

2012 Parts Pricing – CN Parts

Part Number	DESCRIPTION	2012 Price
CN-0001-01	Air Filter Base (Hepa)	\$ 82.69
CN-0002-01	Base Air Filter	\$ 95.81
CN-0005-01	Serial Cable (5 Ft)	\$ 107.63
CN-0006-01	Short Sample Probe	\$ 185.06
CN-0007-01	Long Sample Probe	\$ 240.19
CN-0008-01	Sample Tube Holder (Large, 1.5 mL)	\$ 70.88
CN-0009-01	Sample Tube Holder (Small, 1.2 mL)	\$ 80.06
CN-0010-01	Sheath Filter w/Quick Disconnect	\$ 140.44
CN-0011-01	Sheath Bottle	\$ 273.00
CN-0012-01	Waste Bottle	\$ 257.25
CN-0013-01	Syringe Cylinder w/Seal	\$ 255.94
CN-0014-01	Syringe Seal (Qty 4)	\$ 124.69
CN-0015-01	Sample Needle Height Alignment Kit	\$ 135.19
CN-0016-01	Alignment Tube (XYP)	\$ 115.50
CN-0017-01	Heater Block (XYP)	\$ 173.25
CN-0018-01	USB Cable	\$ 102.38
CN-0019-01	Fuse 2AMP 250V Fast Acting (Qty 10)	\$ 72.19
CN-0022-01	Cleansing Reservoir (XYP)	\$ 112.88
CN-0023-01	V-Housing Shield	\$ 99.75
CN-0025-01	Allen Wrench (3/32)	\$ 76.13
CN-0027-01	Intake Air Filter (Qty 5)	\$ 70.88
CN-0028-01	Sheath Filter	\$ 82.69
CN-0031-01	Fan Guard (XYP)	\$ 66.94
CN-0032-01	DeBubbler Vented Cap	\$ 84.00
CN-0037-01	Solvent Inlet Filter (10 Micron)	\$ 114.19
CN-0048-01	Inline Sheath Filter (Quick Replace)	\$ 57.75
CN-0051-01	3A Fuse (250V, Fast Acting)	\$ 66.94
CN-0053-01	Air Filter (Front Mount)	\$ 59.06
CN-0092-01	SD Tubing Kit	\$ 287.44
CN-0096-01	PM Kit (6 Month)	\$ 135.19
CN-0118-01	Quad Ring Seals, Q4003, 70D EPDM (Set of 8)	\$ 36.75
CN-0119-01	Lee Valve to Injection Needle Tubing	\$ 110.25
CN-0144-01	Foam Lined Crate (LX100 HTS)	\$ 662.81
CN-0145-01	LX100 Shipping Box	\$ 216.56
CN-0146-01	XYP Shipping Box	\$ 242.81
CN-0147-01	HTS Shipping Box	\$ 220.50
CN-0148-01	SDS Shipping Box	\$ 114.19
CN-0151-01	O Ring, Quick Disconnect Coupling Replacement Seal (Qty 14)	\$ 74.81
CN-0152-01	CAN Bus Cable	\$ 87.94

CN-0154-01	Sample Line Shipping Plug	\$ 45.94
CN-0158-01	Total System Installation/Evaluation Form, Quantity of 50	\$ 64.31
CN-0162-01	LED Housing with Quick Disconnect Fitting	\$ 84.00
CN-0163-01	Probe Holder Assembly (with Thumbscrews)	\$ 112.88
CN-0172-01	2 Amp Fuse, 250 Volt, SloBLo, (Qty 2)	\$ 23.63
CN-0183-01	Installation Service, Luminex 100/200 and Luminex 100/200 IS (Europe)	\$ 1,312.50
CN-0184-01	Installation Service, Luminex 100/200 and Luminex 100/200IS (North America)	\$ 1,312.50
CN-0199-01	Washable Teflon Coated Tips	\$ 1,253.44
CN-0208-01	De-installation Service for LX100/200 systems (North America)	\$ 1,312.50
CN-0209-01	De-installation Service for LX100/200 systems (EU)	\$ 1,312.50
CN-0215-01	6 Month PM Kit (FM3D, Increased Throughput)	\$ 69.56
CN-0221-01	Sample Probe Needle	\$ 102.38
CN-0222-01	Air Filter: Access Door	\$ 38.06
CN-0223-01	Heater Block for 384-Well Microtiter Plates	\$ 140.34
CN-0224-01	Heater Block for 96-Well Microtiter Plates	\$ 96.90
CN-0225-01	Off-Plate Reservoir	\$ 116.81
CN-0226-01	Fuse (6 Amp)	\$ 28.88
CN-0239-01	Kit, 2 Port PCI Serial Card in Dell GX760 PC	\$ 132.56
CN-0257-01	Air Filter, 4.5 x 4.5	\$ 43.31
CN-0258-01	Drive Fluid Filter	\$ 105.00
CN-0259-01	Sample to Valve Tubing Assembly	\$ 70.88
CN-0263-01	Sample Probe Height Adjustment Kit	\$ 35.44
CN-0267-01	Assy, Air Tube, 36"	\$ 53.81
CN-0268-01	Assy, Sheath Out Tube, 36"	\$ 53.81
CN-0269-01	Magnetic Plate Separator	\$ 680.00
CN-P001-01	Power Cord (USA)	\$ 89.25
CN-P002-01	Power Cord (Australia)	\$ 84.00
CN-P003-01	Power Cord (NEMA 5-15P to IEC-60320-C13)	\$ 84.00
CN-P004-01	Power Cord (Denmark)	\$ 84.00
CN-P005-01	Power Cord (Germany, Sweden, France, Belgium, and Spain)	\$ 84.00
CN-P006-01	Power Cord (Switzerland)	\$ 84.00
CN-P007-01	Power Cord (Israel)	\$ 103.69
CN-P008-01	Power Cord (UK)	\$ 95.81
CN-P009-01	Power Cord (Italy)	\$ 84.00
CN-P010-01	Power Cord (Japan)	\$ 84.00
CN-P012-01	Power Cord (China)	\$ 87.94
CN-P013-01	Power Cord (India, South Africa)	\$ 78.75