



Product Information

RNA MARKER 0.28-6.6 KB

Product No. R 7644

TECHNICAL BULLETIN

Product Description

Suitable for use as a molecular weight marker for formaldehyde agarose gel electrophoresis.

Usage: 3 μ l per lane. See suitability assay for details. Sufficient for 25 uses.

Storage: -70 °C

Storage Buffer: 10 mM Tris-HCl, pH 8.0, 1 mM EDTA

10X MOPS Electrophoresis buffer: 400 mM MOPS, 100 mM sodium acetate, 10 mM EDTA

SUITABILITY ASSAY

RNA Marker sample solutions were prepared for electrophoresis as follows:

- 3 μ l RNA Marker
- 2 μ l RNA Sample Loading Buffer (Product No. R 4268) 62.5% Deionized Formamide, 1.14 M Formaldehyde, 1.25X MOPS-EDTA-Sodium Acetate Buffer (Product No. M 5755, diluted 1:8) 200 μ g/ml Bromphenol Blue 200 μ g/ml Xylene Cyanole 50 μ g/ml Ethidium Bromide.

The RNA marker sample solution was incubated at 65 °C for 5 minutes and immediately cooled on ice.

The entire 5 μ l of RNA Marker solution was run with appropriate RNA markers on a 10 mm thick denaturing (formaldehyde) agarose gel. Electrophoresis was performed in a mini submarine-type apparatus at 100 V for 2 hours in 1x MOPS electrophoresis buffer with buffer recirculation. The gel was stained in 5 μ g/ml ethidium bromide for 15 minutes and destained 1 hour with shaking in water. Nine bands were resolved and the band pattern was consistent with the sizes listed below.

FRAGMENT SIZES (bases)

6583	1383
4981	955
3638	623
2604	281
1908	

References

1. Sambrook, J., and Russell, D. W., *Molecular Cloning, A Laboratory Manual*, (Cold Spring Harbor Laboratory 2001) p. 7.31-7.34
2. Fasman, G.D., ed., *Practical Handbook of Biochemistry and Molecular Biology*, (CRC Press, 1986) p. 464.

JLL/JWM 2/04

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.