

Traditional Gel Clot Limulus Amebocyte Lysate (LAL) Assay Procedure Quick Guide

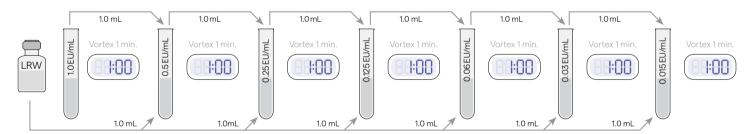
This is a step by step guide depicting how to perform a traditional gel clot LAL assay. Prior to initiating the assay procedure, allow reagent vials to equilibrate to room temperature.

Step 1 Step 2 Step 3 Reconstitute Control Vortex for 15 minutes. Dilute endotoxin with LRW to a Standard Endotoxin (CSE) concentration of 1.0 EU/mL using the 15:88 with LAL Reagent Water endotoxin potency identified on the (LRW). EU/mL CSE Certificate of Analysis (CoA). 0. LRW CSE CSE ĹRW

Step 4

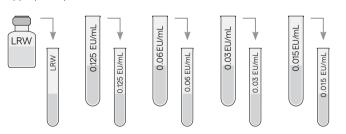
Label the tubes with the appropriate endotoxin concentration and add 1.0 mL of LRW to each. Using the 1.0 EU/mL solution, prepare serial 2-fold dilutions to bracket lysate sensitivity.

(Example based on a test using 0.06 EU/mL lysate sensitivity.)



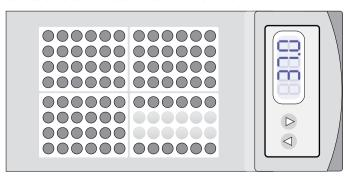
Step 5

Add 100 μ L of endotoxin standards and negative control (LRW) into each appropriately labeled 10 × 75 mm reaction tube.



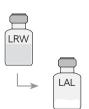
Stan 8

Place tubes in 37°C ± 1°C non-circulating water bath or heating block for 60 minutes. Note: Water or dry bath should be located in the lab away from all sources of vibration that could disturb clot formation.



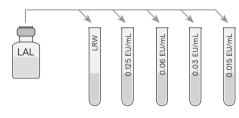
Step 6

Immediately prior to use, reconstitute LAL and gently swirl.



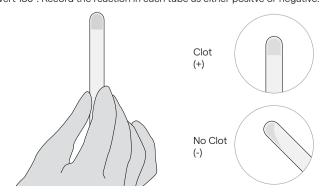
Step 7

Add 100 µL reconstituted LAL into each reaction tube.



Step 9

After 60 minutes (± 2 minutes) of incubation, carefully remove each tube and invert 180°. Record the reaction in each tube as either positive or negative.



Materials, equipment and documents needed

Reagents

- Limulus Amebocyte Lysate (LAL) Reagent (PYROGENT® Reagent)
- Control Standard Endotoxin (CSE)
- LAL Reagent Water (LRW) (# W50-640, W50-100, W50-500)

Kits are available in a wide range of sizes and configurations.

Please contact your local sales representative for additional information.

Accessories

- Glass dilution tubes, 13 × 100 mm (# N207)
- Glass reaction tubes, 10 × 75 mm (# N201, N205)
- Individually wrapped serological pipettes (optional)
- Tips

Use pyrogen-free accessories that have been qualified for endotoxin testing.

Equipment and software

- Heating block or non-circulating hot water bath
- Pipettors
- Timer
- Vortex mixer
- Test tube rack

Supporting documents

- Certificate of Analysis (CoA), www.lonza.com/coa
- Limulus Amebocyte Lysate (LAL) PYROGENT® and PYROGENT® Plus Package Insert

▲ Points to consider

- Use matched LAL and CSE reagents
- Plastic tubes are not recommended for making endotoxin dilutions
- Follow all suggested endotoxin vortexing times
- Use pyrogen-free accessories that have been qualified for endotoxin testing
- Equilibrate reagents to room temperature before use
- Do not vortex the LAL
- Avoid contaminating samples, standards, LRW and accessories
- Equipment should be installed, validated and maintained appropriately
- Avoid disturbing the clot

Contact us

North America

Customer Service: +1800 638 8174 (toll free)

order.us@lonza.com

Scientific Support: +1800 521 0390 (toll free)

scientific.support@lonza.com

Europe

Customer Service: + 32 87 321 611

order.europe@lonza.com

Scientific Support: + 32 87 321 611 scientific.support.eu@lonza.com

International

Contact your local Lonza Distributor
Customer Service: +13018987025
Fax: +13018458291
scientific.support@lonza.com

Lonza Walkersville, Inc. - Walkersville, MD 21793

All trademarks belong to Lonza, registered in USA, EU or CH or to third party owners and used only for informational purposes. The information contained herein is believed to be correct and corresponds to the latest state of scientific and technical knowledge. However, no warranty is made, either expressed or implied, regarding its accuracy or the results to be obtained from the use of such information and no warranty is expressed or implied concerning the use of these products. The buyer assumes all risks of use and/or handling. Any user must make his own determination and satisfy himself that the products supplied by Lonza Group Ltd or its affiliates and the information and recommendations given by Lonza Group Ltd or its affiliates are (i) suitable for intended process or purpose, (ii) in compliance with environmental, health and safety regulations, and (iii) will not infringe any third party's intellectual property rights. The user bears the sole responsibility for determining the existence of any such third party rights, as well as obtaining any necessary licenses. For more details: www.lonza.com/legal. ©2021 Lonza. All rights reserved. RT-DS014 01/22

www.lonza.com/gelclot