46117 Landing Parkway, Fremont, CA U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Thiazole Green, 10,000X in DMSO

CATALOG NUMBER: 40086-0.5mL/40086-1mL

MOLECULAR $C_{32}H_{37}ClN_4S$ **INFORMATION:** Mwt: 545.19

CAS#: 163795-75-3

PROPERTIES:

Color & Form Orange solution
Purity > 95% by HPLC

 $\lambda_{\text{ex}} \setminus \lambda_{\text{em}}$ 497/520nm (with DNA)

STORAGE AND HANDLING:

Store at -20 °C and protect from light.

APPLICATION: 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. 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 has the identical spectral properties as SYBR® Green I

TOXICITY: Unknown

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on

clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.