

A DATA CATALOG FOR BUSINESS INTELLIGENCE & ANALYTICS

Spend Less Time Gathering Data and More Time on Analytics

Analysts may have self-service BI, but still spend far too much time provisioning the right data for BI and advanced analytics tools. Forrester Research estimates that "...80% of any Business Intelligence initiative is spent on data integration." Given such a slow and inefficient data discovery process, many BI users continue to burden IT with multiple requests to gain access to additional data sets. Alternatively, they develop analyses based on incomplete information. **This is not agile BI.**

Attvio's Semantic Data Catalog provides a fast and easy way to find and understand data, wherever it lives. Attvio delivers the right information, in context, directly into BI and analytics tools so analysts can spend less time looking for data sources and more time developing data-driven insights.



Immediate Discovery

Do you spend too much time gathering data sources for analysis? With Attvio, you can instantly find the right data sets from all of your enterprise information, so you can identify the data that is most relevant. It's as easy as Google, and takes only minutes to find the right data sources.

The Attvio Difference

Attvio's foundational technology powers four unique differentiators.

BIG DATA SEARCH & DISCOVERY

Provides immediate visibility into the right information

SEMANTIC ENRICHMENT

Automatically discovers the meaning and relationships in your information ecosystem

DATA UNIFICATION

Unifies disparate information across silos to deliver unique insights

AGILE ENTERPRISE PLATFORM

Secure, scalable foundation for delivering data-driven applications

Attvio's Semantic Data Catalog

The Attvio Semantic Data Catalog helps analysts, data scientists, and data managers find, understand, unify, and provision the right information for analytics. With an easy shopping cart experience, data consumers can get the data they need, instantly.



360° View of All Relevant Information

What if you could instantly see the connections between your data sources?

Attivio gives you the power to build a unified model from disparate data sources. With both recommended and user-defined connections between data sources, you control the model. You then provision the model to your BI or analytics tool. It's that easy.



Self-Service Access

Do you have all of the data you need for analysis at your fingertips? The process of getting the best data for analysis involves time-consuming back-and-forth cycles between analysts and IT. Attivio gives you instant access to all of your enterprise information – structured and unstructured – so you have complete confidence that you are using the right data, in context.

If you spend too much time gathering data sources, and not enough time analyzing data, **you'll want to see how Attivio can help you get the right data in a fraction of the time.**

Deepen Your Insights with Unstructured Content

Whether it's customer comments, field notes, social posts, or other unstructured information, companies in any industry can benefit from insights available in text-based content. With its Text Analytics technology, Attivio adds structure to unstructured content so that disparate sources can be correlated, modeled, and provisioned, quickly and easily.

About Attivio

Attivio provides software that empowers its customers to get their hands on the right data and to work with it to quickly get it to the point of informing decisions. Visit us online at www.attivio.com.

About PerkinElmer

PerkinElmer, Inc. is a global leader committed to innovating for a healthier world. Our dedicated team of over 9,000 employees worldwide are passionate about providing customers with an unmatched experience as they help solve critical issues especially impacting the diagnostics, discovery and analytical solutions markets. Our innovative detection, imaging, informatics and service capabilities, combined with deep market knowledge and expertise, help customers gain earlier and more accurate insights to improve lives and the world around us. PerkinElmer is the exclusive distributor for both Attivio® and TIBCO Spotfire® for scientific and clinical R&D applications.