

---

## CERTIFICATE OF ANALYSIS

**Catalog Number** C048  
**Product Name** LysoBeacon™ Green, 1 mM in DMSO  
**Appearance** yellow solution  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption<sup>1</sup></b> Optical Density	0.24 at 458 nm	0.25 ± 0.03 at 457 ± 3 nm

1. Solvent: Methanol. Method: Optical Density. 100x dilution