



TIBCO SPOTFIRE X: NOW WITH EVEN MORE INTELLIGENCE

Around the world, scientists and engineers are leveraging the power and possibilities of TIBCO Spotfire® analytics to uncover deeper insights into their data. And with the new TIBCO Spotfire® X platform, you and your organization can better seize on new opportunities and mitigate risks with maximum flexibility. From interactive dashboards and visualizations to predictive analytics, TIBCO Spotfire® software helps you explore data and reach actionable insights quickly, without having to rely heavily on IT. You can analyze disparate data from multiple sources and create a better, more comprehensive picture of what's happening right now.

THE POWER OF X

With TIBCO Spotfire® X, we've harnessed the power of artificial intelligence (AI) to boost productivity even further – so you can reach those powerful insights that much faster. And together with machine learning capabilities, this latest version of the platform enables you to quickly access, explore, and analyze data from screening, preclinical, translational, and even pharmacovigilance research with anyone, anywhere. Visual analytics; data wrangling; and predictive, location, and streaming analytics: it's all possible with the TIBCO Spotfire® X platform.



KICKSTART YOUR JOURNEY TO BETTER ANALYTICS

With artificial intelligence and machine learning capabilities, the TIBCO Spotfire® X platform is helping scientists around the globe to access and analyze data from a whole host of sources, faster than ever before. Here are some resources to help you better understand this remarkable platform – and utilize it in your research.

TIBCO Spotfire® X Overview

This video provides a quick overview of the platform's AI and machine learning updates, which give you the ability to quickly access, explore, and analyze data from a wide variety of sources.

Gartner Magic Quadrant 2019

A Magic Quadrant delivers a graphical competitive positioning of four types of technology providers in high-growth markets with differentiated providers. TIBCO Spotfire® technology was designated a Magic Quadrant Visionary.

Latest Platform Features and Benefits

Our webinar shows you all the features and benefits of the TIBCO Spotfire® X platform.

TIBCO Spotfire® X for Macs

This webinar demonstrates TIBCO Spotfire® X on Macs, with tips and tricks about research and data analysis with the TIBCO Spotfire® Mac client, ad hoc exploration of early-phase clinical-trial data on the Mac, and much more.

Life Sciences Use Cases

TIBCO Spotfire® X is the ideal platform for visualizing results in a wide array of life sciences applications. Get the inside story in this 30-minute presentation.

TIBCO Spotfire® Community

Read forum posts (or contribute your own posts) on our TIBCO Spotfire® Wiki.

TIBCO Spotfire® Training

Looking to acquire new skills with the TIBCO Spotfire® X platform? Sign up for a training course.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

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

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

