116-5Pack DCIS / 1000z: 910-0145-5Pack	DCIS / 1000z: 910-0110-5Pack

Fluxion Biosciences, Inc.

Toll Free (USA): +1 (866) 266-8380

Main: +1 (650) 241-4777

Quotes: sales@fluxionbio.com

Orders: orders@fluxionbio.com

www.fluxionbio.com

