

# DATA SECURITY AND PRIVACY IN THE CLOUD

## Introduction

PerkinElmer Signals™ is an agile data warehouse designed for storing assay and clinical research data in a common, cloud-based warehouse. The core of Signals is a highly flexible and scalable Measurement Store that users and administrators can adapt to store experimental results from everything from protein-based assays and cellular imaging assays to test results from clinical trials. Easy-to-use query interfaces empower scientists with access to a wide range of assay data without having to rely on data managers.

## Security and Privacy

A recent report by Gartner titled “Clouds Are Secure: Are You Using Them Securely?” makes the point that cloud-based infrastructure tends to be at least as secure as infrastructure located on-premise. However, the report goes on to say that the safe use of cloud infrastructure “requires new organizational policies, skills and activities”.

We take our responsibility towards the safekeeping of our customer’s data very seriously. As a cloud-based, hosted



SaaS solution, PerkinElmer Signals™ has been designed from the beginning with a strong emphasis on maintaining data safety and privacy. This starts with the use of strong encryption for all data, regardless of whether they are in rest or transit and ends with a single tenant deployment model using a Virtual Private Cloud (VPC) for secure access and separation between customers. When dealing with sensitive research data, and in particular medical data, the physical location of the data is often of paramount importance. By offering PerkinElmer Signals™ deployment in multiple geographic regions we help our customers meet regulatory and legal requirements.

Please contact us as [Signals@PerkinElmer.com](mailto:Signals@PerkinElmer.com) for further information.

Requirement	Details
Hosting Service	PerkinElmer Signals™ is hosted in the Amazon Elastic Compute Cloud® (EC2) and is offered as a Software as a Service solution by PerkinElmer (with Amazon acting as an Infrastructure as a Service provider).
Data Segmentation	PerkinElmer Signals™ is a single tenant solution with strict separation between customer instances. There is no sharing of resources and/or components between customer instances.
Data Localization	PerkinElmer Signals™ utilizes Amazon EC2® Regions to provide optimal performance with low latency and ensure that data are confined to specific regions and helps meet regulatory and legal requirements.
Data Encryption	PerkinElmer Signals™ leverages Amazon encryption to protect all customer data using strong encryption (AES 256) when at rest. The encryption uses customer-specific encryption key to guarantee access and privacy.
Network Configuration	All communication with PerkinElmer Signals™ utilizes secure HTTPS connections. A Virtual Private Cloud is set up for every customer to ensure secure access from the customer’s internal networks.
User Management	PerkinElmer Signals™ can be configured to use Single Sign-On (SSO) systems for authentication and authorization (incl. SAML 2.0, OpenID 2.0, and OAUTH 2.0).
Information Security	Amazon EC2® is certified to comply with ISO 27001; PerkinElmer generally follows the guidelines of ISO 27001 but is not certified at this time.

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