



**Applied BioProbes**  
Reagents for Life Sciences

info@abpbio.com  
www.abpbio.com

## CERTIFICATE OF ANALYSIS

**Catalog Number** N100  
**Product Name** GreenView DNA Gel Stain  
**Appearance** orange solution  
**Lot Number** A300LY170

	LOT DATA	SPECIFICATION
<b>Absorption</b> Optical Density <sup>1</sup>	0.7 at 500 nm	0.65 ± 0.05 at 500 ± 3 nm

1. Solvent: water. Method: Optical Density, 4000× dilution.