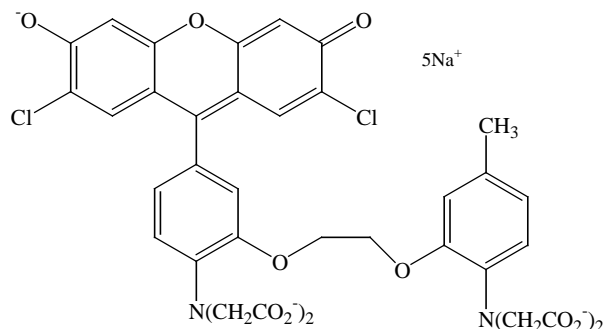


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Fluo-3, pentasodium salt**CATALOG #:** 50012**MOLECULAR INFORMATION:** C₃₆H₂₅Cl₂N₂Na₅O¹³
MWt: 879**PROPERTIES:**

Color & Form Orange red solid
Purity ≥ 95% by HPLC
Solubility Soluble in water or DMSO
Absorption/Emission 506nm/526nm (low or high [Ca²⁺])
Extinction Coefficient 86,000 M⁻¹cm⁻¹ (506 nm)

STORAGE AND HANDLING:

Store desiccated at 4 °C upon receipt. Protect from light, especially when in solution.

APPLICATION:

Fluo-3 has its absorption maximum at 506 nm, thus making it excitable by the argon-ion laser. Unlike fura-2 and indo-1, neither the excitation nor the emission maximum of the sensor shifts significantly before and after Ca²⁺ binding. As a result, the ratioing technique is not applicable to fluo-3. Fluo-3 is essentially nonfluorescent without Ca²⁺ present, but the fluorescence increases at least 40 times on Ca²⁺ binding. Also, because fluo-3 binds Ca²⁺ more weakly (higher K_d) than do fura-2 and indo-1, it is more useful for measuring high transient Ca²⁺ concentration during Ca²⁺ spikes.

Fluo-3 pentasodium salt is membrane-impermeant but can be loaded into cells via microinjection or scrape loading.

Ref: 1) Zucker, R.S., et al. *Cell Calcium*. **13**, 29(1992); 2) Merritt, J.E., et al. *Biochem. J.* **269**, 513(1990); 3) Lattanzio, F.A., et al. *Biochem. Biophys. Res. Comm.* **171**, 102(1990); 4) Jaffe, L. et al. *Proc. Natl. Acad. Sci. USA* **88**, 9883(1991). 5) Minta, A., et al. *J. Biol. Chem.* **264**, 8171(1989); 6) Kao, J.P.Y., et al. *J. Biol. Chem.* **264**, 8179(1989).

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: 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.
