

Product Information

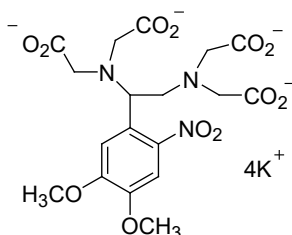
DMNP-EDTA, tetrapotassium salt (Caged Calcium)

Catalog Number: 50050

Unit Size: 5 mg

Molecular Information:

Molecular formula: $C_{18}H_{19}K_4N_3O_{12}$
Molecular weight: 626



Storage and Handling

Store DMNP-EDTA tetrapotassium salt at -20°C , protected from light. Stock solutions may be prepared in water.

Product Application

DMNP-EDTA (also known as DM-Nitrophen®) is a photolyzable calcium chelator that can be used to release a pulse of calcium upon light exposure. Upon photolysis, the K_d of DMNP-EDTA for calcium increases from 5 nM to 3 mM. DMNP-EDTA may also be useful for caged magnesium studies, with a K_d for magnesium of 2.5 μM .

References

Chem Rev (2008) 108:1603; Nat Methods (2007) 4:619.

DM-Nitrophen is a registered trademark of Calbiochem-Novabiochem Corp.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

Other Related Products

Biotium offers a wide selection of calcium indicator dyes and related products. Please visit our website at www.biotium.com for details.