

## Certificate of Analysis

### CAT# 307-L Aequorin Photoprotein in liquid

Date of Test:	23 Jan 2019
Compound ID:	Calcium Activated Aequorin Photoprotein Fully formed & ready to make light from Ca <sup>+2</sup> binding
Quantification:	Supplied as 0.5 mg in 250 µl (2mg/ml) in 1.2 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub> ABS <sub>280</sub> of 3.00 = 1.0 mg/ml of Aequorin
Production:	recombinant protein from E.coli purified via multi-step chromatography
EC Number:	1.13.12.5 (same as Renilla luciferin 2 monooxygenase)
CAS#	Not Registered
Storage:	Store in the dark at -20°C or for long term storage at -80°C. avoid repeated freeze/thaw cycles
Appearance:	dense yellow liquid
Maximal Turnover:	1.2 X 10 <sup>16</sup> QPS/mg using excess Calcium Chloride

#### SDS Gel Electrophoresis:

