TMRE, 2 mM in DMSO

TMRE (Tetramethylrhodamine ethyl ester, perchlorate) is used to quantify changes in mitochondrial membrane potential in live cells.



Product Description

TMRE (Tetramethylrhodamine ethyl ester, perchlorate) is a cell permeable, cationic, red-orange fluorescent dye that accumulates active mitochondria. It is used to detect changes in mitochondrial membrane potential of live cells via microplate fluorometry, fluorescence microscopy and flow cytometry. Inactive or depolarized mitochondria have decreased membrane potential and thus fail to retain the TMRE dye.

- Red-orange fluorescence, $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 549/574 nm
- Suitable for staining mitochondria in live cells only. It does not work with fixed cells
- TMRE is used for quantitative measurement of membrane potential using the Nernst equation
- Store the solution at 4°C and protect from light
- C₂₆H₂₇CIIN₂O₇
- MWt: 515

TMRE and TMRM (70017) are preferred dyes for quantitative measurements of membrane potentials using the Nernst equation. TMRE can be 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).

References

1. J Histochem Cytochem doi: 10.1369/0022155413520404

This datasheet was generated on November 15, 2025 at 02:36:40 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/tmre-2-mm-dmso/

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

Product attributes

Probe cellular localization	Mitochondria
CAS number	115532-52-0
For live or fixed cells	For live/intact cells
Assay type/options	No-wash staining, Real-time imaging
Cell permeability	Membrane permeant
Apoptosis/viability marker	Mitochondrial potential
Potential dependence	Mitochondrial potential-dependent
Colors	Red
Excitation/Emission	549/574 nm