

## CERTIFICATE OF ANALYSIS

**Catalog Number** C007  
**Product Name** propidium iodide  
**CAS Number** 25535-16-4  
**Molecular Formula** C<sub>27</sub>H<sub>34</sub>I<sub>2</sub>N<sub>4</sub>  
**Molecular Weight** 668.4  
**Appearance** purple solid  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup>		
Maximum	494 nm	493 ± 3 nm
Extinction	6100 cm <sup>-1</sup> M <sup>-1</sup>	6000 ± 300 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup>		
Emission Maximum	632 nm	635 ± 4 nm
<b>HPLC</b>		
Purity	95 % at 254 nm	≥ 92 % at 254 nm
<b>NMR</b>		
Result	meets specification	consistent with structure

1. Solvent: Buffer, 50 mM potassium phosphate, pH 7.