

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: Propidium Iodide, 1.0mg/ml solution in water

**CATALOG #** 40017

**MOLECULAR**  $C_{27}H_{34}I_2N_4$  **INFORMATION:** MWt: 668.4

[25535-16-4]

H<sub>2</sub>N — NH<sub>2</sub>

N<sub>+</sub> CH<sub>2</sub>CH<sub>3</sub>

CH<sub>2</sub>CH<sub>2</sub>CH<sub>2</sub>N — CH<sub>3</sub>

2I CH<sub>2</sub>CH<sub>3</sub>

PROPERTIES:

Color & FormOrange red liquidPurity $\geq 95\%$  by HPLC

**Soluble** in water or DMSO

**Absorption/Emission** 535/617nm (DNA-bound); 493/636nm (no DNA)

**Extinction** 5400 (DNA-bound); 5900 (no DNA)

Coefficient

STORAGE AND Stored d
HANDLING:

Stored desiccated at 4  $^{\circ}\text{C}$  and protect from light, especially when in solution.

**APPLICATION:** Propidium Iodide is a membrane-impermeant nucleic acid intercalator. The dye is

commonly used to selectively stain dead cells in a cell population and also used as a

counterstain in multicolor fluorescent imagings.

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.