

SYNCT™

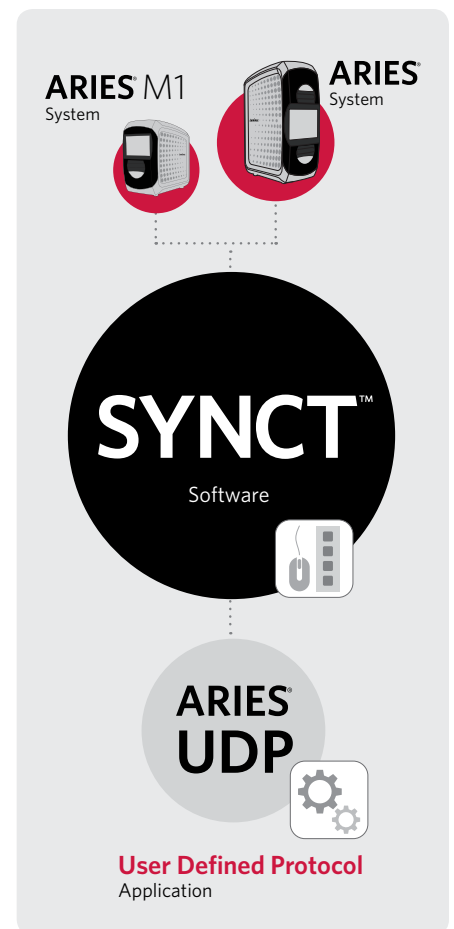
SYNCT™ Software | The Power Behind Your Lab

SYNCT™ ARIES® UDP App

The ARIES® User Defined Protocol (UDP) Application enables laboratories to create user defined protocols for ARIES® Systems. The application is an optional utility that may be installed on SYNCT™ Software.

Create User Defined Assay Protocol Files in Four Simple Steps

- 1. Define Targets:** Define targets on any of the 1-6 channels, including Sample Processing Controls (SPC).
- 2. Set Thermal Cycling:** Set thermal cycling conditions easily with drag and drop functionality, or leave the default Luminex universal cycling conditions to run laboratory developed tests (LDTs) in the same batch as ARIES® IVD assays. (See Figure 1)
- 3. Gather Data and Set Call Parameters:** Run your LDT to gather data needed to graphically set the Ct and Tm windows used in making call determinations.
- 4. Save and Run:** Save the new assay protocol file, and then export the file to an ARIES® instrument to begin running LDTs.



SYNCT™ Software provides access to all ARIES® data in one place, and offers novel applications to expand functionality.

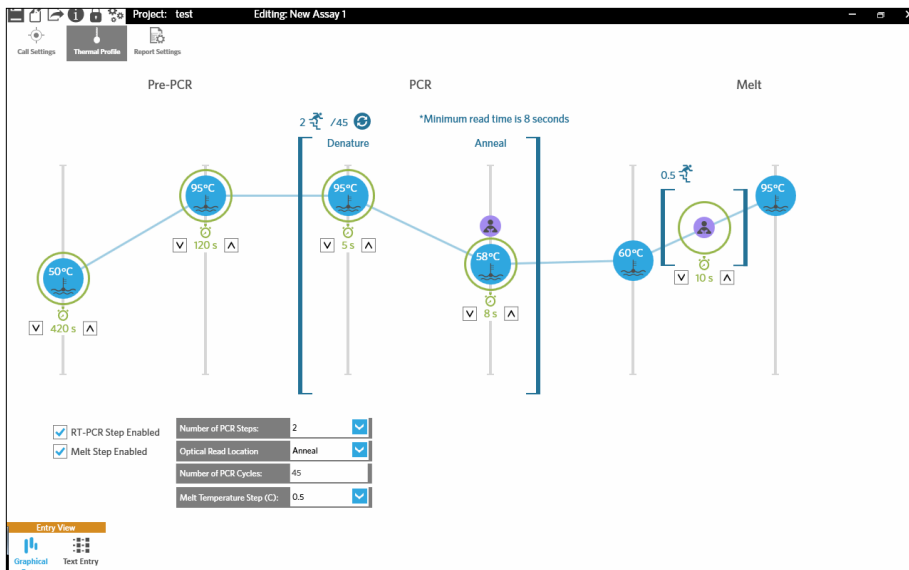


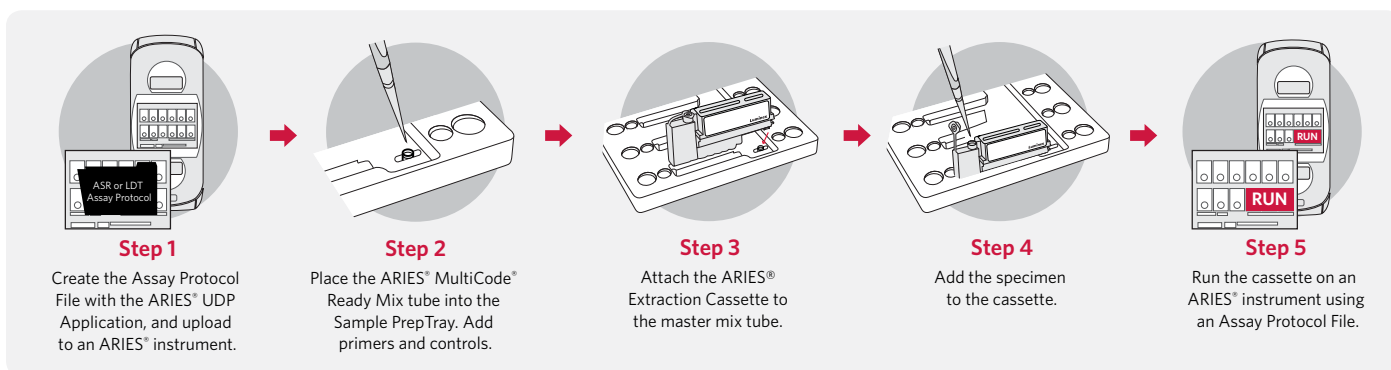
Figure 1: Image from SYNCT™ thermal cycling condition screen.

Learn More About Our UDP Tools

In addition to SYNCT™ Software tools, we offer other items to support your testing needs.

For more information, please contact your local Luminex representative.

Workflow



Ordering Information

Product Name	Part Number
ARIES® UDP App	CN-SW52-01
SYNCT™ (IVD)*	CN-SW47

*SYNCT™ is required to run the ARIES® UDP (User Defined Protocol) App.

Luminex
complexity simplified.

orders@luminexcorp.com or support@luminexcorp.com

ARIES® Systems are cleared by the FDA only when used with ARIES IVD assays which have been cleared by the FDA. Products are region specific and may not be approved in some countries/regions. Please contact Luminex at support@luminexcorp.com to obtain the appropriate product information for your country of residence. ARIES Systems are class 1(I) laser products.

©2016-2017 Luminex Corporation. All rights reserved. Luminex and ARIES 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

+86.21.802.31150

info@cn.luminexcorp.com

JAPAN

+81.3.5545.7440

info@jp.luminexcorp.com

5540846