

## 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.

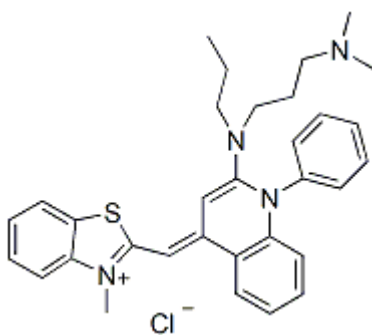


### Product attributes

|                             |                                     |
|-----------------------------|-------------------------------------|
| 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 nm (with DNA)               |

### 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 June 6, 2026 at 02:55:32 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/thiazole-green-10000x-in-dmso/>