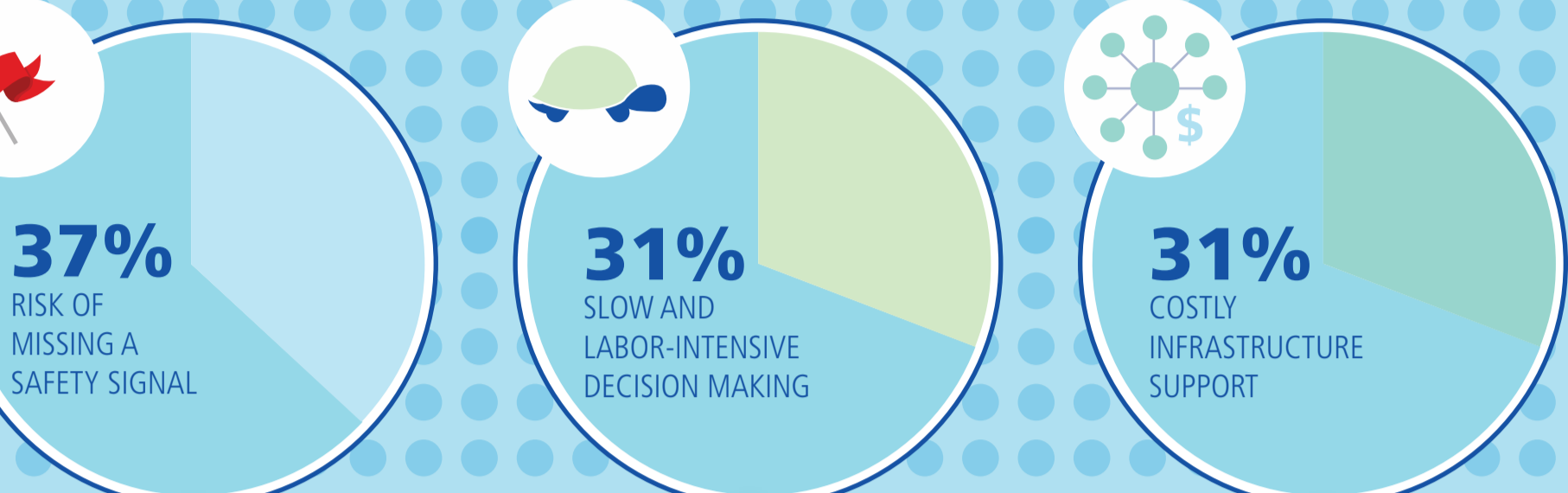


Never miss a safety signal.

As the complexity of clinical trial data exponentially grows, so does the pressure on clinical development teams not to miss a safety signal. To help teams meet this challenge, PerkinElmer, a leading innovator in life sciences introduces Signals Medical Review to provide medical monitors with fast, reliable, and actionable insights through a combination of advanced data technologies and purpose-built workflows.

What are the Greatest Challenges in Medical Review Today?

In a recent poll of clinical development professionals, the top 3 challenges in Medical Review are:



Signals Medical Review addresses all three pain points and more, through the combination of streamlined workflows and advanced analytics.

LEADING TO THREE KEY OUTCOMES:

GET SUBMISSION-READY FASTER

Medical monitors endure sluggish safety data reviews due to the lack of visibility of data from different systems. Signals Medical Review solves this by delivering personalized landing pages, alerts on specific cohorts of interest, curated analysis visualizations, and the ability to track line listing review – effectively trimming up to 5 days per month / per trial on medical data review.

FACT:

Decreasing time to market by 1 month can save a company up to \$50 million in a year.

REDUCE REJECTED SUBMISSIONS OR MARKET WITHDRAWAL

Powered by TIBCO Spotfire® visual analytics, Signals Medical Review allows safety teams to find immediate answers to safety questions from the FDA. This will get you way ahead of competitors who still rely on SAS listings, without compromising patient safety.

FACT:

Data shows that drug recalls can cost companies \$650 million or more

STREAMLINE DATA PROVISIONING, AGGREGATION AND REVIEW

Data aggregation is an incredibly taxing process, especially due to multiple systems, different formats, and high costs. Signals Medical Review unifies all these steps through its transformative, end-to-end cloud solution – from data analysis to action. It automates over 50% of all the manual steps involved, boosting productivity by 25% while ensuring higher accuracy and lower risk of human error.

FACT:

Simplifying study setup can save companies \$3-8 million a year

THE RESULTS

PERKINELMER SIGNALS MEDICAL REVIEW COULD SET THE BAR IN THE MEDICAL REVIEW INDUSTRY, AND THESE FIGURES CAN PROVE IT:

Biostatistics receive **20-40%** less ad-hoc requests

With the platform, days per month per trial is reduced by **2-5 DAYS**

Time in database unlocks is reduced by around **50%**

Time to database lock reduced by **10-30%**

IT and customer support costs are reduced by around **67%**