

## CERTIFICATE OF ANALYSIS

**Catalog Number** C042  
**Product Name** MitoBeacon™ Red, 1 mg/mL in DMSO  
**CAS Number** 167095-09-2  
**Molecular Formula** C<sub>32</sub>H<sub>32</sub>Cl<sub>2</sub>N<sub>2</sub>O  
**Molecular Weight** 531.52  
**Appearance** purple solution  
**Lot Number** A300LY209

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup> Maximum Extinction	578 nm 111,000 cm <sup>-1</sup> M <sup>-1</sup>	578 ± 3 nm 117,000 ± 8,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup> Emission Maximum	600 nm	600 ± 4 nm
<b>HPLC</b> Purity	95 % at 578 nm	≥ 90 % at 578 nm
<b>NMR</b> Result	meets specification	consistent with structure

1. Solvent: Methanol.