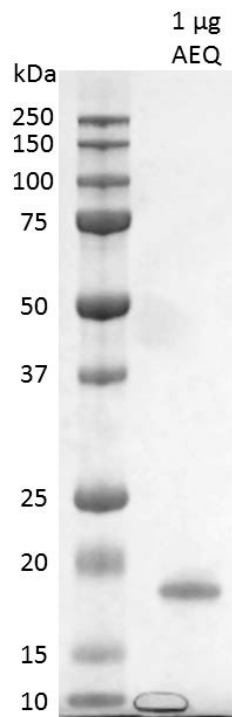


Certificate of Analysis

CAT#307-P Aequorin Protein lyophilized

Aequorin, a photoprotein originating from the jellyfish *Aequorea victoria* emits light in the presence of a trace amount of Ca^{2+} without the requirement of any other cofactor. Aequorin was "charged" with unmodified (native) Coelenterazine and will emit light at 465 nm upon Ca^{2+} contact.

Compound:	<i>Aequorea victoria</i> photoprotein (untagged)
Source:	recombinantly produced in <i>E. coli</i> , purified via multi-step chromatography
Sequence:	MTSKQYSVKLTSDFDNPRWIGRHKHMFNFLDVNHNGKISLDEMVKASDIVINN LGATPEQAKRHKDAVEAFFGGAGMKYGVETDWPAYIEGWKKLATDELEKYAKNE PTLIRIWGDALFDIVDKDQNGAITLDEWKAYTKAAGIIQSSDCEETFRVCDIDES QLDVDEMTRQHLGFWYTMDPACEKLYGGAVP
Quantity:	100 µg per vial
Reconstitute:	in 100 µl of ddH ₂ O (18.2 MΩ, Ca^{2+} free) as 1 mg/ml solution
MW:	22.3 kDa (theoretical)
Storage:	The protein is freeze dried and sealed under vacuum in a 10 fold excess of lyophilization stabilizer. We recommend storage at $\leq -20^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles after reconstitution.
Purity:	>95% by Coomassie staining 1 µg was applied on a 12% SDS-Gel and Coomassie stained



Activity of Aequorin

1. Aequorin was reconstituted in 100 μ l distilled water resulting in a 1 mg/ml solution
2. dilutions of Aequorin (in 10 mM K_2HPO_4 , 50 mM NaCl, pH 7.2) were pipetted into a black 96-well plate with a total volume of 50 μ l
3. 100 μ l of 10 mM $CaCl_2$ in dilution buffer was injected
4. luminescence was integrated over 10 sec with 0.0 sec delay after injection

