

## Y1BFP, Protein

Recombinant full-length protein expressed in *E. coli* cells

### Catalog # B620-30BH

Lot # H2680-6P1

### Product Description

Recombinant full-length vibrio fischeri Y1BFP was expressed in *E. coli* cells using an N-terminal His tag. The Y1BFP protein accession number is [BAG71497](#).

### Gene Aliases

BFP; Y1-blue fluorescent protein

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, and 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

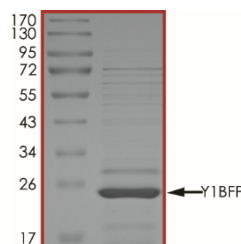
### Scientific Background

Y1BFP, a blue fluorescence protein from *Vibrio fischeri* Y1, was characterized to be responsible for the blue bioluminescence emission from *Photobacterium phosphoreum* (1). The Y1BFP, as a bound fluorophore, can bind with riboflavin, 6, 7-dimethyl-8-ribityllumazine (lumazine), and 6-methyl-7-oxo-8-ribityllumazine without riboflavin-synthase activity (2).

### References

1. Petushkov V.N., et al.: Purification and characterization of flavoproteins and cytochromes from the yellow bioluminescence marine bacterium *Vibrio fischeri* strain Y1. *Eur J Biochem.* 245(3):790-6, 1997.
2. <http://www.uniprot.org/uniprot/P80893>.

### Purity



The purity of Y1BFP was determined to be **>80%** by densitometry, approx. MW **25 kDa**.

## Y1BFP, Protein

Recombinant full-length protein expressed in *E. coli* cells

Catalog #	B620-30BH
Lot #	H2680-6P1
Purity	>80%
Concentration	0.2 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**