46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Fluo-4, pentapotassium salt

CATALOG #: 50019

**MOLECULAR**  $C_{36}H_{25}F_2K_5N_2O_{13}$ MWt: 927.09 **INFORMATION:** 

CH<sub>3</sub>  $N(CH_2CO_2)_2$ 

PROPERTIES:

Color & Form Orange solid **Purity** > 95% by HPLC

**Solubility** Soluble in water or DMSO

494nm/516nm (low or high [Ca<sup>2+</sup>]) **Absorption/Emission** 

**Extinction** 82,000 M<sup>-1</sup>cm<sup>-1</sup> (494 nm) Coefficient

Kd for Ca<sup>2+</sup> ~335 nM

STORAGE AND **HANDLING:** 

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

APPLICATION:

Fluo-4 has its absorption maximum at 494 nm, thus making it excitable by the argonion laser. Fluo-4 is essentially nonfluorescent without Ca<sup>2+</sup> present, but the fluorescence increases at least 100 times on Ca<sup>2+</sup> binding. Also, because fluo-4 binds Ca<sup>2+</sup> more weakly (higher K<sub>d</sub>) than do fura-2 and indo-1, it is more useful for measuring high transient Ca<sup>2+</sup> concentration during Ca<sup>2+</sup> spikes.

Fluo-4 pentapotassium salt is membrane-impermeant but can be loaded into cells via microinjection or scrape loading.

Ref: 1) Gee, K.R., et al. Cell Calcium. 27, 97(2000).

Unknown **TOXICITY:** 

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on FIRST AID:

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.