



Applied BioProbes
Reagents for Life Sciences

info@abpbio.com
www.abpbio.com

CERTIFICATE OF ANALYSIS

Catalog Number	N101
Product Name	GreenView Plus DNA Gel Stain
Appearance	orange solution
Lot Number	A300LY171

	LOT DATA	SPECIFICATION
Absorption Optical Density ¹	0.46 at 500 nm	0.45 ± 0.05 at 500 ± 3 nm

1. Solvent: water. Method: Optical Density, 1000× dilution.