

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, AM Ester

CATALOG #: 50043/50044

**MOLECULAR**  $C_{47}H_{51}N_3O_{22}$  **INFORMATION:** MWt: 1010

[112926-02-0]

## **PROPERTIES:**

Color & FormLight yellow solidPurity $\geq 96\%$  by HPLCSolubilitySoluble in DMSO

**Absorption/Emission** 356 nm/478 nm (no  $Ca^{2+}$ ); 330 nm/400 nm (high  $Ca^{2+}$ ) Extinction 33,000 M  $^{-1}$  cm  $^{-1}$  (no  $Ca^{2+}$ ); 33,000 M  $^{-1}$  cm  $^{-1}$  (high  $Ca^{2+}$ ) Coefficient

STORAGE AND

HANDLING:

Store desiccated at -20 °C. Protect from light, especially when in solution

**APPLICATION:** Membrane-permeant

Membrane-permeant form and thus can be loaded into cells via incubation. Because of the relatively low water solubility of the AM ester, Pluronic F-127 (See **59000** and refs. therein), a mild detergent, is often used as a dispersing agent to facilitate the loading. Indo-1/AM itself does not bin d Ca<sup>2+</sup>, but it is readily hydrolyzed to indo-1 by endogenous esterases once the dye is inside the cells.

Ref.: Grieson, J.P., et al. J. Neurophysiol. 67, 704(1992); Babcock, D.R., et al. J. Biol. Chem. 262, 15041(1987); Lazzari, K.G., et al. J. Biol. Chem. 261, 9710(1986).

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.