

## RFP, Protein

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

### Catalog # R620-30H

Lot # H2726-7

### Product Description

Recombinant full-length Sea anemone RFP was expressed in *E. coli* cells using an N-terminal His tag. The RFP protein accession number is [AAF03369](#).

### Gene Aliases

drFP583; DsRed

### 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.

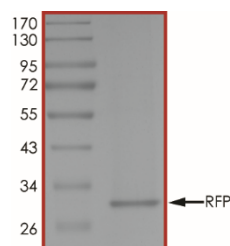
### Scientific Background

RFP (red fluorescent protein drFP583) has the protection role from the UV light and the conversion of suitable light for the microalgal symbionts in the coral reefs (1). The RFP proteins can be used in multicolor tagging experiments, avoiding natural cellular autofluorescence, and the resonance energy transfer-based experiments (2). The broad fluorescence emission spectrum peaks is at 583 nm (3).

### References

1. Matz, M.V., et al.: Fluorescent proteins from non-bioluminescent Anthozoa species. *Nat. Biotechnol.* 17 (10), 969-973, 1999.
2. Yarbrough D., et al.: Refined crystal structure of DsRed, a red fluorescent protein from coral, at 2.0-Å resolution. *Proc Natl Acad Sci USA.*; 98(2):462-7, 2001.
3. <http://www.uniprot.org/uniprot/Q9U6Y8>.

### Purity



The purity of RFP was determined to be **>90%** by densitometry, approx. MW **29 kDa**.

## RFP, Protein

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

Catalog #	R620-30H
Lot #	H2726-7
Purity	>90%
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.**