

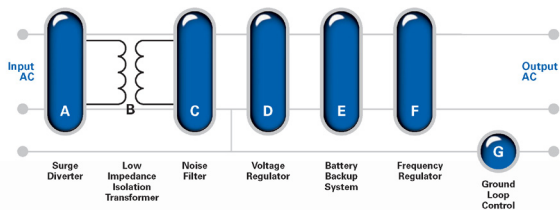


Power Conditioners and UPS Systems

Avoid all of the unnecessary costs – financial and otherwise – of equipment downtime.

Whatever your business, equipment that crashes can cause real problems for your entire operation.

Damaged or compromised components, disrupted processes, lack of reliability – they all add up to frustration, broken schedules and costly downtime. The problem is caused by a host of power disturbances – some visible and many invisible – that threaten your equipment’s operation every year. Power “disturbances” are simply those large and small variations in the quality of the electric power you use day in and day out. Some come from your local utility company, but most are created within your own facility as a result of the distribution and use of electrical power. These variations include high-energy voltage transients, sags and swells, electrical noise, and common mode voltage, as well as the power outages everyone witnesses when the lights go out.



The ABC's of power conditioning.

FEATURES & BENEFITS

- Conditioners provide clean power at 120, 208, or 240 VAC for loads up to 62 amps
- Dual-output models can protect both high-voltage system loads and low-voltage peripherals at less expense and effort than installing two conditioners
- Protects against: high-voltage surges, common-mode voltage, electrical noise, voltage swells and sags, power outages, unstable AC frequency, ground loops
- All models include POWERVAR®'s five-year warranty on parts and labor



Power Conditioners

Instrument Model	Description	Frequency (Hz)	Part No.
ATOMIC ABSORPTION			
AAAnalyst™ 100/200/300/400	520 VA Power Conditioner	60	N9307504
AAAnalyst 100/200/300/400/PinAAcle™ 900 F	750 VA Power Conditioner	50	N9307521
AAAnalyst 600/800	5.8 kVA Power Conditioner	60	N9307511
AAAnalyst 600/800	6.0 kVA Power Conditioner	50	N9307523
AAAnalyst 700	3.8 kVA Power Conditioner	60	N9307509
AAAnalyst 700	3.6 kVA Power Conditioner	50	N9307522
PinAAcle 900 H/T/Z	5.8 kVA Power Conditioner	60	N9307760
PinAAcle 900 H/T/Z	6.0 kVA Power Conditioner	50	N9307523
PinAAcle 900 F	720 VA Power Conditioner	60	N9307515
ICP-OES			
Optima™ 2x00/4x00/5x00/7x00/8x00	3.8 kVA Power Conditioner	60	N9307512
Optima 2x00/4x00/5x00/7x00/8x00	3.6 kVA Power Conditioner	50	N9307522
ICP-MS			
ELAN® 6x00/9000 Controller Side	3.8 kVA Power Conditioner	60	N9307519
ELAN 6x00/9000 Controller Side	3.6 kVA Power Conditioner	50	N9307522
AAAnalyst 600/800/ELAN 6x00/9000 RF Generator Side	5.8 kVA Power Conditioner	60	N9307511
ELAN 6x00/9000/NexION® RF Generator Side	6.0 kVA Power Conditioner	50	N9307523
NexION	5.0 kVA Power Conditioner	60	N0777690
TITAN MPS™ SYSTEMS			
Titan MPS	2.0 kVA Power Conditioner	60	N9306758
Titan MPS	2.0 kVA Power Conditioner	50	N9306755
GAS CHROMATOGRAPHY			
Clarus® 580/680 GC	1920 VA Power Conditioner	60	N9307500
Clarus 580/680 GC	2.0 kVA Power Conditioner	50	N9306755
Clarus SQ 8 (MS Portion)	840 VA Power Conditioner	60	N9307517
Clarus/SQ 8 Complete System	3.8 kVA Power Conditioner	60	N9307510
Clarus/SQ 8 Complete System	3600 VA Power Conditioner	50	N9307522
LIQUID CHROMATOGRAPHY			
Flexar™, Series 200, Flexar FX-10, Flexar FX-15	1440 VA Power Conditioner	60	N9307502
Flexar, Series 200, Flexar FX-10, Flexar FX-15	2.0 kVA Power Conditioner	50	N9306752
Flexar SQ 300/AxION® TOF	2.0 kVA Power Conditioner	60	N9306753
Flexar SQ 300 MS/AxION TOF	2.0 kVA Power Conditioner	50	N9306755
THERMAL AND ELEMENTAL			
DSC 7/TGA 7	240 VA Power Conditioner	60	N9307507
TGA 7	1920 VA Power Conditioner	60	N9307501
PYRIS™ 1 DSC, Elemental 2400	480 VA Power Conditioner	60	N9307513
DSC 6/Diamond DMA	360 VA Power Conditioner	60	N9307514
TMA 7	720 VA Power Conditioner	60	N9307515
2400 Series II®	1920 VA Power Conditioner	60	N9307503
UV/VISIBLE			
LAMBDA™	520 VA Power Conditioner	60	N9307505
LAMBDA	480 VA Power Conditioner	60	N9307513

On-Line Conditioned Uninterruptable Power Supply Systems

With the Security Plus Series, you get much more protection and a higher comfort level than you get with most other UPS systems. The Security Plus Series also provides complete power conditioning and, because the Security Plus Series features on-line inverter design, added peace of mind. And regardless of input fluctuations, the Security Plus Series ensures that the output remains continuous and regulated.

UPS Systems



Instrument Model	Description	Frequency (Hz)	Part No.
ATOMIC ABSORPTION			
PinAAcle 900F/AAnalyst 100/200/300/400	800 VA UPS	60	N0777681
PinAAcle 900F/AAnalyst 100/200/300/400	800 VA UPS	50	N0777689
AAnalyst 700/HGA 900	5.2 kVA UPS	50/60	N9308150
PinAAcle 900 T/H/Z	10 kVA UPS	50/60	N9306757
AAnalyst 600/800	10 kVA UPS	50/60	N0777613
ICP-OES			
Optima 2x00/4x00/5x00/7x00/8x00	5.2 kVA UPS	50/60	N0777511
ICP-MS			
NexION	8.0 kVA UPS	50/60	N0777682
ELAN	12 kVA UPS	50/60	N0777719
GAS CHROMATOGRAPHY			
Clarus 580/680 GC	3000 VA Power Conditioned UPS	60	N9302556
Clarus 580/680 GC/SQ 8 MS	5.2 kVA On-Line Conditioned UPS	50/60	N9308150
LIQUID CHROMATOGRAPHY			
Flexar, Series 200, Flexar FX-10, Flexar FX-15	1440 VA Power Conditioned UPS	60	N9306750
Flexar SQ 300 MS/AxION TOF	3.0 kVA Power Conditioned UPS	50/60	N9306751
AxION, Flexar SQ 300 MS	5.2 kVA On-Line Conditioned UPS	50/60	N0777511
UV/VISIBLE			
LAMBDA	800 VA Power Conditioned UPS	60	N0777681
LAMBDA	800 VA Power Conditioned UPS	50	N0777689
Spectrum® 65/100/100N FT-IR	420 VA Power Conditioned UPS	60	N9306756
ACCESSORIES			
Power Probe	0-250 V input	50/60	N3151391
Opto-Isolator			N9306754

Every day, you can count on PerkinElmer to provide you with solutions that deliver reliable performance, control operating costs and maximize operational time. Our complete portfolio of consumables, parts, supplies, training and service helps you meet both routine and demanding measurement challenges. We invest heavily in testing and validating our products to ensure you receive guaranteed compatibility and performance – on-time, for every instrument in your laboratory.

For a complete listing of consumables, please visit www.perkinelmer.com/supplies

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2011-2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

009595C_01