Thiazole Green (SYBR® Green I), 10,000X in DMSO

Thiazole Green, which is structurally identical to SYBR® Green I Nucleic Acid Gel Stain, is one of the most sensitive stains available for detecting double-stranded DNA (dsDNA) in agarose, polyacrylamide gels and qPCR.

Blotum

Product attributes

| Probe cellular localization | Nucleus & cytoplasm |
|-----------------------------|-------------------------------------|
| For live or fixed cells | For live/intact cells |
| Assay type/options | No-wash staining, Real-time imaging |
| Cell permeability | Membrane permeant |
| Colors | Green |
| Excitation/Emission | 498/522 pm (with DNA) |

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

Product Description

Thiazole Green, is structurally identical to the well known qPCR and DNA gel stain SYBR® Green I. Thiazole Green is one of the most sensitive stains available for detecting double-stranded DNA (dsDNA) in agarose, polyacrylamide gels and qPCR. Thiazole Green can also be used to detect ssDNA and RNA in denaturing agarose/formaldehyde and polyacrylamide/urea gels without any prewashing steps, although the sensitivity is lower. Thiazole Green may also be used as a green nuclear stain for all cells in live cultures, but will lose nuclear specificity after fixation. Thiazole Green has the identical spectral properties as SYBR® Green I (Ex/Em 498/522 nm with DNA).

SYBR is a registered trademark of Thermo Fisher Scientific.

This datasheet was generated on November 11, 2025 at 07:18:42 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/thiazole-green-10000x-in-dmso/