

# LM Agarose, L pack

Low Melting (LM) Agaroses are derivatized by organic synthesis which generates methoxylate groups from the basic agarose structure. The main properties of these agaroses are their low melting and gelling temperatures when compared with standard agaroses. LM Agarose is a recommended agarose for electrophoresis of DNA fragments  $\geq 10000$  bp, tissue/cell culture and viral plaque assays. LM Agarose is characterized by lower gel strength and highest clarity compared to standard agaroses. The low melting temperature allows the recovery of undamaged nucleic acids below the denaturation temperature. The low gelling temperature ensures that the agarose will be in a liquid state at a temperature range where in-Gel manipulations can be performed without prior extraction of the DNA from the gel slice.

## Ordering Information

[Order Online »](#)

JBS-PCR-277L	5x50g
--------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      Low melting Agarose

Appearance      white fine homogeneous powder

**Technical Info / Product Notes**      For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)