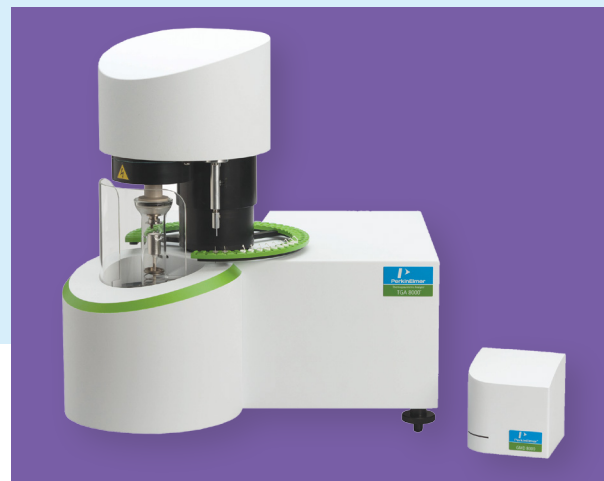


## Thermal Analysis

GMD 8000  
Gas Mixing Device

The GMD 8000 allows mixing of up to three gases, delivering a custom gas environment to the TGA 8000™ or other devices. The GMD 8000 has three gas inputs, each with its own mass flow controller (MFC). Gas mixing is customized via a Windows® PC-based software program. The customized mixed gas exits from a single outlet and is delivered to the sample through the TGA 8000 reactive gas line or a non-controlled input port of other devices.

## GMD 8000 Specifications

Design		Compact stand-alone controller for delivering a custom gas environment for analytical instruments. Includes three gas inputs, each with its own mass flow controller (MFC), and a single outlet for delivery of a user defined custom gas mixture.
Gases		Air, Argon, Carbon dioxide, Helium, Nitrogen, Oxygen <sup>1</sup>
Gas Management	Gas Input	3 mass flow-controlled inputs (software controlled input ratio)
	Gas Output	1 output
	Input Pressure	15 - 90 PSI
	Output Flow	Up to 600 ml/min (1 ml/min increments) <sup>2</sup>
	Accuracy	2% of set value or 2 ml/min (using most common gases)
User Interface	Repeatability	0.5%
	Software	Microsoft® Windows® 7; English, Japanese, Chinese languages
	Features	Single program mixing; Switch from one gas mixture to a second; Ramp between two gas mixtures
Communication	Display	LED status indicator provides setup and runtime monitoring
	PC interface	USB
System	External interface	SeaLINK® (trigger) interface to control ON/OFF of gas flow
	Dimensions (H x W x D)	100 mm x 100 mm x 100 mm
	Weight	1 Kg (2 Lbs)
	Power	100 - 240 Volt, 50/60 Hz



1. Input gas must be dry and filtered. Not designed for humid gases. Hydrogen gas or corrosive gases cannot be used.
2. With input ratio 33%/33%/33% and each input flow is greater than 200 ml/min (200 sccm (20 °C, 1 atm))

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

  
**PerkinElmer**<sup>®</sup>  
For the Better

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2015 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.