DMNP-EDTA, Tetrapotassium Salt (Caged Calcium)

DMNP-EDTA (also known as DM-NitrophenTM) is a caged calcium chelator. Upon photolysis, the Kd for Ca²⁺ increases from 5 nM to 3 mM, resulting in a pulse of free Ca²⁺.



Product attributes

Call us: 800-304-5357

117367-86-9

Email: btinfo@biotium.com

CAS number Cell permeability Membrane impermeant

Product Description

DMNP-EDTA (also known as DM-NitrophenTM) is a caged Ca²⁺ chelator. Upon photolysis, the Kd for Ca²⁺ increases from 5 nM to 3 mM, resulting in a pulse of free Ca²⁺.

- Light yellow solid soluble in water (pH >6) and DMSO
- Store desiccated at -20°C and protect from light
- Aqueous solutions stable for at least three months when stored at 4°C, protected from light
- C₁₈H₁₉K₄N₃O₁₂
- MW: 626

Biotium also offers a membrane-permeant DMNP-EDTA (Caged Calcium), AM Ester that may be used for live cell studies. Once inside the cell, intracellular esterases will hydrolyze the AM group. The resulting AM ester compound is then contained inside the cell and can accumulate.

DM-Nitrophen is a trademark of MilliporeSigma.

This datasheet was generated on November 16, 2025 at 10:49:50 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dmnp-edta-tetrapotassium-salt-caged-calcium/