

## ReadiPrep™ Lipopolysaccharide (LPS) Isolation Kit

Catalog number: 60201  
Unit size: 10 Preps

Component	Storage	Amount (Cat No. 60201)
Component A: LPS Isolation Buffer	Refrigerated (2-8 °C), Minimize light exposure	1 Bottle (100 mL)
Component B: Proteinase K (20 mg/mL)	Freeze (< -15 °C)	1 Vial (0.6 mL)

### OVERVIEW

The ReadiPrep™ Lipopolysaccharide (LPS) Isolation Kit provides a safe and effective way to isolate LPS from the outer membrane of Gram-negative bacteria. It utilizes a bacterial membrane lysis buffer and enzymatic protein digestion to extract pure LPS from bacterial cultures, which can be quantified using carbohydrate detection methods. Unlike conventional approaches that rely on hazardous organic chemicals like chloroform or phenol, this kit provides a safer and more time-efficient alternative for isolating high-purity LPS for research applications. Lipopolysaccharides, or LPS, are found in the outer membrane of Gram-negative bacteria. It is a carbohydrate with a low molecular weight of 10-20 kDa. It is a complex molecule made of heterogeneous components comprising O antigen, core oligosaccharide, Lipid A, and non-carbohydrate components such as phosphate and amino acids groups. Lipid A, characterized by its multiple fatty acid chains, anchors the LPS into the bacterial membrane and contributes to the toxicity of the Gram-negative bacteria. Known as endotoxins, LPS is notorious for inducing potent inflammatory responses and sepsis when consumed by animals.

### AT A GLANCE

#### Protocol Summary

1. Grow bacteria on the LB agar plate.
2. Add LPS isolation buffer to the bacterial pellet. Sonicate the mixture, then incubate it on ice for 10 minutes.
3. Add proteinase K to the bacterial lysate and incubate the mixture at 60°C for one hour.

#### Important Note

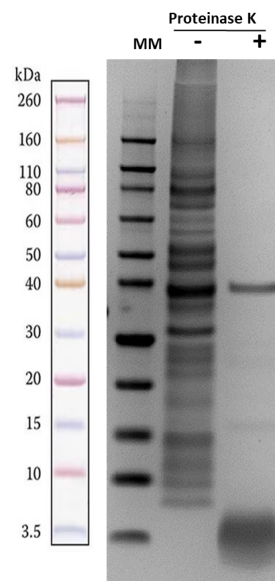
Thaw all the kit components at room temperature before starting the experiment.

### SAMPLE EXPERIMENTAL PROTOCOL

1. Grow the bacteria on an LB agar plate overnight at 37°C.
2. The next day, collect bacterial colonies in ice-cold PBS. Measure the optical density (OD) at 600 nm using a spectrophotometer. Ensure the OD at 600 nm is greater than 0.6.
3. Centrifuge the tube at 2500g for 10 minutes to collect the pellet. After removing the supernatant, centrifuge again at 2500g for 10 minutes to ensure all the supernatant is removed.
4. Measured the weight of the pellets.  
  
**Note:** For an E. coli suspension with OD > 0.6, the pellet weight should be greater than 10 mg.
5. Add 200 µL of isolation buffer to the LPS pellet. Use 200 µL for a 20 mg pellet.
6. Sonicate the pellet three times for 30 seconds each at 10 watts. Then, incubate the tube on ice for 10 minutes.

7. Centrifuge the tube at 2500g for 10 minutes at 4°C.
8. Transfer the lysate into a new, clean tube.
9. Add Proteinase K to the lysate at a concentration of 0.1 mg/mL.  
  
**Note:** For example, if you have 200 µL of lysate, add 1 µL of 20 mg/mL Proteinase K solution.
10. Incubate the tube at 60 °C for one hour.
11. Centrifuge the tube at 2500g for 10 minutes at 4°C.
12. Collect the supernatant.
13. LPS can be measured using carbohydrate detection methods. Alternatively, its purity can be assessed by running it on an SDS-PAGE gel and then staining it with Coomassie blue.

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Bacterial lysate, prepared using the ReadiPrep™ Lipopolysaccharide (LPS) Isolation Kit, was boiled at 85 °C for 3-5 minutes in 1X loading buffer. The sample was then loaded onto a 4-12% SDS-PAGE gel and run for 70 minutes at 110V. Afterward, the gel was stained with Coomassie Blue to visualize the proteins.



**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.