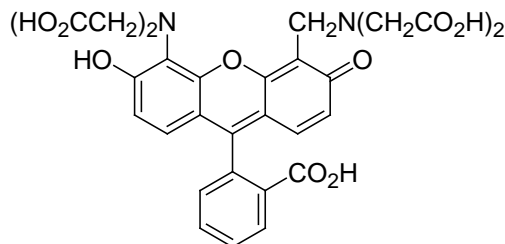


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Calcein**CATALOG #:** 80013**MOLECULAR INFORMATION:** C₃₀H₂₆N₂O₁₃
MW: 623**PROPERTIES:**

Color & Form	Yellow-orange solid
Purity	≥ 95% by TLC
Solubility	Soluble in DMSO, DMF or water (pH>6)
Absorption/Emission	494/517 nm (pH 8)
Extinction Coefficient	60,000

STORAGE AND HANDLING: Store desiccated at 4°C. Protect from light.

APPLICATION: 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 the membrane-permeant version of calcein: calcein AM (**80011**), which hydrolyzes into calcein once inside cells.

TOXICITY: Unknown.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--