Mag-Fluo-4, Tetrapotassium Salt

Membrane-impermeant form of Mag-Fluo-4, a fluorescent magnesium and calcium ion indicator that is an analog of Fluo-4.



Product attributes

Cell permeability	Membrane impermeant
Indicator type	Non-ratiometric
Excitation/Emission	493/517 nm (high Mg2+)

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

Product Description

Mag-Fluo-4 is a fluorescent ion indicator and analog of Fluo-4. The indicator has an affinity for magnesium (Kd = 4.7 mM) and low-affinity for calcium (Kd = 22 uM) with absorbance/emission at 493/517 nm (high Mg²⁺). The cell membrane-impermeant tetrapotassium salt form can be introduced into cells by microinjection.

- λ_{Ex}/λ_{Em} (high Mg²⁺): 493/517 nm
- Orange solid soluble in water
- Store 4°C and protect from light
- C25H13F2K4NO10
- MW: 682

BAPTA-based ion indicators like Mag-Fluo-4 have been shown to be fixable in situ by EDC/EDAC (cat# 59002). The fixation of indicator dyes is useful for downstream immunofluorescence and IHC studies (Cell Calcium 1997, 21(3), 175).

As the indicator does not covalently bind to cellular components, it may be actively effluxed from the cell by organic anion transporters. The rate of efflux increases with temperature, and may vary between cell types, resulting in variable retention times of a few minutes to hours. Experiments using indicators in cells usually are performed within one or two hours of loading, but it may be possible to re-load cells with indicator if needed. The organic anion transporter inhibitor Probenecid (#50027) can be used to slow the rate of indicator efflux from cells.

References

- 1. Biophys J. (2009) Oct 7; 97(7):1864 doi: 10.1016/j.bpj.2009.07.021 2. Biochem J. (2002) Oct 1; 367(Pt1):137 doi: 10.1042/BJ20020305
- 3. Biochima et Biophys Acta Mol Cell Res. (2017) June; 1864(6):977 doi.org/10.1016/j.bbamcr.2016.11.026
- 4. Biochem & Biophys Res Comm. (2014) November 28; 454(4):572
- doi.org/10.1016/j.bbrc.2014.10.125 5. Methods Cell Biol, 99, 113, (2021), DOI: 10.1016/B978-0-12-374841-6.00005-0

This datasheet was generated on November 9, 2025 at 08:50:12 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mag-fluo-4-tetrapotassium-salt/