

## CERTIFICATE OF ANALYSIS

**Catalog Number** C013-1  
**Product Name** NucBeacon™ Green Live nucleic acid stain, 1 mM in DMSO  
**Appearance** orange solution  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption<sup>1</sup></b> Optical Density	0.63 at 494 nm	0.60 ± 0.03 at 494 ± 4 nm

1. Solvent: Buffer, 50 mM potassium phosphate, pH 7. Method: Optical Density. 100× dilution.