

# Preparation of X-Gal/IPTG LB Agar Plates for Blue/White Colony Screening

This protocol is for the Preparation of X-Gal/IPTG LB Agar Plates for Blue/White Colony Screening.

For individual LB (Luria Broth) agar plates:

1. Pour sterile warm LB agar (about 25 mL) into a Petri dish.
2. Dry opened LB plates at room temperature under UV light for about 30 minutes.
3. Add 40  $\mu$ L of the Thermo Scientific X-Gal Solution (20 mg/mL), ready-to-use (Cat #R0941).
4. Add 40  $\mu$ L of 100 mM Thermo Scientific IPTG Solution, ready-to-use (Cat #R1171)
5. Spread evenly on the plate with a sterile spatula.

For batch use, add the following directly per 1 mL of the liquid LB agar (kept at about 50 °C):

6. 1  $\mu$ L of X-Gal Solution (20 mg/mL), ready-to-use.
7. 1  $\mu$ L of 100 mM IPTG Solution, ready-to-use.
8. Mix well.
9. Pour 25 mL of prepared LB agar into each Petri dish.
10. Dry opened LB plates at room temperature under UV light for about 30 minutes.

[thermoscientific.com/onebio](http://thermoscientific.com/onebio)

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

#### Europe

Customer Service  
cs.molbio.eu@thermofisher.com

Technical Support  
ts.molbio.eu@thermofisher.com

Tel 00800 222 00 888  
Fax 00800 222 00 889

#### United States

Customer Service  
cs.molbio@thermofisher.com

Technical Support  
ts.molbio@thermofisher.com

Tel 800 235 9880  
Fax 800 292 6088

#### Canada

Customer Service  
cs.molbio@thermofisher.com

Technical Support  
ts.molbio@thermofisher.com

Tel 800 340 9026  
Fax 800 472 8322

**Thermo**  
S C I E N T I F I C

Part of Thermo Fisher Scientific