

# Product Information

## Propidium Iodide Buffer, 50 µg/ml

<b>Catalog Number:</b>	40048
<b>Packaging Size:</b>	2 ml, filtered (0.45 micron)
<b>Composition:</b>	50 µg/ml propidium iodide in phosphate buffered saline (PBS)
<b>Molecular Information:</b>	C <sub>27</sub> H <sub>34</sub> I <sub>2</sub> N <sub>4</sub> MW: 668.4 [25535-16-4]
<b>Spectral Properties:</b>	Abs/Em (nm): 535/617 (with DNA) Abs/Em (nm): 493/636 (w/o DNA)
<b>Extinction Coefficient:</b>	5400 (with DNA) 5900 (w/o DNA)
<b>Storage:</b>	Store at 2-8°C, protect from light
<b>Caution:</b>	Propidium iodide is toxigenic and mutagenic. Handle with care.

## Product Description

Propidium iodide (PI) is a nucleic acid intercalator. It is cell impermeable and is excluded from viable cells. PI can be used in apoptotic studies to distinguish necrotic cells or cells in the late stages of apoptosis from viable cells or cells in early apoptosis where cell membranes are still intact. PI is also utilized as a counterstain in multicolor fluorescent imaging.

## Application

This product can be used in flow cytometry in conjunction with our series of Annexin V products: Annexin V, CF488A conjugate (cat. #29005), Annexin V, biotin conjugate (cat. #29013), and 5X Annexin V Binding Buffer (cat.#99902).

## Related Products

Cat.#	Product Name	Unit Size
29013	Annexin V, biotin conjugate, 50 µg/ml	0.5 ml
29005	Annexin V, CF488A conjugate, 50 µg/ml	0.5 ml
29001	Annexin V, FITC conjugate (FITC Annexin V), 50 µg/ml	0.5 ml
30017	Apoptosis & Necrosis Quantification Kit	50 assays
30018	Apoptotic, Necrotic & Healthy Cells Quantification Kit	50 assays
30060	CF488A Annexin V and 7-AAD Apoptosis Kit	100 assays
30061	CF488A Annexin V and PI Apoptosis Kit	100 assays
99902	5X Annexin V Binding Buffer	15 ml
10400	NucView™ 488 Caspase-3 Enzyme Substrate 1 mM in 1X PBS	100 µl
30030	NucView™ 488 Caspase-3 substrate and sulforhodamine 101-Annexin V Dual Apoptosis Kit	50-250 assays
30027	Viability/Cytotoxicity Assay Kit for Bacteria Live and Dead Cells	100-1000 assays

**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.