

Certificate of Analysis

CAT#315 Pt-GFP-HIS

Date: April 14, 2019
 Compound ID: Ptilosarcus sp. Green Fluorescent Protein-HIS
 Batch Number: 42005
 Manufacturer: Prolume Ltd
 Genbank Link: <http://www.ncbi.nlm.nih.gov/protein/AAG54097.1>
 Theoretical pI/Mw: 5.91 / 27047.72

Molecular Formula/Structure:

```

      10      20      30      40      50      60
MNRNVLKNTG LKEIMSAKAS VEGIVNNHVF SMEGFGKGNV LFGNQLMQIR VTKGGPLPFA

      70      80      90      100     110     120
FDIVSIAFQY GNRTFTKYPD DIADYFVQSF PAGFFYERNL RFEDGAIVDI RSDISLEDDK

      130     140     150     160     170     180
FHYKVEYRGN GFPSNGPVMQ KAILGMEPSF EVVYMNSGVL VGEVDLVYKL ESGNYYSCHM

      190     200     210     220     230
KTFYRSKGGV KEFPEYHFH HRLEKTYVEE GSFVEQHETA IAQLTTIGKP LGSLEHWV
  
```

Test Items	Specification	Results	Test method
Appearance	Green Solid	Conforms	Visual
Purity	>75%	89%	PAGE
Identification	Emission Spectrum	Conforms	Ex/Em
	Absorption @480nm/280nm >2	Conforms	By HPLC
Storage	Store dry and dark at -20°C or for long term storage at -80°C.		