

Clarus Gas Chromatograph









The Clarus GC Family – Performance you can rely on from the leader in Gas Chromatography and sample-handling solutions




With more than 50 years' experience in Gas Chromatography (GC), the PerkinElmer® family of Clarus® GC instruments can be depended upon to meet the most demanding performance and throughput needs for today's analytical, process monitoring and quality control facilities.

Glass Inlet Liners for Injector Systems

Programmed Temperature Split/Splitless (PSS) Injector Liners

Description	Dimensions (mm) i.d. x o.d. x length	Pkg.	Part No.
 Quartz Split Liner PerkinElmer Standard (ships with instrument)	2 x 4 x 86.2	1	N6121004
 Quartz Split Liner with Silanized Glass Wool	2 x 4 x 86.2	1	N6121008
 Quartz Split Liner with Silanized Glass Wool	2 x 4 x 86.2	5	N6121009
 Quartz Splitless Liner Excellent for low-volume analysis (ships with instrument)	1 x 4 x 86.2	1	N6121006
 On-Column Glass Liner	2.4 x 4 x 86.2	1	N6101539
 Borosilicate Glass Split Liner This liner is loosely packed with wool for the specific use of analyzing fuel in oil	2 x 4 x 86.2	5	N9302949
Silanized Glass Wool (2 oz.)		1	03300905



Capillary Split/Splitless Injector Liners

Description	Dimensions (mm) i.d. x o.d. x length	Pkg.	Part No.
 Quartz Liner for Split Operation Good for large volume injections	4 x 6.2 x 92.1	1	N6121001
 Quartz Liner for Split Operations with Silanized Glass Wool Good for large volume injections	4 x 6.2 x 92.1	1	N6121010
 Quartz Liner for Split Operations with Silanized Glass Wool Good for large volume injections	4 x 6.2 x 92.1	5	N6121011

Capillary Split/Splitless Injector Liners, *continued*

	Description	Dimensions (mm) i.d. x o.d. x length	Pkg.	Part No.
	Quartz Liner for Splitless Operation Standard Injector Liner (ships with instrument)	2 x 6.2 x 92.1	1	N6121002
	Glass Liner for Split Operation Universal liner for general purpose analysis	4 x 6.2 x 92.1	1	N6101052
	Glass Liner for Splitless Operation Universal liner for general purpose analysis	2 x 6.2 x 92.1	1	N6101372
	Precision Split Glass Liner Deactivated Surface with Wool Wool can be adsorptive if fibers are broken	4 x 6.2 x 92.1	5	N6121020
	Zero Dilution Glass Outer Liner Ideal for trace headspace work	2.7 x 6.3 x 69	1	N1011445
	Zero Dilution Inner Liner Use with Outer Liner (N1011445)	1 x 2 x 73	1	N1011446

Packed Column Injector Liners

	Description	Dimensions (mm) i.d. x o.d. x length	Pkg.	Part No.
	Wide-Bore Column On/Off Quartz Liner	6 x 4 x 92.1	1	N6121003
	Wide-Bore Column Glass Liner	6 x 4 x 92.1	1	N6101375
	Drilled Uniliner (hole on bottom) Good for PPC equipped GCs. Recommended for analysis in which compounds of interest could be affected by a tailing solvent peak	4 x 6.2 x 92.1	5	N6502013
	Gooseneck Drilled Uniliner (hole on top) Use for trace, active samples. Gives high recovery and linearity	4 x 6.2 x 92.1	5	N6502014

O-rings

Description	Pkg.	Part No.
For Injector Split/Splitless		
O-ring Silicone – Max. Temp. 250 °C	10	N6101374
O-ring Kalrez – Max. Temp. 450 °C	1	N9302782
For Injector PSS		
O-ring Silicone – Max. Temp. 250 °C	10	N6101747
O-ring Kalrez – Max. Temp. 450 °C	1	09921004

Capillary Column Ferrules

Description	Pkg.	Part No.
Ferrules for Injector		
Graphite Ferrule 1/16 in, for Capillary Column from 0.18 mm to 0.32 mm	10	09903700
Graphite Ferrule 1/8 in, for Capillary Column from 0.32 mm to 0.53 mm	10	09920141
Ferrules for Detector		
Graphite Ferrule 1/8 in, for Capillary Column from 0.18 mm to 0.32 mm	10	09903981
Graphite Ferrule 1/8 in, for Capillary Column from 0.32 mm to 0.53 mm	10	09903394
Ferrules for Split/Splitless and PSS		
Column Nut 1/16 in for Reversed Ferrule	5	09903392

Autosampler Syringes



Syringes from PerkinElmer are individually inspected for accuracy and performance. Recommended auto-sampler syringes are available in 0.5, 5 and 50 μL capacities. For routine analyses, the metal plunger in barrel with PTFE-tipped seal is the standard syringe as shipped with each Clarus GC instrument. Alternative syringes to use are the metal plunger in barrel or the 0.53 mm on-column injection.

Description	Pkg.	Part No.
50 μL Syringe, Metal Plunger 0.63 mm o.d. Needle	1	N6101760
5 μL Syringe, Metal Plunger PTFE-tipped Seal 0.63 mm o.d. Needle	1	N6101390
5 μL Syringe, Metal Plunger PTFE-tipped Seal 0.63 mm o.d. Needle BLUE Barrel	5	N6103240
5 μL Syringe, Metal Plunger PTFE-tipped Seal 0.63 mm o.d. Needle BLUE Barrel	10	N6103241
5 μL Syringe, Metal Plunger 0.63 mm o.d. Needle	1	N6101251
0.5 μL Low Injection Volume Syringe, Metal Plunger 0.63 mm o.d. Needle	1	N6101252
0.5 μL Low Injection Volume Syringe Metal Plunger 0.63 mm o.d. Needle BLUE Barrel	5	N6103242
0.5 μL Low Injection Volume Syringe Metal Plunger 0.63 mm o.d. Needle BLUE Barrel	10	N6103243

On-Column Injection

5 μL On-column Syringe Metal Plunger 0.47 mm o.d. Needle	1	N6101380
0.5 μL Low Injection Volume Syringe, Metal Plunger 0.47 mm o.d. Needle	1	N6101253



Injector Septa

Description	Pkg.	Part No.
PerkinElmer Green Injection Septum, 11 mm diameter These septum have extremely low bleed over a wide range of inlet temperatures – 100 to 350 $^{\circ}\text{C}$. Easier needle penetration and high puncture tolerance make this septum ideal for autosamplers. This septum is already conditioned and ready to use.	50	N6621028
BTO™ (Bleed Temperature Optimized) Injector Septa Maximum recommended operating temperature 300 $^{\circ}\text{C}$. 11 mm diameter.	25	N9303343
PTFE/Silicone Injector Septa, 11 mm diameter	50	00090652
Green Injection Port Septa, 11 mm diameter Rated to 400 $^{\circ}\text{C}$. The advanced green septum combines significantly longer injection life, low bleed and low injection port adhesion. The result is a general use green septum made of uniquely formulated silicone rubber that can be used for all your daily analyses. Packaged in a pre-cleaned glass screw top jar for high purity.	10 50	N9306218 N9306219
Orange Injection Port Septa Rated to 400 $^{\circ}\text{C}$. Uniquely formulated silicone rubber septa is Bleed and Temperature Optimized (BTO®) for today's most demanding GC and GC/MS applications. Septa BTO® is formulated to extend low-bleed and outstanding mechanical properties of premium GC septa. It retains remarkable softness at high temperatures and has been optimized to reduce injection port adhesion. Packaged in a pre-cleaned glass screw top jar for high purity.	50	N9302972
White Injection Port Septa Rated to 400 $^{\circ}\text{C}$. 125 thick silicone, 9 mm diameter	12	02261144

Injector Kits

Description	Pkg.	Part No.
Split/Splitless Injector Starter Kit – Manual Gas Control	1	N6120101
Kit includes: 2 mm i.d. quartz liner, 4 mm i.d. quartz liner, silicone O-rings Green Injector Port Septa (50), 0.55 mm graphite ferrules (10), 0.8 mm graphite ferrules (10), 1/16 in stainless steel nuts (5), untreated quartz wool, packing rod (1) and water scibes (10)		
PSS Injector Starter Kit	1	N6120102
Kit includes: universal connectors (5), 0.53 mm i.d. deactivated fused silica (5 m), 2 mm i.d. quartz liner, 1 mm i.d. quartz liner, glass hour-glass liner, 1/16 in stainless steel nuts (5), 0.5 mm graphite ferrules (10), 0.8 mm graphite ferrules (10), Viton O-rings (5), graphite O-rings (5), Green Injector Port Septa (50), untreated quartz wool, water scibes (10)		

Vials, Caps and Septa Kits

We have expanded our product selection to cover additional test applications and to make it easier for our customers to order prepackaged kits for their daily production requirements.

Tested to meet your application requirements and guaranteed for fit and compatibility.

Autosampler Crimp and Screw Top Vials, Caps and Septa. Our screw thread vials are custom designed for PerkinElmer Clarus GCs. Manufactured from Type 1 borosilicate glass, the vials have high-temperature tolerance and are chemically inert. Color variety crimp caps offer easy identification within the laboratory on 2 mL crimp top vials. Microsampling with a maximum capacity of 200 µL can be accomplished with the use of a low-volume insert and support foot.

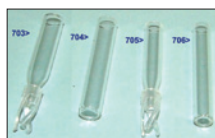
Vials, Caps and Septa Kits

Description	Pkg	Part No.
Crimp Top Vials (also see Convenience Kits for Additional Vials and Caps Sold Separately)		
2 mL 8 mm Crimp Top Clear Glass Vials	200	N9301069
2 mL 11 mm Crimp Top Clear Glass Vials	100	N9301385
2 mL 11 mm Crimp Top Amber Glass Vials	100	N9302680
2 mL 11 mm Crimp Top Clear Wide Mouth Glass Vials with Write-On Patch and Fill Lines	100	N9306223
2 mL 11 mm Crimp Top Amber Wide Mouth Glass Vials with Write-On Patch and Fill Lines	100	N9302679
Crimp Top Septa		
8 mm Silver Crimp Cap with PTFE/Rubber Septa	1000	03300806
11 mm Silver Crimp Caps with PTFE/Silicone Septa	100	N9306228
11 mm Silver Crimp Caps with PTFE/Silicone/PTFE Septa	100	N9306229
11 mm Silver Crimp Caps with Red Rubber Septa	100	N9306230
11 mm Blue Crimp Caps with Teflon®/Rubber Septa	100	N9302686
11 mm Green Crimp Caps with Teflon®/Rubber Septa	100	N9302684
11 mm Red Crimp Caps with Teflon®/Rubber Septa	100	N9302685
11 mm Silver Crimp Caps with Teflon®/Rubber Septa	100	N9306015
11 mm Silver Crimp Cap with Black Viton® Septa	1000	N9302784

Vials, Caps and Septa Kits, *continued*

Description	Pkg.	Part No.
Screw Top Vials/Caps/Septa		
2 mL 8 mm Screw Top Clear Glass Vials	100	N9302945
8 mm Screw Caps – no Septa – Phenolic Cap with Hole	100	N9303441
8 mm Screw Black Caps PTFE/Silicone Septa	100	N9303449
8 mm PTFE-Coated Butyl Rubber Septa For 8 mm Screw Caps	100	N9303442
2 mL 9 mm Screw Top Clear Vials	100	N9306201
2 mL 9 mm Screw Top Amber Vials	100	N9306220
2 mL 9 mm Screw Top Clear Vials with Write-On Patch and Fill Lines	100	N9307801
2 mL 9 mm Screw Top Amber Vials with Write-On Patch and Fill Lines	100	N9307802
9 mm Blue Screw Cap PTFE/Silicone Septa	100	N9306202
9 mm Blue Screw Cap Pre-Slit PTFE/Silicone Septa	100	N9306203
2 mL 10 mm Screw Top Clear Vial	100	N9306053
2 mL 10 mm Screw Top Amber Vial	100	N9306057
10 mm Blue Screw Caps Red Rubber Septa	100	N9306200
10 mm Black Screw Caps PTFE/Silicone Septa	100	N9306205
10 mm Black Screw Caps PTFE/Red Rubber Septa	100	N9306206
Snap Top Vials/Caps/Septa		
2 mL 11 mm Clear Snap Wide Opening Vial	100	N9303418
2 mL 11 mm Clear Snap Wide Opening Vials with Write-On Patch and Fill Lines	100	N9306207
2 mL 11 mm Amber Snap Wide Opening Vials with Write-On Patch and Fill Lines	100	N9306208
11 mm Pre-Slit PTFE/Silicone Clear Snap Cap	100	N9303416
11 mm Pre-Slit PTFE/Silicone/PTFE Clear Snap Cap	100	N9303417
11 mm PTFE/Silicone Clear Snap Cap	100	N9303419
Waste and Wash Vials/Caps		
Cap for Waste and Wash Vials (15 mm)	1 Each	9923032
Clear Waste and Wash Vial 4 mL (15 mm)	1 Each	9923031
Clear Waste and Wash Vial 4 mL (15 mm)	100	N9306247
Septa for Waste Wash Vials (For use with Viscous or Toxic Samples)	50	N9302780
Diffusers for Waste and Wash Vials	1	N6101276
<i>Diffusers are sold exclusively by PerkinElmer. The conical shape allows the needle to penetrate into solvent or waste. Reduces spillage and large losses through evaporation.</i>		
Inserts		
200-µL Low-Volume Clear Glass Insert	1000	N9302681
Vial Support for Insert	500	N9302682

Vials, Caps and Septa Kits, *continued*



Description	Pkg.	Part No.
Low Volume Vial Glass Inserts		
6 mm x 29 mm Clear Glass Insert, Rimless Pulled Point with Bottom Spring, 250 µL Usable Volume	100	N9300703
6 mm x 31 mm Clear Glass Flat Bottom Insert, 400 µL Usable Volume	100	N9300704
Use With:		Vial #
10 mm 2 mL Amber Screw Top Vial with White "P" Icon with Patch		N9306057
10 mm 2 mL Clear Screw Top Vial with White "P" Icon with Patch		N9306053
11 mm Clear Large Opening Snap Ring Vial with Black "P" Icon		N9303418
11 mm Clear Large Opening Crimp Vial with Black "P" Icon		N9306231
9 mm 2 mL Amber Screw Top Vial with Black "P" Icon		N9306220
9 mm 2 mL Clear Screw Top Vial with Black "P" Icon		N9306201
5 mm x 29 mm Clear Glass Insert Rimless Pulled Point with Bottom Spring, 150 µL Usable Volume	100	N9300705
5 mm x 31 mm Clear Glass Flat Bottom Insert, 200 µL Usable Volume	100	N9300706
Use With:		Vial #
11 mm Standard Amber Crimp Vial with Black "P" Icon		N9302680
11 mm Standard Clear Crimp Vial with Black "P" Icon		N9301385
8 mm 2 mL Clear Screw Top Vial with Black "P" Icon		N9302945

Vial Holder Racks

36-Vial Capacity 20 mm Vial Holder Rack	1	N9301304
50-Vial Capacity 11 mm Vial Holder Rack	1	N9301303



NEW! Vials with Fused Glass Sample Inserts

2 mL 11 mm Large Opening Crimp Top Clear 0.3 mL Insert	100	N9300709
2 mL 11 mm Large Opening Crimp Top Amber 0.3 mL Insert	100	N9300710
2 mL 11 mm Large Opening Snap Top Ring Clear 0.3 mL Insert	100	N9300711
2 mL 11 mm Large Opening Snap Top Ring Amber 0.3 mL Insert	100	N9300712
2 mL 8 mm Screw Top Clear 0.1 mL Insert	100	N9300713
2 mL 8 mm Amber Screw Top 0.1 mL Insert	100	N9300714
2 mL 9 mm Large Opening Screw Top Clear 0.3 mL Insert	100	N9300715
2 mL 9 mm Large Opening Screw Top Amber 0.3 mL Insert	100	N9300716
2 mL 10 mm Large Opening Screw Top Clear 0.3 mL Insert	100	N9300717
2 mL 10 mm Large Opening Screw Amber 0.3 mL Insert	100	N9300718

Crimping Tools

Whatever your need may be, PerkinElmer offers a wide range of crimping tools for your convenience. Our universal voltage, precision control, power crimpers with adjustable settings are designed to deliver hundreds of crimps on a single battery charge.



Description	Qty.	Part No.
Ergonomic Hand Crimper		
11 mm Ergonomic Hand Crimper	1	N6621035
11 mm Ergonomic Hand Decapper	1	N6621036
20 mm Ergonomic Hand Crimper	1	N6621037
20 mm Ergonomic Hand Decapper	1	N6621038

Crimping Tools, *continued*



Description	Qty.	Part No.
Hand Held and Bench-top Crimpers		
8 mm Hand Crimper	1	N9306127
11 mm Hand Crimper	1	00090699
20 mm Hand Crimper	1	N9302785
11 mm Crimper Jaws	1	N6621008
20 mm Crimper Jaws	1	N6621009
Bench-top Vial Crimper	1	N6621006
Electronic Crimpers		
20 mm Electronic Universal Crimper with Global Mains Plug Pack	1	N9304501
20 mm Electronic Universal Decapper with Global Mains Plug Pack	1	N9304503
11 mm Electronic Universal Crimper with Global Mains Plug Pack	1	N9304500
11 mm Electronic Universal Decapper with Global Mains Plug Pack	1	N9304502
Rechargeable Battery 6.4 Volt Lithium Ion	1	N9304504

Autosampler Vials/Caps/Septa Kits

PerkinElmer understands your challenges and offers a variety of kits so that you can easily order and restock your laboratory supplies.

Description	Qty.	Part No.
Autosampler Starter Kit	1	N6120105
Crimper Tool (Pkg/1)	1	00090699
Syringe 9000 5.0 µL 0.63 mm o.d. (Pkg/1)	1	N6101390
Vial Rack for 50 Vials-12 mm (Pkg/1)	1	N9301303
2 mL 11-mm Crimp Top Vial, Clear (Pkg/100)	1	N9301385
Decapper 11 mm Tool (Pkg/1)	1	N9301390
2 mL Crimp Top Vial, Amber (Pkg/100)	1	N9302680
11 mm Crimp Cap, Green with Septa	1	N9302684
11 mm Crimp Cap, Red with Septa (Pkg/100)	1	N9302685
11 mm Crimp Cap, Blue with Septa (Pkg/100)	1	N9302686
Septa for Waste/Wash Vials (Pkg/50)	2	N9302780
11 mm Crimp Cap Silver with Red Rubber/PTFE Septa (Pkg/100)	1	N9306015
GC Educational Consumables Kit with Electronic Crimper	1	N6500570
Crimper Tool 11-mm Electronic Hand Crimper with Battery (Pkg/1)	1	N9304500
Syringe 5.0 µL 0.63 mm o.d. (Pkg/1)	2	N6101390
Rack for 50 Vials 12 mm (Pkg/1)	1	N9301303
11-mm Crimp Top Vial, 2 mL (Pkg/100)	2	N9301385
11 mm Crimp Cap Silver with Red Rubber/PTFE Septa (Pkg/100)	2	N9306015
Elite™ Column – 5-30 M x 0.25 µm x 0.25 mm (Pkg/1)	1	N9316076

Autosampler Vials/Caps/Septa Kits, *continued*

Description	Qty.	Part No.
GC Educational Consumables Kit with Manual Crimper	1	N6500571
Hand Crimper Tool (Pkg/1)	1	00090699
Syringe 5.0 µL 0.63 mm o.d. (Pkg/1)	2	N6101390
Rack for 50 Vials, 12 mm (Pkg/1)	2	N9301303
11-mm Crimp Top Vial, 2 mL (Pkg/100)	2	N9301385
11 mm Crimp Cap Silver with Red Rubber/PTFE Septa (Pkg/100)	2	N9306015
Elite Column – 5-30 M x 0.25 µm x 0.25 mm (Pkg/1)	1	N9316076
Screw Top Autosampler Vial Convenience Kit – Pkg/100	1	N9301945
2 mL 8 mm Clear Screw Top Vials (Pkg/100)		
8 mm Black Caps with Preassembled PTFE/Rubber Septa (Pkg/100)		
<i>Packed in an Environmentally Clean, Re-sealable Clamshell Pack</i>		
Crimp Top Autosampler Vial Convenience Kit – Pkg/100	1	N9300654
2 mL 11 mm Clear Crimp Top Vials (Pkg/100)	1	N9301385
11 mm Silver Caps with Preassembled PTFE/Rubber Septa (Pkg/100)	1	N9306015
<i>Packed in an Environmentally Clean, Re-sealable Clamshell Pack</i>		
Easy-to-Order Convenience Kits		
11 mm PTFE Red Rubber Crimp Convenience Kit – Pkg/100	1	N9300502
11 mm PTFE Red Rubber/Silver Crimp Assembly (Pkg/100)	1	N9306230
11 mm Clear 2 mL Vial Kit (Pkg/100)	1	N9306231
11 mm PTFE Red Rubber Crimp Convenience Kit – Pkg/500	1	N9300503
11 mm PTFE Red Rubber/Silver Crimp Assembly (Pkg/100)	5	N9306230
11 mm Clear 2 mL Vial Kit (Pkg/100)	5	N9306231
11 mm PTFE/Sil Crimp Convenience Kit – Pkg/100	1	N9300500
11 mm PTFE/Sil Crimp Assembly (Pkg/100)	1	N9306228
11 mm Clear 2 mL Vial Kit (Pkg/100)	1	N9306231
11 mm PTFE/Sil/PTFE Crimp Convenience Kit – Pkg/100	1	N9300501
11 mm PTFE/Sil/PTFE Crimp Assembly (Pkg/100)	1	N9306229
11 mm Clear 2 mL Vial Kit (Pkg/100)	1	N9306231
9 mm PTFE Red Rubber/Screw Cap Convenience Kit – Pkg/100	1	N9300699
9 mm PTFE Red Rubber/Screw Cap Assembly (Pkg/100)	1	N9306200
9 mm Clear 2 mL Vial Kit (Pkg/100)	1	N9306201
9 mm PTFE/Sil Screw Cap Convenience Kit – Pkg/100	1	N9300700
9 mm PTFE/Sil Screw Cap Assembly (Pkg/100)	1	N9306202
9 mm Clear 2 mL Vial Kit (Pkg/100)	1	N9306201



3-Cartridge Gas Purification System

This system provides for high capacity contaminant removal for GC supply gases, and allows for individual cartridge replacement. Cartridge replacement is recommended after processing 19 cylinders of 8 m³ high purity gas (99.997%).

Description	Part No.	System	Capacity	Efficiency
Purification System, 3 Head All Steel System	N9306135	Oxygen	850 cc	< 1 ppb
Cartridge Set, Gas Purification All System Cart (Replacement)	N9306136	Water	12 g	< 10 ppb
Cartridge, Gas Purification Moisture	N9306137	Hydrocarbons	8 g	< 1 ppb
Cartridge, Gas Purification Hydrocarbon	N9306138			
Cartridge, High Capacity O ₂ Replacement	N9306004			
Hardware, Gas Purification Manifold/Mtg	N9306140			
Hardware, Gas Purification Manifold/Mtg for Bench or Wall	N9306014			

Features and Benefits

- All stainless steel cartridges are easy to replace
- High capacity
- Wall bracket gets the system up and out of the way



Advanced Filter System

The Advanced Filter System has high capacity and efficiency levels for oxygen, water and hydrocarbons. The recommended maximum flow rate is 2.0 L/min and 200 psi maximum operating pressure.

Features and Benefits

- Two indicators, for oxygen and moisture
- Easy cartridge replacement with on/off knob
- Double seal construction for safety
- Check valves protect gas lines during cartridge replacement

A polycarbonate shield surrounding the glass indicator section of the filter is sealed, unlike other gas filters, the gas flow is secure even if the glass should break. This redundant sealing system and robust construction provides a new level of security in gas filtration. Maximum flow rate is 2.0 L/min with 200 psi max operating pressure.

Description	Part No.
Advanced Filter System (manifold and mounting hardware included)	N9303963
Replacement Cartridge	N9303964
Manifold and Mounting Hardware	N9306139

Click-On Inline Super Clean™ Gas Purifiers

Stainless Steel Trap Kits



Description	Connector (Qty)	Part No.
Combination: Oxygen/Moisture Trap	1/8-in Brass (2)	N9306108
Combination: Oxygen/Moisture Trap	1/8-in SS (2)	N9306109
Combination: Moisture/Hydrocarbons Trap	1/8-in Brass (2)	N9306117
Combination: Moisture/Hydrocarbons Trap	1/8-in SS (2)	N9306118
Triple: Oxygen/Moisture/Hydrocarbons Trap	1/8-in Brass (2)	N9306110
Triple: Oxygen/Moisture/Hydrocarbons Trap	1/8-in SS (2)	N9306111
Triple Gas Specific (He): Oxygen/Moisture/Hydrocarbons	1/8-in Brass (2)	N9306112
Triple Gas Specific (He): Oxygen/Moisture/Hydrocarbons	1/8-in SS (2)	N9306113

Stainless Steel Traps

Description	Part No.
Moisture Trap	N9306100
Oxygen Trap	N9306101
Hydrocarbons Trap	N9306102
Combination: Oxygen/Moisture Trap	N9306103
Combination: Moisture/Hydrocarbons Trap	N9306105
Triple: Oxygen/Moisture/Hydrocarbons Trap	N9306104
Triple Gas Specific (He): Oxygen/Moisture/Hydrocarbons	N9306106



High-Purity Brass Regulators

PerkinElmer regulators are constructed of high-purity brass barstock and have stainless steel diaphragms and metal-to-metal seals. They are suitable for use with high-purity (>99.995% pure) non-corrosive gases. Regulators terminate in a 1/4 in NPT Swagelok fitting.

Key Benefits

- Barstock body construction
- Stainless steel diaphragms
- Metal-to-metal seals
- Use with high-purity carrier gas

High-Purity Brass Regulators (Dual Stage)

CGA Fitting	Delivery Pressure Delivery Pressure Use	Cylinder Pressure Range (psig)	Gauge (psig)	Gauge (psig)	Part No.
CGA-350	H ₂ and Ar/CH ₄	4 – 100	0 – 200	0 – 4000	09907128
CGA-580	He, Ar, N ₂	4 – 100	0 – 200	0 – 3000	09907127

High-Purity Brass Regulators, *continued*

High-Purity Brass Regulators (Single Stage)

CGA Fitting	Delivery Pressure Delivery Pressure Use	Cylinder Pressure Range (psig)	Gauge (psig)	Gauge (psig)	Part No.
CGA-350	CO, H ₂ and Ar/CH ₄ Mixes	4 – 100	0 – 150	0 – 4000	00230091
CGA-350	CO, H ₂	10 – 200	0 – 400	0 – 4000	00230253
CGA-590*	Air	10 – 200	0 – 400	0 – 4000	00230090

*Supplied with 590-580 Adapter

GC Startup Kits

1/8 in Tubing and Fitting for (3) Gases

N9306304

Description	Qty.
1/8-in Tubing x 50 ft Coil Copper Special Cleaning	1
1/8-in Compression Brass Tee Two Piece Ferrule Brass	3
Tee 1/8-in Com x 1/8-in Comp x 1/4-in fnpt Brass	3
Adjustable Safety Relief Valve Brass 50 – 150 PSI	3
1/4-in fnpt x 1/8-in Comp Fitting Brass	3
1/8-in Port Connector Brass	3
1/8-in Ferrule Brass	3
1/8-in Compression Brass Nut	3
1/8-in Compression Brass Fitting Cap	3
Tubing Cutter 1/8-in Tubing	1
Teflon® Tape	1

1/8 in Tubing and Fitting with One Single Stage Regulator

N9306304

Description	Qty.
1/8-in Tubing x 50 ft Coil Copper Special Cleaning	1
1/8-in Compression Brass Tee Two Piece Ferrule Brass	3
Tee 1/8-in Com x 1/8-in Comp x 1/4-in fnpt Brass	3
Adjustable Safety Relief Valve Brass 50 – 150 PSI	3
1/4-in fnpt x 1/8-in Comp Fitting Brass	3
1/8-in Port Connector Brass	3
1/8-in Ferrule Brass	3
1/8-in Compression Brass Nut	3
1/8-in Compression Brass Fitting Cap	3
Tubing Cutter 1/8-in Tubing	1
Teflon® Tape	1
Single Stage Analytical 0 – 100 psig delivery, CGA 350 (H ₂ , Carbon Monoxide, Ethylene)	1

1/8 in Tubing and Fitting with One Dual Stage Regulator

N9306304

Description	Qty.
1/8-in Tubing x 50 ft Coil Copper Special Cleaning	1
1/8-in Compression Brass Tee Two Piece Ferrule Brass	3
Tee 1/8-in Com x 1/8-in Comp x 1/4-in fnpt Brass	3
Adjustable Safety Relief Valve Brass 50 – 150 PSI	3
1/4-in fnpt x 1/8-in Comp Fitting Brass	3
1/8-in Port Connector Brass	3
1/8-in Ferrule Brass	3
1/8-in Compression Brass Nut	3
1/8-in Compression Brass Fitting Cap	3
Tubing Cutter 1/8-in Tubing	1
Teflon® Tape	1
Dual Stage Analytical 0 – 100 psig delivery, CGA 580 (N ₂ , Argon, He)	1

Velocity (The Economical Choice) and Elite (Premium, Low Bleed, Rugged) Series Capillary Columns

Velocity-1, Velocity-5, and Velocity-Wax, Elite-1, Elite-5, Elite-Wax, Elite-624

PerkinElmer Velocity columns are excellent for standard daily test applications. They combine quality and affordability with reproducible results.

Velocity-1 and Elite-1 – 100% Dimethyl Polysiloxane

General purpose columns with a highly versatile phase that is extremely rugged, exhibiting long column lifetime, and high operating temperatures. Ideal for the analysis of non-polar petrochemical samples, such as detailed hydrocarbon analysis, hydrocarbon gases, petroleum oxygenates, petroleum aromatics, fuels, waxes, oils, sulfur compounds, mercaptans, and carbon disulfide. It also is an excellent phase for solvents, chemicals, flavors, fragrances, essential oils, air toxins, chlorofluorocarbons, arson analysis, pesticides, and hydrocarbons. Thermal stability to 350 °C. Visit our website for additional products.

Description	Dimensions	Part No.
Velocity-1	15 M x 0.25 mm x 0.25 µm	N9306319
Velocity-1	30 M x 0.25 mm x 0.25 µm	N9306312
Velocity-1	30 M x 0.25 mm x 1.00 µm	N9306323
Velocity-1	30 M x 0.32 mm x 0.25 µm	N9306318
Velocity-1	60 M x 0.25 mm x 0.25 µm	N9306320
Elite-1	15 M x 0.53 mm x 0.15 µm	N9316045
Elite-1	30 M x 0.32 mm x 3.00 µm	N9316025
Elite-1	30 M x 0.25 mm x 0.25 µm	N9316010
Elite-1	60 M x 0.25 mm x 0.25 µm	N9316013

Velocity-5 and Elite-5 – 5% Diphenyl and 95% Dimethyl Polysiloxane

This column is ideal for general purpose analysis of drugs, pesticides, hydrocarbons, essential oils and semi-volatiles and solvent impurities. Low polarity phase with thermal stability to 350 °C. (Elite – 5HT 400 °C.)

Description	Dimensions	Part No.
Velocity-5	15 M x 0.32 mm x 0.25 µm	N9306325
Velocity-5	30 M x 0.25 mm x 0.25 µm	N9306311
Velocity-5	30 M x 0.32 mm x 0.25 µm	N9306313
Velocity-5	30 M x 0.32 mm x 1.00 µm	N9306316
Elite-5	30 M x 0.25 mm x 25 µm	N9316076
Elite-5	30 M x 0.32 mm x 25 µm	N9316086
Elite-5HT	15 M x 0.32 mm x 0.10 µm	N9316274

Velocity-Wax and Elite-Wax – Polyethylene Glycol

This column is ideal for intermediate to high polarity compounds. Thermal stability to 250 °C.

Description	Dimensions	Part No.
Velocity-Wax	30 M x 0.32 mm x 0.25 µm	N9306314
Velocity-Wax	30 M x 0.25 mm x 0.25 µm	N9306315
Elite-Wax	30 M x 0.25 mm x 0.25 µm	N9316403
Elite-Wax	30 M x 0.32 mm x 0.25 µm	N9316412
Elite-Wax	30 M x 0.32 mm x 0.50 µm	N9316413

Specialty Phase for Analysis of Volatiles

Description	Dimensions	Part No.
Elite-624	30 M x 0.53 mm x 3.0 µm	N9316207
Elite-624	30 M x 0.25 mm x 1.4 µm	N9316201
Elite-624	30 M x 0.32 mm x 1.80 µm	N9316203

PerkinElmer – The Clear Choice in Chromatography Consumables

- Genuine PerkinElmer consumables products
- 100% Tested and Approved
- Guaranteed compatibility

Portable Gas Leak Detector

The new PerkinElmer compact handheld electronic gas leak detector is the ideal solution for detecting gas leaks in your Gas Chromatography systems. Leaks in your system waste gas and can cause detector noise, baseline instability, and shorter column life. This portable



unit detects minute leaks of any gas with thermal conductivity different from air. The reference gas inlet draws in ambient air for comparison to air drawn into the sample probe. A leak is detected by both LED bar graph display and audible alarm.

N9306089

MiniTemp MT4 Non-contact Temperature Measurement with Laser Sighting

The MiniTemp MT4 is capable of displaying thermal measurement readings using easy point and shoot infrared technology in a pocket size configuration. This high accuracy unit is great for instrument thermal test confirmation, including GC injector port and detector measurements,



thermostatted LC vials, and enzymatic hydrolysis baths. The MT4 is only 6 inches in height, light weight, and fits in your pocket. It provides easy resolution to common temperature measurement requirements in the lab including difficult to reach locations, with a red laser dot sight indicator!

N9306074

Digital Bubble Flow Meter

The PerkinElmer Model 520 is a volumetric flow meter. It can measure the flow rate of any gas or combination of gases, such as air, without adjustment. The flow meter has a digital display



and a single push-button input. It is made of stainless steel and anodized aluminum. It comes with a certificate of calibration, and is accurate to $\pm 3\%$. The Model 520 can measure flow rates of 0.5 to 500 mL/min.

N9302974

PerkinElmer FlowMark Electronic Flowmeter

PerkinElmer's FlowMark™ flowmeter is specifically designed for use with GC instruments. The probe is applied directly to the gas flow stream and the measured flow rate is presented on the LCD screen. Units of flow are measured in mL/min. This unit provides continuous



real-time measurements of gas streams ranging from 0.50 mL/min to 500 mL/min. Because the technology uses volumetric flow measurement, the unit is compatible with all laboratory gases. The flowmeter is designed to measure clean, dry non-corrosive gases.

N9307086

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 more details, call your local representative or visit www.perkinelmer.com/gcsupplies.

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

009737A_01