

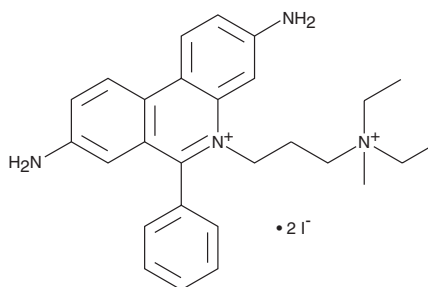
# PRODUCT INFORMATION



## Propidium Iodide Solution

Item No. 10008351

|                          |   |
|--------------------------|---|
| <b>Synonym:</b>          | PI Solution   |
| <b>Formulation:</b>      | 1 mg/ml solution in PBS, pH 7.4,<br>5 ml total                      |
| <b>Purity:</b>           | ≥95%  |
| <b>Application:</b>      | Nuclei counter stain for fluorescent<br>immunocytochemical staining |
| <b>Working Dilution:</b> | 1:200-1:1,000   |
| <b>Storage:</b>          | 4°C   |
| <b>Stability:</b>        | ≥2 years  |



### Description

Propidium Iodide (PI) binds to double-stranded DNA and RNA after cells have been permeabilized. Once bound to nucleic acids, the nuclei appear red in color as observed by fluorescence microscopy with a rhodamine filter. PI is useful for multi-color fluorescent staining analysis. The bright red staining of nuclei makes this an excellent counter stain to be used simultaneously with FITC, PE, or Cy5. For analysis by flow cytometry, PI staining can be monitored in the FL2 channel. PI can also be used to measure cell viability since this dye is excluded from viable cells.

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/22/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM