



LE Agarose

general use Agarose

Cat. No.	Amount
PCR-269S	2 x 100 g
PCR-269L	2 x 500 g

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 24 months

Form: fine homogeneous powder

Color: white

Applications:

- Nucleic acid analytical and preparative electrophoresis
- Blotting
- Protein electrophoresis such as radial immunodiffusion

Description:

LE Agarose is a general use agarose for routine nucleic acid analytical and preparative electrophoresis of fragments between 200 - 20000 bp, blotting and protein electrophoresis. This agarose has a low electroendosmosis value (EEO), ensuring high electrophoretic mobility of nucleic acids. Because of the high gel strength, easy-to-handle gels can be prepared with this agarose at concentrations as low as 0.4 %. LE Agarose has an exceptionally low fluorescence background.

Specifications:

Moisture	≤ 10 %
Ash	≤ 0.4 %
Electroendosmosis	0.05 - 0.13
Sulphate	≤ 0.1 %
Clarity 1.5 % (NTU)	≤ 3
Gel Strength 1 %	≥ 1200 g/cm ²
Gel Strength 1.5 %	≥ 2500 g/cm ²
Gelling Temperature 1.5 %	36 ± 1.5 °C
Melting Temperature 1.5 %	88 ± 1.5 °C

Functional tests

DNase/RNase activity	none detected
DNA resolution > 1000 bp	finally resolved
Gel background	very low

Recommended Agarose Concentrations:

Size Range (base pairs)	Final Agarose Concentration [%] 1x TAE Buffer
1000 - 20000	0.60
500 - 12000	0.80
300 - 8000	1.00
200 - 6000	1.20
100 - 3000	1.50
100 - 2000	2.00