

## Certificate of Analysis

### CAT# 316 Renilla reniformis Green Fluorescent Protein

Date: 17 July 2019  
 Compound ID: Renilla reniformis Green Fluorescent Protein (RR-GFP)  
 Batch Number: 62005  
 Manufacturer: Prolume Ltd  
 Genbank Link: <http://www.ncbi.nlm.nih.gov/protein/AAK54757.1>  
 Theoretical pI/Mw: 6.31 / 27123.19

#### Molecular Formula/Structure:

```

      10      20      30      40      50      60
MDLAKLGLKE VMPTKINLEG LVGDHAFSME GVGEGNILEG TQEVKISVTK GAPLPFAFDI
      70      80      90     100     110     120
VSVAFSYGNR AYTGYPEEIS DYFLQSFPEG FTYERNIRYQ DGGTAIVKSD ISLEDGKFIV
      130     140     150     160     170     180
NVDFKAKDLR RMGPVMQQDI VGMQPSYESM YTNVTSVIGE CIIAFKLQTG KHFTYHMRTV
      190     200     210     220     230
YKSKKPVETM PLYHFIQHRL VKTNVDTASG YVVQHETAIA AHSTIKKIEG SLP
  
```

Test Items	Specification	Results	Test method
<b>Appearance</b>	Green Solid	Conforms	Visual
<b>Purity</b>	>75%	76%	PAGE
<b>Identification</b>	Emission Spectrum	Conforms	Ex/Em
	Absorption @480nm/280nm >2	Conforms	By HPLC
<b>Storage</b>	Store dry and dark at -20°C or for long term storage at -80°C.		