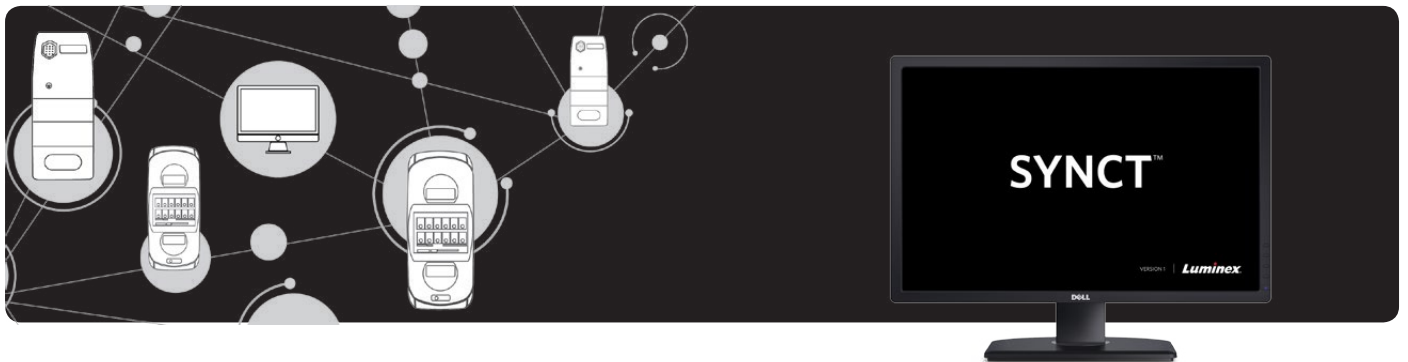




SYNCT™

Software



Welcome to the New Way to Work

SYNCT™ Software empowers laboratorians to manage data from any place in the lab. It is a molecular middleware that provides flexible, customizable workflows by connecting up to four ARIES® or NxTAG®-enabled MAGPIX® Systems.

SYNCT Software enables labs to manage orders from laboratory information systems (LIS), view test results, prepare test reports, perform administrative functions, and maintain security control.

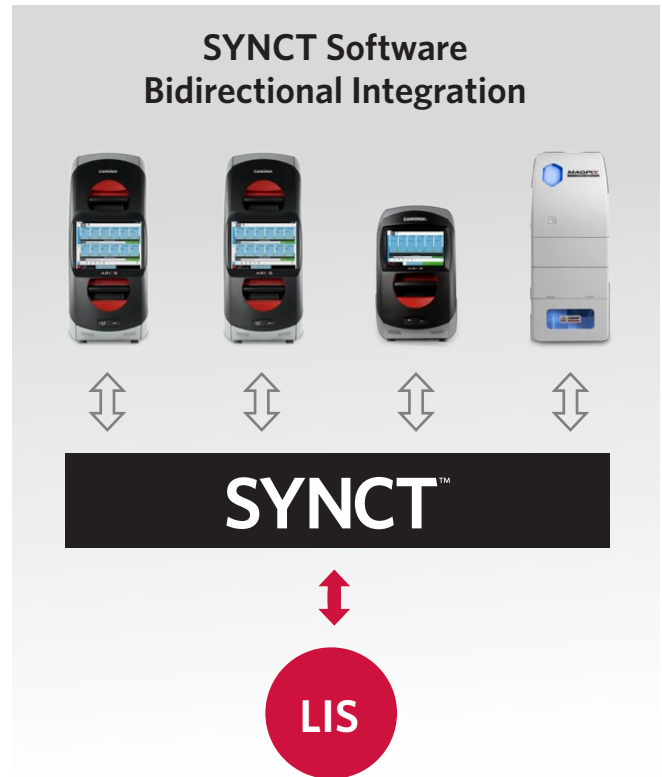
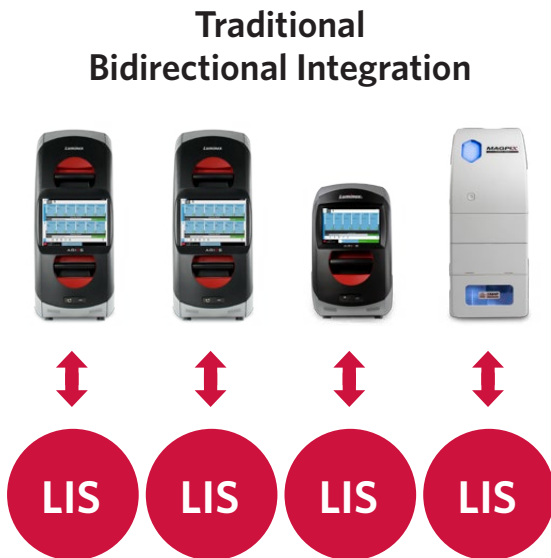
This software offers:

- Bidirectional interface, enabling the transfer of both orders and results
- One connection between SYNCT Software and your LIS, allowing labs to easily connect multiple instruments to the LIS
- Results monitoring and management from anywhere in the lab
- The ability to associate cassettes to orders from anywhere in the lab

To learn more about the benefits of SYNCT Software, please contact your Luminex sales representative today.

How to Connect the ARIES® System to Your LIS

One connection through SYNCT Software to your LIS enables your laboratory to easily connect multiple instruments to your network. It's never been easier.



Luminex will provide you with both an integration guide and access to a LIS expert to guide you through the setup process. To get started, it only takes three easy steps:

1. Locate the assay codes in the provided product insert, and enter the codes into your LIS.
2. Connect the ARIES® System to your laboratory network.
3. Integrate SYNCT to your LIS using Luminex's integration guide once, and then connect all the ARIES® Systems and NxTAG-enabled MAGPIX Systems with a few clicks of a mouse.

Luminex
complexity simplified.

To learn more, please visit: www.luminexcorp.com

For In Vitro Diagnostic Use. Products are region specific and may not be approved in some countries/regions. Please contact Luminex to obtain the appropriate product information for your country of residence. The ARIES® instrument is a Class 1(I) laser product. Validation of the LIS compatibility must be performed by the end user.

©2017 Luminex Corporation. All rights reserved. Luminex ARIES, NxTAG, and MAGPIX are trademarks of Luminex Corporation registered in the U.S. and other countries. SYNCT is a trademark of Luminex Corporation.

www.luminexcorp.com

HEADQUARTERS
UNITED STATES
+1.512.219.8020
info@luminexcorp.com

EUROPE
+31.73.800.1900
europe@luminexcorp.com

CANADA
+1.416.593.4323
info@luminexcorp.com

CHINA
+86.21.8036.9888
info@luminexcorp.com

JAPAN
+81.3.5545.7440
info@luminexcorp.com

FL19258