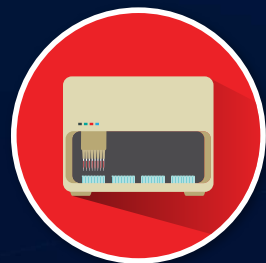
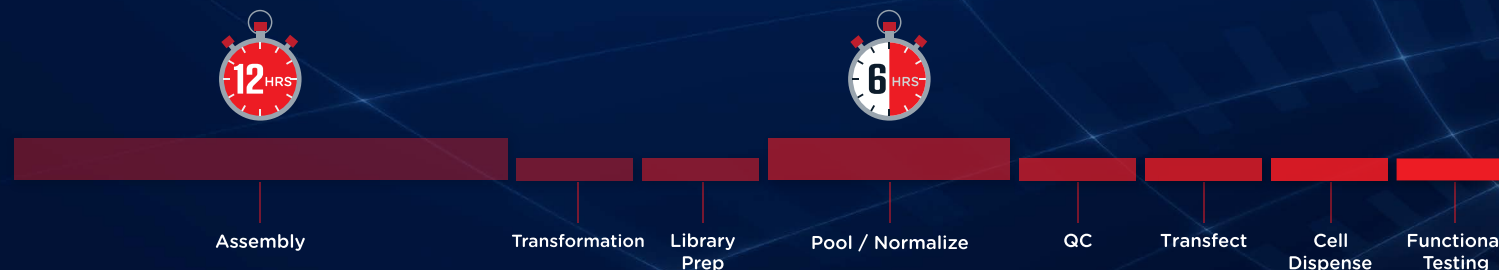


Save 15 Hours on Your Synthetic Biology Workflow

LABCYTE 
The Future of Science is Sound



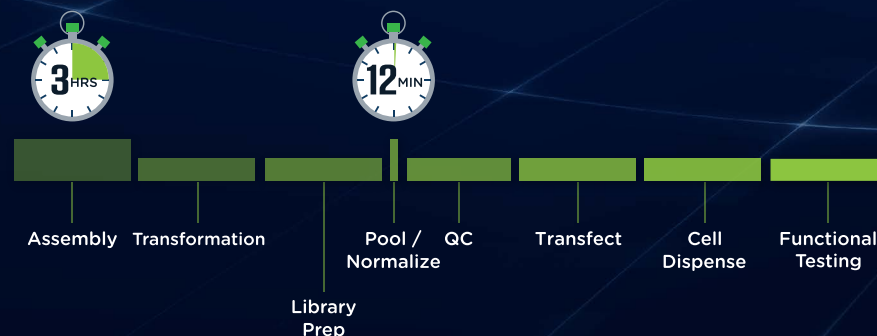
Traditional
LIQUID HANDLER



Advantages of Echo[®] Acoustic Liquid Handling



Echo[®] Acoustic
LIQUID HANDLER




15 HRS
faster



 Echo[®] Acoustic LIQUID HANDLING
for SYNTHETIC BIOLOGY

www.labcyte.com/synbio

info-us@labcyte.com

Follow Us on Twitter:  @LabcyteInc

FOR RESEARCH USE ONLY. Not for use in diagnostic procedures.

© 2017 LABCYTE INC. All rights reserved. Labcyte[®] and Echo[®] are registered trademarks of Labcyte Inc., in the U.S. and/or other countries.