



## Efficient Rodent Anesthesia System with Active Scavenging Capabilities



### RAS-4 Rodent Anesthesia System

Part # CLS146737

The RAS-4 Rodent Anesthesia System enables real-time *in vivo* imaging to monitor and record molecular and genetic activity in mice and rats using isoflurane gas. The RAS-4 Rodent Anesthesia System is compatible with PerkinElmer's suite of *in vivo* imaging systems. The system facilitates efficient gas delivery via four delivery circuits while minimizing excess gas exposure to lab personnel with active scavenging capabilities.

## 360 Degree Surface Mapping



### MVI-2 Multi-View Imager Kit

Part # CLS143807

The MVI-2 Multi-view imager enables complete 360 degree 2D surface mapping of two mice simultaneously for the identification, localization, and quantification of whole body imaging of probes.

Please visit our website for a complete list of accessories <https://www.perkinelmer.com>.

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)



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

Copyright © 2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

69324 (57955A) PKI

**DRAFT**