



Proven Methods. Superior Support

Whatever the scale or nature of your purchasing needs, an e-enabled procurement solution from Sigma-Aldrich can result in Significant savings, and total peace of mind for your organization.

The future of e-commerce is being provided by Sigma-Aldrich – and that means empowerment is here to stay.

FAQ's

What are Sigma-Aldrich's e-commerce capabilities?

Sigma-Aldrich offers various solutions that accommodate B2B/ERP platforms, "custom" websites, EDI, catalogue files and electronic invoicing.

What efficiencies can be gained by implementing an Electronic Solution with Sigma Aldrich?

- Streamlined ordering process-eliminating paper transactions
- Time savings
- Order Tracking
- Product availability and contract pricing
- Reduced error rates

Does the Sigma-Aldrich Website support work-flow-approval process?

Our customized PIPELINE solution will support workflow based on various transaction levels

What support is available for implementing an e-solution?

You will have a dedicated implementation Analyst available for pre and post implementation, who will help you seamlessly, manage the implementation process.

What systems does Sigma-Aldrich integrate with?

Sigma-Aldrich will evaluate any integration request and access if we are able to integrate with the technology used. We currently integrate using the following technologies: CXML, XCBL, OAG, OCI, SAP, iDOC XML, EDI and custom XML(s).