

## CERTIFICATE OF ANALYSIS

**Catalog Number** C047  
**Product Name** LysoBeacon™ Blue, 1 mM in DMSO  
**Appearance** colorless solution  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption<sup>1</sup></b> Optical Density	0.48 at 373 nm	0.48 ± 0.04 at 373 ± 3 nm

1. Solvent: Methanol. Method: Optical Density. 26x dilution