



Applied BioProbes
Reagents for Life Sciences

info@abpbio.com
www.abpbio.com

CERTIFICATE OF ANALYSIS

Catalog Number N102
Product Name RedView DNA Gel Stain
Appearance purple solution
Lot Number A300LY172

	LOT DATA	SPECIFICATION
Absorption Optical Density ¹	0.19 at 500 nm	0.19 ± 0.02 at 500 ± 3 nm

1. Solvent: water. Method: Optical Density, 500× dilution.