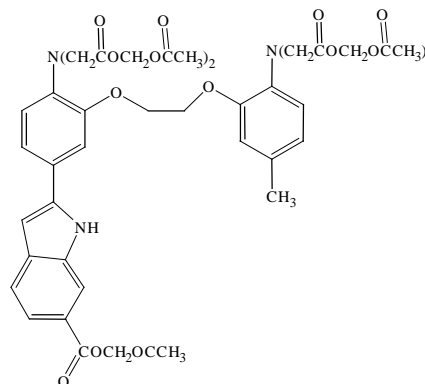


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Indo-1, AM Ester**CATALOG #:** 50043/50044**MOLECULAR INFORMATION:** C₄₇H₅₁N₃O₂₂
MWt: 1010
[112926-02-0]**PROPERTIES:**

Color & Form	Light yellow solid
Purity	≥ 96% by HPLC
Solubility	Soluble in DMSO
Absorption/Emission	356 nm/478 nm (no Ca ²⁺); 330 nm/400 nm (high Ca ²⁺)
Extinction Coefficient	33,000 M ⁻¹ cm ⁻¹ (no Ca ²⁺); 33,000 M ⁻¹ cm ⁻¹ (high Ca ²⁺)

STORAGE AND HANDLING:

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

APPLICATION:

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 bind Ca²⁺, 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.
