



Fluxion Biosciences
Ordering Information
 2009

To place an order:
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	900-0011	
	Vapor Trap, BioFlux	900-0019	
	Power Cord, UK, 8 ft, CE	400-0063	
	Power Cord, Schuko, 8ft, CE	400-0064	
	Power Cord, Swiss, 8ft, CE	400-0065	
Well Plate Microfluidic™ Devices	BioFlux Plate, 24-well, 20dyne, 1/pack	910-0009	
	BioFlux Plate, 24-well, 20dyne, 5/pack	910-0006	
	BioFlux Plate, 24-well, 20dyne, 10/pack	910-0008	
	BioFlux Plate, 24-well, 20dyne, 25/pack	910-0007	
	BioFlux Plate, 24-well, 20dyne, 50/pack	910-0005	
	BioFlux Plate, 48-well, 20dyne, 1/pack	910-0004	
	BioFlux Plate, 48-well, 20dyne, 5/pack	910-0001	
	BioFlux Plate, 48-well, 20dyne, 10/pack	910-0003	
	BioFlux Plate, 48-well, 20dyne, 25/pack	910-0002	
	BioFlux Plate, 48-well, 20dyne, 50/pack	910-0000	
	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	