



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

EMBER™ Ultra RNA Gel Kit

The most sensitive RNA gel stain available in convenient pre-coated agarose format. Includes formamide-containing RNA loading dye.



Product attributes

| | |
|--------------------|--|
| Storage Conditions | Components have different storage temperatures, open kit on receipt. |
| Assay type/options | RNA gel staining |

Product Description

The EMBER™ Ultra RNA Gel system provides the ultimate in sensitivity and resolution for RNA gel electrophoresis in a fast and simple agarose gel format. Along with being much more sensitive than any other RNA gel stains, EMBER™ Ultra also does not require a post-electrophoresis staining step. Simply cast your agarose gel using the EMBER™ Ultra Precoated Agarose, and prepare your samples in the denaturing EMBER™ Ultra RNA Loading Dye. Bands can be imaged immediately after electrophoresis.

Note: The number of gels per kit is based on a 9×11 cm, 1% agarose mini gel (0.5 g agarose in 50 mL buffer per gel). The actual number may vary depending on gel size.

Simple, Quick, & Highly Sensitive RNA Gel Staining

EMBER™ Ultra Precoated Agarose is ultra-pure molecular biology grade agarose with low electroendosmosis that is precoated with Biotium's fluorescent EMBER™ Ultra Dye. This novel gel system using EMBER™ Ultra Precoated Agarose with EMBER™ Ultra Loading Dye is more sensitive than currently available RNA dyes, allowing detection of as little as ≤5 ng RNA. The gel system does not require a post-electrophoresis staining step or a special gel system, for convenient and fast results. EMBER™ Ultra Precoated Agarose gives excellent results at percentages between 1% and 5% agarose, for analyzing different sizes of RNA.

EMBER™ Ultra Dye has green fluorescence with excitation/emission at 500/530 nm when bound to nucleic acids. For best results, we recommend imaging with a Biotium's [Gel-Bright™ Laser Diode Gel Illuminator](#) or a blue light illuminator. Gels can also be imaged using a UV transilluminator equipped with a SYBR® filter (recommended) or EtBr filter (with lower signal). The EMBER™ Ultra RNA Loading Dye is a formamide-containing denaturing RNA loading dye specifically formulated to work with EMBER™ Ultra Precoated Agarose to provide high sensitivity fluorescent staining of RNA in a fast and convenient agarose gel. The loading dye contains blue and orange tracking dyes that migrate at ~1500 bp (blue) and ~50 bp (orange) in a 1% agarose gel.

Note: DNA will also stain on EMBER™ Ultra RNA Gels, so any contaminating DNA in your RNA sample will also be detected. For best results with DNA, we recommend our [EMBER™ Ultra DNA Agarose Gel Kit](#).

[Learn more about Biotium's selection of unique solutions for RNA workflows, including reagents for RNA extraction, quantitation, and RNase decontamination.](#)

SYBR is a registered trademark of Thermo Fisher Scientific.

This datasheet was generated on January 12, 2026 at 09:20:16 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ember-ultra-rna-gel-kit/>