



## PERKINELMER SOLUTIONS FOR TRIAL OPERATIONS



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

### Achieve Clinical Trial Operations Excellence

PerkinElmer's innovative solution gives clinical operations organizations the unprecedented ability to reduce the time, cost, and risk of running clinical trials while accelerating time to market for products across the entire portfolio of clinical candidates. The solution provides timely information, enabling clinical study teams to make decisions based on live, interactive scorecards, and gain insights that enable optimal study startup, budget, and milestone tracking.

The solution provides a strategic transformation for clinical operations professionals who may rely on inflexible BI tools and standard reporting processes, or spend vast amounts of time compiling spreadsheets and manipulating data to create charts and graphs.

Easily rendered visualizations enable the rapid determination of the current state of a program, trial or site status which helps teams gain insight into areas of potential issues, allow the understanding of the source cause of issues, and improve the management of clinical performance through metrics and KPIs.

### AT A GLANCE

PerkinElmer's Trial Operations solution is built on TIBCO Spotfire®, which is recognized as the "de facto end user of choice in the strategic assessment of clinical data."<sup>1</sup> It is the life science's leading analytics and data discovery software, used by top pharmaceutical, biotech, medical device, and contract research organizations globally.

### BENEFITS

#### Fastest to Actionable Insight

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

#### Visibility into the Unknown

Uncover trends and patterns that represent risks or opportunities through intuitive visualizations, analytic dashboards, and applications.

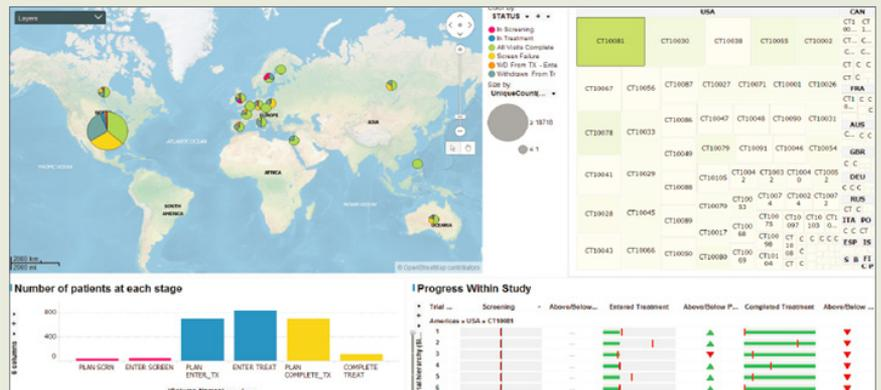
#### Self-Service Discovery

Empower a broad spectrum of Program Managers, Country Managers, Study Managers,

CRA's, and others with no advanced technical, computer, or statistical skills to go directly against the data, freely explore, and answer their own questions.

#### Universal Adaptability

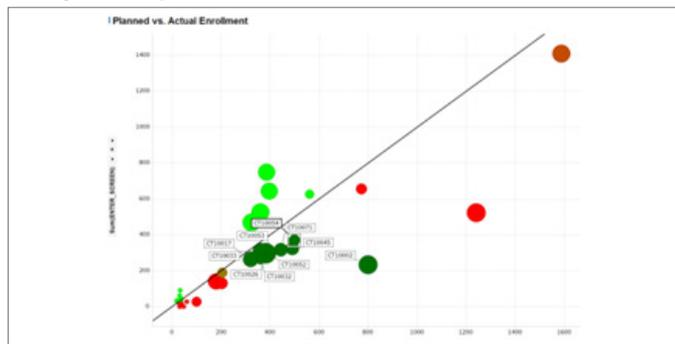
Empower a broad spectrum of users to freely explore operational data, identify trends representing risks and opportunities earlier, and be able to dig deep into the data. This allows organizations to gain a faster review of operational data and provide better study oversight.



Progress within Trials Across Clinical Development Portfolio

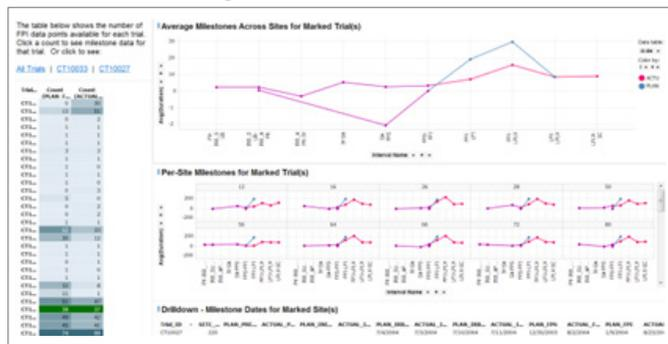
# ATTRIBUTES AND CAPABILITIES

## Study Startup



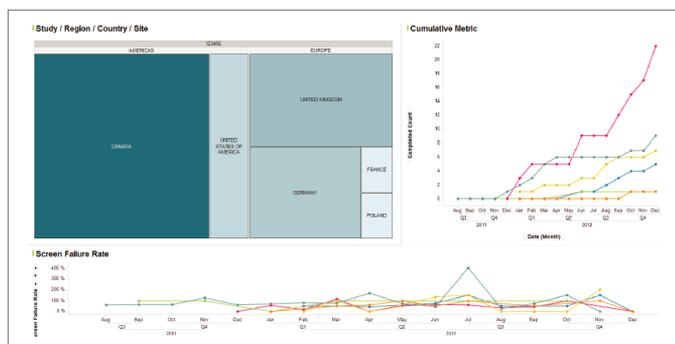
Tasks associated with study startup including site identification and feasibility and planning for patient recruitment can be enhanced.

## Milestone Tracking



Companies can see how trials are progressing compared to plan for key milestones such as Investigator Review Board approvals, first patient in, last patient visit, etc.

## Recruitment



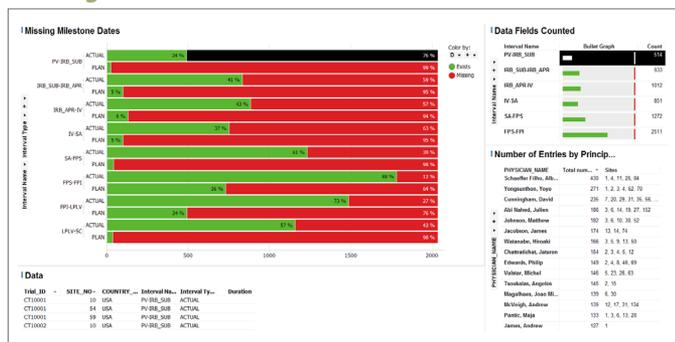
Patient recruitment pace can be tracked and explored in detail, enabling shortened or on-time study timelines.

## Withdrawal/Failure Rates



Withdrawal rates vs. screen failure rates for each site across all trials can easily be viewed and further investigated.

## Missing Data



The percentage of milestone data points present versus missing can easily be viewed along with the distribution of duration for each milestone interval.

*It is the responsibility of the end user to install and validate PerkinElmer solutions within the end users quality systems in accordance with their policies and standard operating procedures.*

## Budget Tracking

Planned versus actual budgets can be tracked at the study, country, and site level showing the cumulative budget versus the actual budget as well as a drill down to each payment made to each site.

## References

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

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

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.