Glowing products for science

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

Call us: 800-304-5357 Email: btinfo@biotium.com

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.

EMBERTM 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-BrightTM 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 EMBERTM Ultra RNA Loading Dye is a formamide-containing denaturing RNA loading dye specifically formulated to work with EMBERTM 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 EMBERTM Ultra RNA Gels, so any contaminating DNA in your RNA sample will also be detected. For best results with DNA, we recommend our EMBERTM Ultra DNA Agarose GelKit.

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 November 17, 2025 at 10:28:31 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ember-ultra-rna-gel-kit/