

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: TMRE, 2mM in DMSO

CATALOG NUMBER: 70005

MOLECULAR $C_{26}H_{27}ClIN_2O_7$ **INFORMATION:** Mwt: 515

PROPERTIES:

Color & FormRed solutionPurity $\geq 95\%$ by HPLC $\lambda_{ex} \mid \lambda_{em}$ 549/574nm

STORAGE AND HANDLING:

Store at 4 °C and protect from light.

APPLICATION: TMRE (Tetramethylrhodamine ethyl ester, perchlorate) and TMRM (70017) are

preferred dyes for quantitative measurements of membrane potentials using the Nernst equation. TMRE has been used to separate viable and non-apoptotic cells from cell sorting samples (1). TMRE staining is reversible and does not affect cell proliferation

and viability. TMRE is also available in solid form (70016).

Ref.: 1) J Histochem Cytochem doi: 10.1369/0022155413520404

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.

PSF006 Page 1 of 1