



Glowing products for science

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.



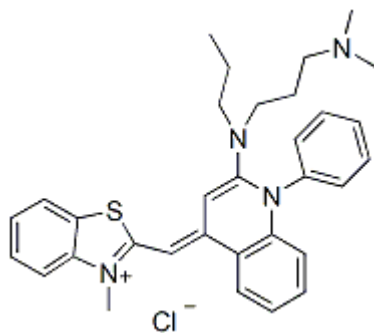
Call us : [800-304-5357](tel:800-304-5357)

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 January 14, 2026 at 10:17:13 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/thiazole-green-10000x-in-dmsol/>