



LM Agarose

Low melting Agarose

Cat. No.	Amount
PCR-277S	50 g
PCR-277L	5 x 50 g

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 12 months

Form: fine homogeneous powder

Color: white

Description:

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 ≥ 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.

Specifications:

Moisture	$\leq 10 \%$
Ash	$\leq 0.4 \%$
Electroendosmosis	≤ 0.12
Sulphate	$\leq 0.12 \%$
Clarity 1.5 % (NTU)	≤ 4
Gel Strength 1.5 %	$\geq 500 \text{ g/cm}^2$
Gelling Temperature 1.5 %	24 - 28 °C
Melting Temperature 1.5 %	$\leq 65 \text{ °C}$

Functional tests

DNase/RNase activity	none detected
DNA Binding	very low
DNA Resolution	finally resolved
Gel Background	very low

Recommended Agarose Concentrations:

Size Range (base pairs)	Final Agarose Concentration [%] 1x TAE Buffer
5000 - 20000	0.60
300 - 15000	1.00
250 - 10000	1.25
200 - 5000	1.50
100 - 2000	1.75
80 - 1000	2.00