Calcein (High Purity)

Calcein is a highly negatively charged, water-soluble fluorescein derivative. Its fluorescence is nearly independent of pH in the range between 6.5 and 12.



Product attributes

CAS number	154071-48-4
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green
Excitation/Emission	494/517 nm (pH 8)

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

Product Description

Calcein is a highly negatively charged, water-soluble fluorescein derivative. Its fluorescence is nearly independent of pH in the range between 6.5 and 12. The dye can be introduced into cells by microinjection and is useful for cell tracing and for the study of cell membrane integrity and fusion. Biotium also offers calcein AM, the membrane-permeant version of calcein that hydrolyzes into calcein once inside cells.

- $\lambda_{Ex}/\lambda_{Em} = 494/517 \text{ nm (pH 8)}$
- $\varepsilon = 75,000$
- Yellow-orange solid soluble in DMSO, DMF or pH>6 water
- Store at 4°C and protect from light
- \bullet C₃₀H₂₆N₂O₁₃
- MW: 623
- [154071-48-4]

This datasheet was generated on November 12, 2025 at 12:09:52 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/calcein-high-purity/