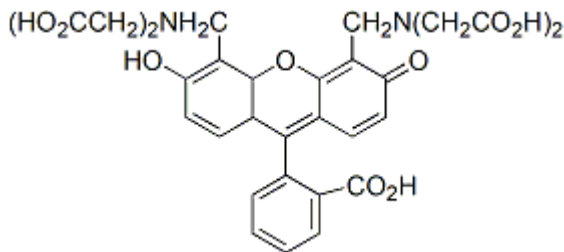


Calcein is a highly negatively charged, water-soluble fluorescein derivative. Its fluorescence is nearly independent of pH in the range between 6.5 and 12.



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.

- $\lambda_{Ex}/\lambda_{Em} = 494/517 \text{ nm (pH 8)}$
- $\epsilon = 75,000$
- Yellow-orange solid soluble in DMSO, DMF or pH>6 water
- Store at 4 °C and protect from light
- $C_{30}H_{26}N_2O_{13}$
- MW: 623
- [154071-48-4]



This datasheet was generated on January 9, 2026 at 03:23:39 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/calcein-high-purity/>

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)