



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

DMNP-EDTA, Tetrapotassium Salt (Caged Calcium)

DMNP-EDTA (also known as DM-Nitrophen™) is a caged calcium chelator. Upon photolysis, the K_d for Ca^{2+} increases from 5 nM to 3 mM, resulting in a pulse of free Ca^{2+} .



Product attributes

| | |
|-------------------|---------------------|
| CAS number | 117367-86-9 |
| Cell permeability | Membrane impermeant |

Product Description

DMNP-EDTA (also known as DM-Nitrophen™) is a caged Ca^{2+} chelator. Upon photolysis, the K_d for Ca^{2+} increases from 5 nM to 3 mM, resulting in a pulse of free Ca^{2+} .

- 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
- $C_{18}H_{19}K_4N_3O_{12}$
- 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 January 8, 2026 at 03:46:53 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dmnp-edta-tetrapotassium-salt-caged-calcium/>