

PURE GAS
PURE POWER
PURE WATER

Get 15% off your essential accessories for a limited time only.

Power Conditioners

Features and 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



Product specifications.

Instrument Model	Description	Input Voltage	Frequency (Hz)	Part No.
Atomic Absorption				
AAAnalyst 100/200/300/400	520 VA	120 V	60	N9307504
PinAAcle 500/900 F	840 VA	120 V	60	N9307517
AAAnalyst 100/200/300/400, PinAAcle 500/900 F	750 VA	220/230 V	50	N9307521
AAAnalyst 600/800	5.8 kVA	208/240 V	60	N9307511
AAAnalyst 600/800	6.0 kVA	220/230 V	50	N9307523
AAAnalyst 700	3.8 kVA	208/240 V	60	N9307509
AAAnalyst 700	3.6 kVA	220/230 V	50	N9307522
PinAAcle 900 H/T/Z	5.8 kVA	208/240 V	60	N9307760
PinAAcle 900 H/T/Z	6.0 kVA	220/230 V	50	N9307523

Product specifications continued.

Instrument Model	Description	Input Voltage	Frequency (Hz)	Part No.
Atomic Absorption				
PinAAcle 500/900 F	1440 VA	120 V	60	N9307503
PinAAcle 500/900 F	1500 VA	220/230 V	50	N9306752
Inductively Coupled Plasma-OES				
Optima™ 2x00/4x00/5x00/7x00/8x00, Avio 200/500	3.8 kVA	208/240 V	60	N9307512
Optima 2x00/4x00/5x00/7x00/8x00, Avio 200/500	3.6 kVA	220/230 V	50	N9307522
Inductively Coupled Plasma-MS				
ELAN® 6x00/9000 Controller Side	3.8 kVA	208/240 V	60	N9307519
ELAN 6x00/9000 Controller Side	3.6 kVA	220/230 V	50	N9307522
ELAN 6x00/9000 RF Generator Side	5.8 kVA	208/240 V	60	N9307511
ELAN 6x00/9000, NexION® RF Generator Side	6.0 kVA	220/230 V	50	N9307523
NexION 300/350/1000/2000	5.0 kVA	208/240 V	60	N077690
Titan MPS™				
Titan MPS	2.0 kVA	208/240 V	60	N9306758
Titan MPS	2.0 kVA	220/230 V	50	N9306755

Laboratory Water Purification Systems

Intuitive and innovative system for all your pure water needs.

The PURELAB flex is designed to deliver accuracy, flexibility and ease of use.

Provides perfect water purity for analytical and life science applications which require Type III water up to ultrapure Type I (18.2 MΩ-cm) water.

simple installation and quick purification pack replacement

- UV lamp to reduce total organic carbon (TOC)
- Realtime TOC monitor for critical analytical applications
- Different dispensing options



15% off all Gas generators for GC, GC/MS and LC/MS

FID Gas Station

The FID gas station is a novel system that combines the reliability of the high purity PG or ultrahigh purity NM hydrogen generators with the superior performance of the zero air generators into a single compact instrument. Designed with space saving in mind, it sits directly under the GC to conserve bench space. You can also apply the FID gas station to GCMS systems with the simple addition of a drawer tray for the MS to rest on. If you want to benefit from the space saving design, but only require a H₂ generator that is also an option.

Features and Benefits

- H₂ flow rate from 100-1000 mL/min
- H₂ pressure up to 11 bar
- Zero air flow rates 5 L/min
- Zero air purity <0.1 ppm
- Minimal maintenance
- CE, CSA, FCC certified



And more gas generators to choose from...

Product specifications.

Description	Capability	Flow Rate	Part No.
FID Station PG Plus	H ₂	150 mL/min	N9300023
FID Station PG Plus	H ₂	260 mL/min	N9300024
FID Station NM Plus	H ₂	150 mL/min	N9300020
FID Station NM Plus	H ₂	350 mL/min	N9300021
FID Station NM Plus	H ₂	650 mL/min	N9300022
FID Station PG Plus	H ₂ and Air	150 mL/min + Air 1500 mL/min	N9300018
FID Station PG Plus	H ₂ and Air	260 mL/min + Air 1500 mL/min	N9300019
FID Station NM Plus	H ₂ and Air	150 mL/min + Air 1500 mL/min	N9300015
FID Station NM Plus	H ₂ and Air	350 mL/min + Air 1500 mL/min	N9300016
FID Station NM Plus	H ₂ and Air	650 mL/min + Air 1500 mL/min	N9300017
FID station NM Plus extension drawer			N9308567

ONE-TIME OFFER

**15% OFF discount
on parts listed here.**

To take advantage of this offer, visit our website and use **promotion code: AMSQ419** at checkout.

* Terms and Conditions: 15% discount on Gas Generators, Water Purification systems, and Line Conditioners. Not cumulative with others discounts, promotions & deals. Valid one time per customer. Not cumulable with other promotions or deals or discount. Limit one offer per customer. Offer valid while supplies last. Other restrictions may apply. Void where prohibited. Valid in US, Canada, Central America, Western Europe, South America, India, Greater China, Japan, Korea. Valid until 12/31/19.

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 ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.