

Hydrogen and Nitrogen Gas Generators



Simply a Smarter Choice

Analytical gas generators can remove the requirement for high-pressure cylinder gases and are typically placed next to the instrument they are servicing. This removes any need for extended gas lines and negates any associated problems impacting on purity, cost and convenience.

The latest gas generators from PerkinElmer utilize new technologies in adsorbents, catalysts, and specialist micro dryers to produce a continual supply of ultra-high purity gases to your instrumentation. This all but eliminates the introduction of impurities, which can be reduced further by the installation of in-line gas purifiers.



FID Tower Plus Hydrogen Gas Generator

	NM PLUS					PG PLUS
Part Number	N9300011	N9300012	N9300013	N9300014	N9300026	N9300002
H2 Flow (mL/min)	250	300	500	600	1000	250
Hydrogen Purity	99.99996%					99.99996%
Application	Carrier and Detector gas					Detector Gas
Dew point at 7 Barg (100 psig)	-25 °C (-77 °F)					-25 °C (-77 °F)
Outlet Pressure Barg (psig)	1.4 to 11 (20 to 160)					0.5 to 11 (7 to 160)
Technology	PEM (Proton Exchange Membrane) - 100% Titanium cell					PEM (Proton Exchange Membrane) - 100% Titanium cell
Drying System	No maintenance cold dual dynamic regeneration system					No maintenance cold dual dynamic regeneration system
Deionized Water Quality	Minimum < 1 micro S/cm @ 25 °C - 1 MΩ-cm @ 25 °C Recommended < 0.2 micro S/cm @ 25 °C - 5 MΩ-cm @ 25 °C					Minimum < 1 micro S/cm @ 25 °C - 1 MΩ-cm @ 25 °C Recommended < 0.2 micro S/cm @ 25 °C - 5 MΩ-cm @ 25 °C
Internal Water Tank (L)	5					5
Safety	Automatic shut down - internal/external hydrogen leak, overpressure and low water					Automatic shut down - internal/external hydrogen leak, overpressure and low water
Display	Touch screen with operating parameters, system status and safety alarms					Touch screen with operating parameters, system status and safety alarms
Electrical Supply	110-120 V 60 Hz / 220-240 V 50 Hz					110-120V 60Hz / 220-240V 50Hz
Power Consumption (W)	140	150	200	300	400	115
Cascading Capable	Yes					Yes



FID Tower Plus Hydrogen Gas Generator with Zero Air

Part Number	NM PLUS ZERO AIR					PG PLUS ZERO AIR
	N9300003	N9300004	N9300005	N9300006	N9300007	N9300009
H2 Flow (mL/min)	250	300	500	600	1000	250
Zero Air Flow (mL/min)	5000					5000
Purity - Hydrocarbons + CO	<0.1 ppm					<0.1 ppm
Inlet Pressure Barg (psig)	4.5 to 10 (65 to 145)					4.5 to 10 (65 to 145)
Inlet Air Quality	Clean dry compressed air ISO8573-1:2010 Class 1.2.1					Clean dry compressed air ISO8573-1:2010 Class 1.2.1
Max Outlet Pressure Barg (psig)	5 (73)					5 (73)
Max HC in	100 ppm					100 ppm
Max CO in	50 ppm					50 ppm



FID Tower Plus Nitrogen Gas Generator

N2 TOWER WITH OVEN	
Part Number	N9300008
N2 Flow (mL/min)	750
Nitrogen Purity	99.999%
Hydrocarbon Purity (Measured as Methane)	N/A
Dewpoint	-50 °C
Outlet Pressure Barg (psig)	Up to 5 max (75)
Inlet Pressure Barg (psig)	7 to 10 (100 to 160)
Actual Inlet Air Requirement (L at 8 Barg)	12
Recommended Compressor Air Inlet (at 8 Barg)	24
Pressure Drop Barg (psig)	1.5 (22)
Inlet Air Quality	Clean dry compressed air ISO8573-1:2010 Class 1.2.1
Technology	Carbon molecular sieve
Warm Up Time (min)	60
Electrical Supply	110-120 V 60 Hz / 220-240 V 50 Hz
Consumption (W)	300



FID Station Plus Hydrogen Gas Generator

	NM PLUS			PG PLUS	
Part Number	N9300020	N9300021	N9300022	N9300023	N9300024
H2 Flow (mL/min)	150	350	650	150	260
Application	Carrier and Detector gas			Detector gas	
Hydrogen Purity	99.99996%			99.9996%	
Dew point at 7 Barg (100 psig)	-73 C° (-103 °F)			-25 C° (-77 °F)	
Outlet Pressure Barg (psig)	1.4 to 11 (20 to 160)			0.5 to 11 (7 to 160)	
Technology	PEM (Proton Exchange Membrane) - 100% Titanium cell			PEM (Proton Exchange Membrane) - 100% Titanium cell	
Drying System	Dual Dynamic Regeneration			Regenerative Permeation Membrane	
Deionized Water Quality	Minimum < 1 micro S/cm @ 25 °C - 1 MΩ-cm @ 25 °C Recommended < 0.2 micro S/cm @ 25 °C - 5 MΩ-cm @ 25 °C			Minimum < 1 micro S/cm @ 25 °C - 1 MΩ-cm @ 25 °C Recommended < 0.2 micro S/cm @ 25 °C - 5 MΩ-cm @ 25 °C	
Internal Water Tank (L)	7			7	
Safety	Automatic shut down - internal/external hydrogen leak, overpressure and low water			Automatic shut down - internal/external hydrogen leak, overpressure and low water	
Display	Touch screen with operating parameters, system status and safety alarms			Touch screen with operating parameters, system status and safety alarms	
Electrical Supply	110-120 V 60 Hz / 220-240 V 50 Hz			110-120 V 60 Hz / 220-240 V 50 Hz	
Power Consumption (W)	100	200	350	100	120
Cascading Capable	Yes			Yes	

	NM PLUS ZERO AIR			PG PLUS ZERO AIR	
Part Number	N9300015	N9300016	N9300017	N9300018	N9300019
H2 Flow (mL/min)	150	350	650	150	260
Zero Air Flow (mL/min)	1500			1500	
Purity - Hydrocarbons + CO	<0.1 ppm			<0.1 ppm	
Inlet Pressure Barg (psig)	4.5 to 10 (65 to 145)			4.5 to 10 (65 to 145)	
Inlet Air Quality	Clean dry compressed air ISO8573-1:2010 Class 1.2.1			Clean dry compressed air ISO8573-1:2010 Class 1.2.1	
Max Outlet Pressure Barg (psig)	5 (73)			5 (73)	
Max HC in	100 ppm			100 ppm	
Max CO in	50 ppm			50 ppm	



Hydrogen Gas Generator

NM PLUS

Part Number	N9308581	N9308582	N9308583	N9308584	N9308585
H2 Flow (mL/min)	100	160	250	500	1000
Hydrogen Purity	99.99996%				
Application	Carrier and Detector gas				
Dew Point at 7 Barg (100 psig)	-25 °C (-77 °F)				
Outlet Pressure Barg (psig)	1.4 to 11 (20 to 160)				
Technology	PEM (Proton Exchange Membrane) - 100% Titanium cell				
Drying System	No maintenance cold dual dynamic regeneration system				
Deionized Water Quality	Minimum < 1 micro S/cm @ 25 °C - 1 MΩ-cm @ 25 °C Recommended < 0.2 micro S/cm @ 25 °C - 5 MΩ-cm @ 25 °C				
Internal Water Tank (L)	2.5				
Safety	Automatic shut down - internal/external hydrogen leak, overpressure and low water				
Display	Touch screen with operating parameters, system status and safety alarms				
Electrical Supply	110-120 V 60 Hz / 220-240 V 50 Hz				
Power Consumption (W)	100	120	165	220	385
Cascading Capable	Yes				

PG PLUS

Part Number	N9308577	N9308578	N9308579	N9308580
H2 Flow (mL/min)	100	160	250	500
Hydrogen Purity	99.9996%			
Application	Detector gas			
Dew Point at 7 Barg (100 psig)	-25 °C (-77 °F)			
Outlet Pressure Barg (psig)	0.5 to 11 (7 to 160)			
Technology	PEM (Proton Exchange Membrane) - 100% Titanium cell			
Drying System	Regenerative Permiation Membrane			
Deionized Water Quality	Minimum < 1 micro S/cm @ 25 °C - 1 Mohm-cm @ 25 °C Recommended < 0.2 micro S/cm @ 25 °C - 5 Mohm-cm @ 25 °C			
Internal Water Tank (L)	2.5			
Safety	Automatic shut down - internal/external hydrogen leak, overpressure and low water			
Display	Touch screen with operating parameters, system status and safety alarms			
Electrical Supply	110-120 V 60 Hz / 220-240 V 50 Hz			
Power Consumption (W)	75	95	140	190
Cascading Capable	Yes			

Shop Online Today!
<https://www.perkinelmer.com/buy>

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 ©2021 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.