

# One Shot™ LB Agar Plates with 100 µg/mL Ampicillin

Catalog Number A55802

Pub. No. MAN0029965 Rev. A.0



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

## Product description

One Shot™ LB Agar Plates are pre-poured plates containing Luria-Bertani (LB) agar medium with 100 µg/mL ampicillin. LB agar is a solid bacterial growth medium used in molecular biology studies for the cultivation and maintenance of *Escherichia coli* strains. Added antibiotic allows selection of ampicillin resistant bacteria.

- Composition of the medium is based on formulation developed by Jeffrey H. Miller and contains twice as much sodium chloride as in the Lennox formulation [1].
- LB agar is a nutrient-rich medium containing casein peptone, yeast extract, sodium chloride, and agar. Casein peptone or tryptone is an enzymatic hydrolysate of casein. It supplies nitrogen, carbon, and amino acids. Yeast extract provides trace elements, vitamins (including B group vitamins), amino acids, and peptides. Sodium chloride supplies sodium cations for the membrane transport and maintains osmotic balance. Agar is a solidifying agent of the growth medium.
- Autoclaved glass plating beads are provided for convenient spreading of bacterial cells across the surface of the solid LB agar medium.

## Contents

| Catalog number | Amount    |
|----------------|-----------|
| A55802         | 20 plates |

The product is shipped as a box containing 20 pre-poured LB agar plates with antibiotic. Plates are divided into two stacks wrapped in a single layer of cellophane, containing 10 plates per stack. Each plate is 90 mm in diameter, contains 21 ± 2 mL of LB agar medium.

A 15 mL conical tube with autoclaved 4 mm glass plating beads is included in the box.

## Storage and stability

Store at 2–14°C in the original packaging (cellophane wrapping and box). Protect from the light. Avoid freezing and overheating.

**Note:** Always store agar plates upside down.

Pre-poured LB agar plates are stable for three months when stored as directed.

## Composition

| Component       | Concentration, g/L |
|-----------------|--------------------|
| Casein peptone  | 10.00              |
| Yeast extract   | 5.00               |
| Sodium chloride | 10.00              |
| Ampicillin      | 0.1                |
| Agar            | 15.00              |

## Procedure

### Innoculate culture plate

Follow procedures established in the laboratory for bacterial culture preparation, inoculation and incubation. For more information, refer to relevant sources.

- Prior to inoculation, it is recommended to prewarm agar plates at 37°C in the incubator for 30 minutes. This step will help to dry out excessive moisture from the surface of the medium and provides better recovery of stressed bacterial cells.
- In case of built-up condensation, dry plates in the laminar flow hood with half-opened lids for 10–30 minutes.

### How to use glass plating beads

Pipet bacterial cells onto the agar plate, then add 5–7 beads and vigorously move the plate back and forth. Make a couple of 360° rotations with 90° stops, until the surface of the medium seems dry. Discard the beads into the waste container by flipping the plate upside down.

## Quality control

### Physical and chemical quality control

- Color: Straw yellow
- Final pH: 7.0 ± 0.2 at 25°C

### Microbiological control

- Inoculum level for productivity testing: a practicable range of 100±20 CFU of the target microorganism per plate with minimum number of 50 CFU per plate.
- Inoculum level for selectivity testing: a range of 10<sup>4</sup>–10<sup>6</sup> CFU of the non-target microorganism per plate.
- Aerobic incubation at 37±1°C. Reading, calculation and interpretation of results after 18–24 hours.

| Microorganism                                   | Result                          |
|---|---------------------------------|
| <i>Escherichia coli</i> ATCC® 25922, WDCM 00013 | Inhibited                       |
| <i>Escherichia coli</i> ATCC® 8739, WDCM 00012  | Inhibited                       |
| <i>Escherichia coli</i> DH5α containing pUC19   | Good growth (productivity ≥50%) |

### Test for absence of microbial contamination on plates

Testing is performed by incubation at 20–25°C and 30–35°C for 48 hours followed by visual inspection. Subsequent incubation under the same conditions for another seven days with following inspection.

Result: No growth.

## References

[1] Miller, J.H. (1972) *Experiments in Molecular Genetics*. Cold Spring Harbor Laboratory, Cold Spring Harbor, New York.

## Related products

The following products are available through [thermofisher.com](https://www.thermofisher.com).

Catalog numbers that appear as links open the web pages for those products.

| Product  | Quantity        | Catalog no.               |
|--|-----------------|---------------------------|
| TOPO™ TA Cloning™ Kit for Subcloning, without competent cells  | 25 reactions    | <a href="#">450641</a>    |
| TOPO™ TA Cloning™ Kit for Sequencing, without competent cells  | 25 reactions    | <a href="#">450030</a>    |
| Zero Blunt™ PCR Cloning Kit, without competent cells   | 25 reactions    | <a href="#">450245</a>    |
| Zero Blunt™ TOPO™ PCR Cloning Kit for Sequencing, without competent cells                                | 25 reactions    | <a href="#">450031</a>    |
| Gateway™ LR Clonase™ II Enzyme Mix   | 20 reactions    | <a href="#">11791020</a>  |
| Gateway™ BP Clonase™ Enzyme Mix  | 20 reactions    | <a href="#">11789013</a>  |
| pENTR™/D-TOPO™ Cloning Kit, with One Shot™ Mach1™ T1 Phage-Resistant Chemically Competent <i>E. coli</i> | 20 reactions    | <a href="#">K240020</a>   |
| pENTR™/D-TOPO™ Cloning Kit   | 20 reactions    | <a href="#">K240020SP</a> |
| GeneArt™ Gibson Assembly® HiFi Master Mix  | 50 reactions    | <a href="#">A46628</a>    |
| GeneArt™ Gibson Assembly® EX Master Mix  | 50 reactions    | <a href="#">A46636</a>    |
| One Shot™ TOP10 Chemically Competent <i>E. coli</i>  | 21 × 50 µL/tube | <a href="#">C404003</a>   |
| One Shot™ Stbl3™ Chemically Competent <i>E. coli</i>   | 21 × 50 µL/tube | <a href="#">C737303</a>   |
| MegaX DH10B™ T1 <sup>R</sup> Electrocomp™ Cells  | 5 × 100 µL      | <a href="#">C640003</a>   |
| ElectroMAX™ Stbl4™ Competent Cells   | 5 × 100 µL      | <a href="#">11635018</a>  |
| Ampicillin, sodium salt, irradiated  | 200 mg          | <a href="#">11593027</a>  |
| Kanamycin Sulfate  | 5 g             | <a href="#">11815024</a>  |
| X-Gal  | 1 g             | <a href="#">15520018</a>  |
| IPTG   | 1 g             | <a href="#">15529019</a>  |
| Bluo-Gal, substrate for β-galactosidase  | 1 g             | <a href="#">15519028</a>  |
| S.O.C. Medium  | 10 × 10 mL      | <a href="#">15544034</a>  |
| LB Broth   | 500 mL          | <a href="#">10855021</a>  |
| LB Agar, powder (Lennox L agar)  | 500 g           | <a href="#">22700025</a>  |
| LB Broth Base  | 500 g           | <a href="#">12780052</a>  |
| One Shot™ LB Agar Plates   | 20 plates       | <a href="#">A55800</a>    |
| One Shot™ LB Agar Plates with 50 µg/mL Kanamycin   | 20 plates       | <a href="#">A55803</a>    |

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](https://www.thermofisher.com/support).



Revision history: Pub. No. MAN0029965 A.0

| Revision | Date         | Description  |
|----------|--------------|--|
| A.0      | 7 March 2024 | Initial product insert release for One Shot™ LB Agar Plates with Ampicillin. |

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.