

# **GENERATORS**

PerkinElmer offers a comprehensive range of analytical gas supply solutions that goes beyond simply 'gas generators'. PerkinElmer is committed to providing analytical chemists with gas solutions that enable them to get the most out of their instruments. Whether the requirement is for ultra high purity hydrogen for use as a GC and GC/MS carrier gas or multiple gases to operate sophisticated LC/MS instrumentation, PerkinElmer's offerings always keep the end user in mind.

#### **HYDROGEN GENERATORS**



#### **Features and Benefits**

- Ultra safe operation
- >99.999% purity
- High purity guaranteed, 24/7
- Class leading PEM cell featuring 2-year warranty as standard
- CE, UL & CSA approved
- Low life cycle cost, excellent return on investment

# ULTRA HIGH PURITY HYDROGEN GENERATORS



### **Features and Benefits**

- Ultra safe operation
- >99.9999% purity
- High purity guaranteed, 24/7
- Class leading PEM cell featuring 2-year warranty as standard
- CE, UL & CSA approved
- Low life cycle cost, excellent return on investment

### **ZERO AIR GENERATORS**



## **Features and Benefits**

- High performance heated platinum catalyst
- High efficiency filtration
- Modular and stackable with all PerkinElmer hydrogen generators
- CE, UL & CSA approved
- Low life cycle cost, excellent return on investment
- Superior resolution and lowered limits of detection
- Safe and sustainable

# LC/MS NITROGEN GENERATORS



## **Features and Benefits**

- Continuous automatic operation, 24/7
- Phthalate-free componentry
- Low life cycle cost, excellent return on investment
- CE, UL & CSA approved
- Safe and sustainable
- Field proven Pressure Swing Adsorption (PSA) technology
- Compact and robust design



## **PRODUCT SELECTION**

# Hydrogen Gas Generators for Fuel Gas Applications

Model	Part No.	Flow Rate (mL/min)	Purity (%)	Water Consumption (24/7, full flow) (L/week)	Delivery Pressure (psi/g)	Height (H) (in.)	Width (W) (in.)	Depth (D) (in.)	Weight (Empty) (lbs.)
20H	N9303225	160	>99.999	1.25	5-100	17.9	13.5	17.2	41.9
40H	N9303226	250	>99.999	2	5-100	17.9	13.5	17.2	41.9
60H	N9303227	500	>99.999	4	5-100	17.9	13.5	17.2	41.9

## Ultra High Purity Hydrogen Gas Generators for Carrier Gas Applications

Model	Part No.	Flow Rate (mL/min)	Purity (%)	Water Consumption (24/7, full flow) (L/week)	Delivery Pressure (psi/g)	Height (H) (in.)	Width (W) (in.)	Depth (D) (in.)	Weight (Empty) (lbs.)
20H-MD	N9303201	160	>99.999	1.69	10-100	17.9	13.5	18.5	45.2
40H-MD	N9303202	250	>99.999	2.41	10-100	17.9	13.5	18.5	45.2
60H-MD	N9303203	500	>99.999	4.82	10-100	17.9	13.5	18.5	45.2
110H-MD	N9303204	1100	>99.999	10.60	10-100	17.9	13.5	18.5	51.8

Ambient Temperature Range		41-104 °F			
Water Quality		Deionized, ASTM® II, >1M $\Omega$ , <1 $\mu$ s, filtered to <100 $\mu$ m			
Maximum Water Supply Pressure/Flow*		1.45 psi g / 1 L/min			
Supply Voltage Range		Universal 120/230 V ±10% (60/50 Hz)			
Port Connections	Hydrogen Outlet Water Drain	1/8" Compression Fitting Quick Release Push in Fitting			

<sup>\*</sup>With Auto Water Refill option

### Zero Air Generators

Model	Part No. (120 V)	Part No. (230 V)	Flow Rate (L/min)	Organic Purity (ppm)	Air Inlet @ 58-145 psig (L/m)	Delivery Pressure (psi/g)	Height (H) (in.)	Width (W) (in.)	Depth (D) (in.)	Weight (lbs.)
UHP-35ZA-S	N9303206	N9303205	3.5	<0.1	4.2	58-145	17.9	13.4	16.7	31.3
UHP-50ZA-S	N9303208	N9303207	5.0	<0.1	6.0	58-145	17.9	13.4	16.7	31.3
UHP-150ZA-S	N9303210	N9303209	15	<0.1	18	58-145	17.9	13.4	16.7	31.3
UHP-300ZA-S	N9303212	N9303211	30	<0.1	35	58-145	17.9	13.4	16.7	31.3

Ambient Temperature Rar	nge	41-104 °F			
Required Inlet Air Quality		Clean dry compressed air ISO8573-1:2001 Class 3.2.1			
Supply Voltage Range		120 V/60 Hz or 230 V/50-60 Hz ±10%			
Port Connections	Outlet (N9303205 and N9303206) Inlet (N9303205 and N9303206) Outlet (N9303207-N9303212 Inlet (N9303207-N9303212)	1/8" Compression Fitting 1/8" Compression Fitting 1/4" Compression Fitting 1/4" Compression Fitting			

## LC/MS Nitrogen Generators

Model	Part No. (120 V)	Part No. (230 V)	Flow Rate (L/min)	Purity (ppm)	Delivery Pressure (psi/g)	Height (H) (in.)	Width (W) (in.)	Depth (D) (in.)	Weight (lbs.)	
With integra	ted compresso	or								
LCMS30-1	N9303223	N930224	30	>98	101.5	27.8	20.1	22.9	298	
Ambient Ter	nperature Ran	ge			41-104 °F					
Supply Volta	ge Range				120 V/60 Hz or 230 V/50-60 Hz ±10%					
Port Connec	tions Nitrogen	Outlet			1/4" Compression Fitting					

<sup>\*</sup>Visit www.perkinelmer.com for a complete listing of flow rates.

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

