Calcein AM

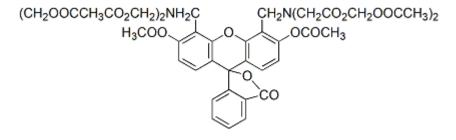
Calcein AM is a widely used green fluorescent cell marker. Calcein AM is membrane-permeant and thus can be introduced into cells via incubation.



Product Description

Calcein AM is a widely used green fluorescent cell marker. Calcein AM is membrane-permeant and thus can be introduced into cells via incubation. Once inside the cells, calcein AM is hydrolyzed by endogenous esterase into the highly negatively charged green fluorescent calcein, which is retained in the cytoplasm. Calcein AM is an excellent tool for the studies of cell membrane integrity and for cell tracing. Supplied in lyophilized form or as a solution in anhydrous DMSO at 4 mM (80011-1) or 1 mg/mL (1 mM) (80011-2). Calcein AM is also offered in our Cell Viability and Cytotoxicity Assay Kit (30002).

- λ_{Ex}/λ_{Em} (calcein) = 494/517 nm (pH 8)
- Colorless solid soluble in DMSO
- Store at -20°C and protect from light
- C46H46N2O23
- MW: 995



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

Product attributes

CAS number	148504-34-1
Probe cellular localization	Nucleus & cytoplasm
Assay type/options	High-throughput assay, Short term staining (<24h)
Cell permeability	Membrane permeant
Apoptosis/viability marker	Live cell stain
Colors	Green
Excitation/Emission	494/517 nm (pH 8)

This datasheet was generated on November 7, 2025 at 06:18:47 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/calcein-am/