



# PERKINELMER SOLUTIONS FOR PHARMACOVIGILANCE AND DRUG SAFETY

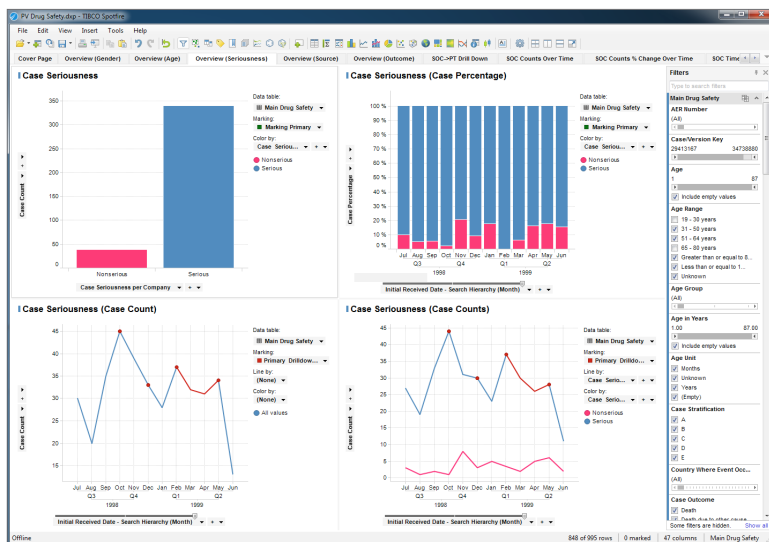
## Strategic Transformation, Time and Costs Savings

Built on the best-in-class scientific analytics and data discovery platform, TIBCO Spotfire®, PerkinElmer's innovative solution for Pharmacovigilance\* enables teams to strategically transform their medical review, safety review, and compliance monitoring processes, gain new insights leading to faster identification of safety risks early in development, and reduce time and costs preparing for safety team reviews. Customers also benefit from streamlined day to day operations and better support reporting activities like PSUR or PBRER creation.

PerkinElmer's solution provides exploratory analysis, so in a fast and easy way, users can freely explore and visually interact with the data, without needing to know the questions they want to ask in advance, as required by traditional reporting out of safety systems. Traditional reporting environments require IT assistance and require you to know what you want to see and then ask for it (query), resulting in time delays and returning only what you asked for in the form of reports, for example Argus with OBIEE. ("Query -> Generate Report -> Next Question -> Repeat.")

PerkinElmer's solution replaces the reporting out of safety systems that leverage standard BI tools, breaking the query, wait, evaluate cycle. Our innovative PV data discovery and analytics solution empowers all users to go directly against the data, intuitively visually interact with it, perform ad hoc analysis, and gain insights previously not possible, requiring no IT assistance and involving no time delays. Users can freely explore certain analysis factors over time, not available in standard safety systems, such as "serial changes" or changes of certain analysis factors over time such as age, race, sex, volume of events over time. Our solution also provides valuable compound case analysis or similar-compound trend analysis.

\*The abbreviation, PV, will be used in place of Pharmacovigilance going forward in this data sheet



PerkinElmer is the exclusive global distributor of the TIBCO Spotfire® platform for certain scientific and clinical R&D applications.

## AT A GLANCE

PerkinElmer's PV solution is built on TIBCO Spotfire®, which is recognized as the "de facto end user of choice for the strategic assessment of clinical data." It is the life science's leading analytics and data discovery software, used by the top pharmaceuticals, biotech, med device companies, and CROs globally.

## BENEFITS

### Fastest to Actionable Insight

Increase agility by providing clinical users with unmatched speed and flexibility to anticipate, answer, and act on answers to PV questions.

### Visibility into the Unknown

Uncover trends and patterns that represent risk or opportunity through intuitive visualizations, analytic dashboards, and applications. Connect to data sources that include industry standard systems like Oracle's Argus®, ArisGlobal's Arisg® and others.

### Self-Service Discovery

Empower a broad spectrum of PV, clinical, and other users with self-service discovery to go directly against the data, answer their own questions, and free Biostatistics and IT groups for higher value projects.

### Universal Adaptability

Empower the broadest spectrum of PV, clinical, and other users to freely explore clinical data, identify trends representing risks and opportunities earlier, and be able to dig deeper into the data without reliance on Biostats and IT groups. This allows organizations to gain a faster review of clinical data, provide better data quality assessment, and make for a drastically improved and quicker process.



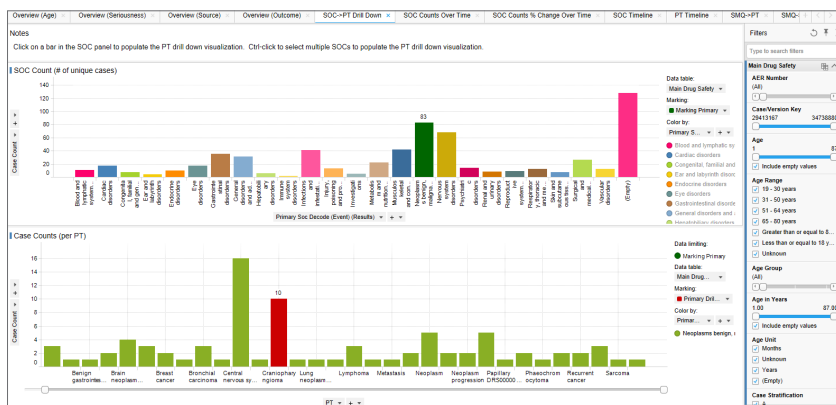
# ATTRIBUTES AND CAPABILITIES

## Hide the Complexity



Easy to understand visualizations such as Serial Changes (changes of certain analysis factors over time such as age, race, sex, volume of events over time) can be leveraged without getting bogged down in data aggregation and statistical modeling.

## SOC and PT Analysis



System Organ Classes can be explored, to populate Patient drill down data visualizations. Further SOC analysis over time, % changes over time, and Timelines can be freely explored and intuitively analyzed in more detail.

**TIBCO Spotfire® is recognized as the “de facto end user of choice for the strategic assessment of clinical data.”**

- IDC Health Insights, “Familiar and User-Friendly: TIBCO Spotfire® Owns the Front End in the Life Sciences”

Leveraging data from industry standard systems like Oracle’s Argus®, ArisGlobal’s Arisg® and others, PerkinElmer’s solution for PV streamlines day to day operations, resulting in faster and enhanced identification of safety issues associated with pharmaceutical, biotech, or medical device products, while minimizing risks that may come to patients and maintaining good standing with regulators.

## RETURN ON INVESTMENT FROM CUSTOMERS:

- Reduce Safety Review Preparation Time from 3 Weeks to 10 Minutes
- Pivot from Case to Event from 1-2 Days to < 1 Minute
- Answer the Next Question from 1 Week to “in the room”

To learn more about our solutions for clinical development, please visit [www.perkinelmer.com/ClinicalInformatics](http://www.perkinelmer.com/ClinicalInformatics)

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.