



Fluxion Biosciences
Ordering Information

2010

To place an order, send purchase order to:

sales@fluxionbio.com
 US Toll Free (866) 266-8380
 FAX (650) 873-3665

Product	Description	Fluxion Part #
Instruments	BioFlux 200 System - 24-well Interface, Controller, Heating Plate, Vapor Trap, BioFlux 200 Software	950-0012
	BioFlux 200 System - 48-well Interface, Controller, Heating Plate, Vapor Trap, BioFlux 200 Software	950-0010
	BioFlux 1000 System	<i>Please inquire</i>
	IonFlux-16 System with Software	950-0013
	IonFlux-HT System with Software	950-0014
Accessories	External Gas Module, BioFlux	950-0035
	Vapor Trap, BioFlux	900-0019
Well Plate Microfluidic™ Devices	BioFlux Plate, 24-well, 20dyne, 1/pack	910-0009
	<i>See pricing sheet or request quotation for quantity discounts</i>	
	BioFlux Plate, 48-well, 20dyne, 1/pack	910-0004
	<i>See pricing sheet or request quotation for quantity discounts</i>	
	BioFlux Plate, 48-well, 200dyne, 1/pack	910-0033
	<i>See pricing sheet or request quotation for quantity discounts</i>	
	Sterile film, BioFlux, 24-well, 50/pack	910-0011
	Sterile film, BioFlux, 48-well, 50/pack	910-0012
	IonFlux 16 Plate, 96 Well, 1 pk	910-0021
	IonFlux 16 Plate, 96 Well, 10 pk	910-0022
	IonFlux 16 Plate, 96 Well, 50 pk	910-0023
	IonFlux 16 Plate, 96 Well, 100 pk	910-0024
	IonFlux 16 Plate, 96 Well, 250 pk	910-0025
	IonFlux 16 Plate, 96 Well, 500 pk	910-0026
	IonFlux HT Plate, 384 Well, 1 pk	910-0027
	IonFlux HT Plate, 384 Well, 10 pk	910-0028
	IonFlux HT Plate, 384 Well, 50 pk	910-0029
	IonFlux HT Plate, 384 Well, 100 pk	910-0030
IonFlux HT Plate, 384 Well, 250 pk	910-0031	
IonFlux HT Plate, 384 Well, 500 pk	910-0032	
Preventive Maintenance and Replacement Parts	Tubing harness, Controller to Vapor Trap	200-0031
	Tubing harness, Vapor Trap to 48-well Interface	200-0032
	Tubing harness, Vapor Trap to 24-well Interface	200-0033
	Vapor Trap Filters, sterile, 16/pack	910-0017
	Gasket, 24 Well Interface	800-0018
	Gasket, 48 Well Interface	800-0019