

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: Indo-1, pentasodium salt

CATALOG #: 50042

MOLECULAR $C_{32}H_{26}Na_5N_3O_{12}$ **INFORMATION:** MWt: 759.5

[132319-56-3]

O₂CCH₃)₂N N(CH₂CO₂-)₂ O O CH₃ NH 5Na⁺

PROPERTIES:

Color & FormPale pink solidPurity $\geq 95\%$ by HPLCSolubilitySoluble in water

Absorption/Emission 349 nm/475 nm (no Ca²⁺); 330 nm/400 nm (high Ca²⁺) **Extinction** 349 nm/475 nm (no Ca²⁺); 330 nm/400 nm (high Ca²⁺) 33,000 M ⁻¹ cm ⁻¹ (high Ca²⁺)

Coefficient

STORAGE AND HANDLING:

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

APPLICATION:

Similar to fura-2, indo-1 is also a UV-excitable fluorescent Ca^{2+} sensor. However, different from fura-2, the fluorescent emission maximum undergoes a large blue shift from 482 nm to 398 nm upon Ca^{2+} binding. Thus, Ca^{2+} concentration can be determined by ratioing the fluorescence intensities at the two wavelengths. As with fura-2, this ratioing technique avoids problems associated with uneven dye distribution, cell or tissue thickness and photobleaching. Indo-1 has been widely used in flow cytometry studies.

Indo-1, pentasodium salt is membrane-impermeant but can be loaded into cells via microinjection or scrape loading.

Ref.: Babcock, D.R., et al. J. Biol. Chem. 262, 15041(1987).

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.

Indo-1, pentasodium salt PSF006

Revised: July 23, 2018

Page 1 of 1