

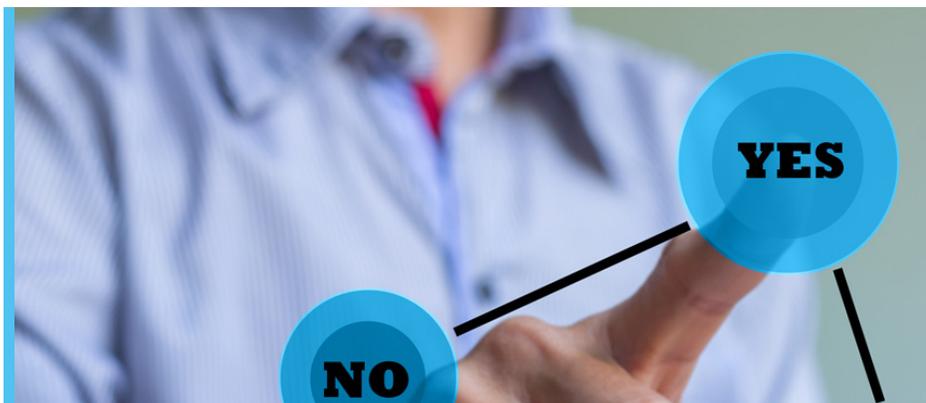
# COGNITIVE SEARCH AND INSIGHT

Drive innovation, operational efficiencies, and improved business outcomes.  
Act with certainty.

It takes a long time to find the right information to make decisions, often too long. According to Findwise, less than 50% of employees have the ability to search across siloed content. In fact, in a recent survey, Capgemini found that 56% of companies cited information silos as the biggest impediment to effective decision-making.

While enterprise architects struggle with current search technologies to tune relevance, add new sources, and enforce security, end users spend more time looking for the right information than actually analyzing it. As a result, decisions typically wind up being made based only on the information that is easy to access – usually just structured data - versus using all the right information from across the enterprise as the basis.

*No one would drive a car if they could see out of only 20% of the windshield, would they? So why drive your business by making decisions using only 20% of relevant information?*



Make better decisions. Act with certainty.

Attivio's cognitive search and insight solutions provide industry-leading natural language processing, machine learning, analytics, and knowledge graphing capabilities at scale to help you drive innovation, operational efficiencies, and improved business outcomes enterprise-wide:

## The Attivio Difference

Attivio's foundational technology powers four unique differentiators.

### INFORMATION INTELLIGENCE

Reveal relationships and insights across structured and unstructured silos through contextual enrichment and semantic understanding

### USER EMPOWERMENT

Anticipate individual information needs and boost productivity by enabling people to find critical information at the point of need

### SECURE ENTERPRISE PLATFORM

Adapt to the most complex data and IT ecosystems, while enforcing rigorous security policies at global scale

### PROVEN EXECUTION

Turn data and content into strategic advantage on the most complete and proven cognitive platform

With Attivio, organizations can quickly gain critical insight and build domain-specific applications, while realizing a rapid return on investment. Attivio's cognitive search and insight far surpasses the expectations of traditional enterprise search by providing game-changing capabilities, including:

- **Natural Language Processing:** Improves human-computer interactions to accelerate knowledge discovery and collaboration.
- **Machine Learning:** Provides improved context for decisions through continuous updating of actions taken.
- **Text Analytics:** Extracts content and metadata from 500+ formats, detect and index entities, and extract key phrases to automatically surface insights in massive volumes of unstructured content.
- **Active Security:** Provides real-time document & field-level security to meet the most complex security needs. A patented, best-in-class approach to security.
- **Predictive Alerting:** Detects outliers in patterns, correlations, and predictions.
- **Relevancy:** Analyzes 100+ signals in data and user behaviors to deliver the most relevant documents and data for a user's query.
- **Recommendations:** Focuses information managers on the most relevant factors and the right answers for decision-making.
- **Simplified UXD:** Optimizes the end user experience with a highly personalized and intuitive UI designed for non-technical users, while masking the sophisticated security calculations behind each query. Responsive design supports all screen sizes and device types.
- **Connectors:** Connects to 100+ enterprise data sources across 500+ file formats – including cloud-based file stores and web crawls that gather and pre-analyze web copy during ingestion – to unify data and uncover insights from across the enterprise.
- **Advanced Linguistics Support:** Supports 45 languages with full linguistic analysis to improve the speed and accuracy of search applications globally. Automatic language detection and entity extraction across all languages.
- **Robust Workflow Capabilities:** Manages data and query transformations through a powerful, flexible, and comprehensive interface. 300+ OOTB stages to transform data and queries, including OCR, entity extraction, key phrase extraction, and entity-level sentiment. Automatically works with 550+ file formats.
- **Query-time JOINS:** Selects relationships across tables, connecting common fields to represent one-to-one, one-to-many, and many-to-many relationships.

## About Attivio

---

Attivio is the leading cognitive search and insight company. Our Fortune 500 clients rely on us to drive innovation, operational efficiencies, and improve business outcomes. Our solutions provide industry-leading natural language processing, machine learning, analytics, and knowledge graphing capabilities at scale. Let Attivio empower you to act with certainty. Visit us online at [www.attivio.com](http://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.