## **TURBOMATRIX™ HEADSPACE TRAP SAMPLERS**

## Up to 100 times lower detection limits



The clear choice in gas chromatography. www.perkinelmer.com/GC



For a complete listing of our global offices, visit www.perkinelmer.com/lasoffices

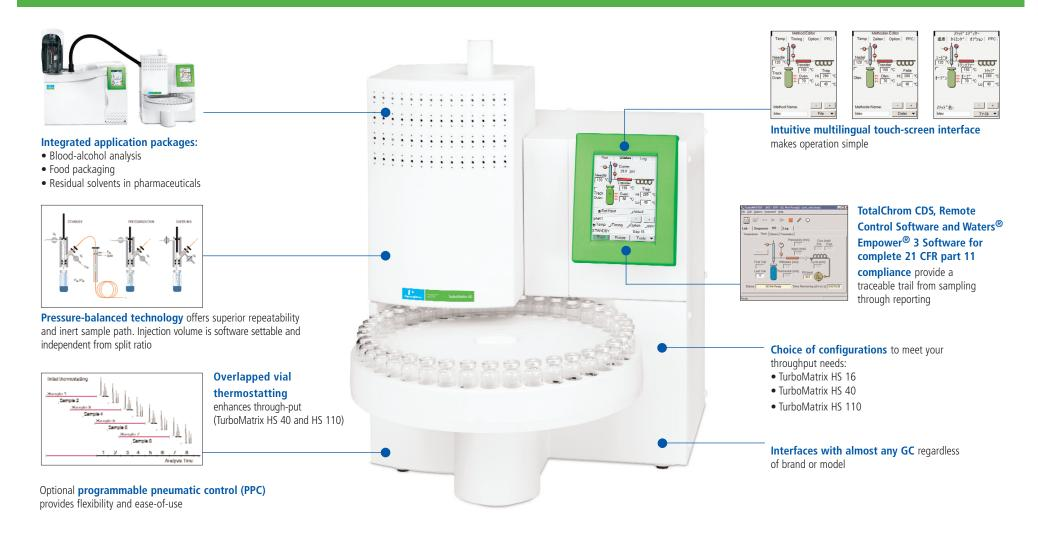
©2007-2016 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.

PerkinElmer, Inc. 940 Winter Street Waltham, MA 02451 USA Phone: (800) 762-4000 or

(+1) 203-925-4602 www.perkinelmer.com

## **TURBOMATRIX™ HEADSPACE SAMPLERS**

The clear choice for any GC or GC/MS volatile-analysis system



The clear choice in gas chromatography. www.perkinelmer.com/GC

## PerkinElmer, Inc. 940 Winter Street



For a complete listing of our global offices, visit www.perkinelmer.com/lasoffices

©2007-2016 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.

Waltham, MA 02451 USA Phone: (800) 762-4000 or

(+1) 203-925-4602 www.perkinelmer.com