



# LB Agar (Premade)

(Cat. # L011)

## ITEMS INCLUDED

Cat. #	Description	Size
L011	LB Agar (Premade)	100ml

## STORAGE CONDITION

Shipped at ambient temperature. Upon receiving, store at room temperature

## PROTOCOL

1. Use microwave to dissolve the LB Agar (Premade) media completely.
2. Once the LB Agar has cooled to hand hot temperature (about 45°C), add appropriate antibiotics and pour a ~0.5cm/ ¼" layer of agar into Petri dishes. This requires approximately 20-25ml volume of LB agar per Petri dish.
3. Leave the Petri dish on a flat surface for 30-40 minutes until completely solidified.

## Approximate Formula/ Liter

Casein Digest Peptone (Tryptone): 10g

Yeast Extract: 5g

Sodium Chloride: 5g

Agar: 15g

Final pH: 7.0 ± 0.2 at 25°C

