

BFP, Protein

Synthetic construct expressed in *E. coli* cells

Catalog # B620-30H

Lot # H2620-8P3

Product Description

Synthetic construct BFP (228-end) was expressed in *E. coli* cells using an N-terminal His tag. The BFP gene accession number is [DQ399412](#).

Gene Aliases

pBI-SS(Tom)(TP)101-EGFP

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°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.

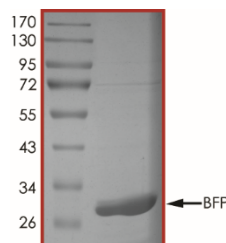
Scientific Background

BFP contains tomato signal sequence and Thr-Pro synthetic glycomodule fused to EGFP (1). When it is expressed in plants, many of the repetitive proline residues are 4-hydroxylated and O-glycosylated (1). The resulting proteoglycan is secreted into the cell wall (2).

References

1. Tan, L., et al: Glycosylation motifs that direct arabinogalactan addition to arabinogalactan-proteins *Plant Physiol.* 132 (3), 1362-1369, 2003.
2. Estevez, J.M., et al: Characterization of synthetic hydroxyproline-rich proteoglycans with arabinogalactan protein and extensin motifs in *Arabidopsis*. *Plant Physiol.* 142 (2), 458-470, 2006.

Purity



The purity of BFP was determined to be **>90%** by densitometry, approx. MW **28 kDa**.

BFP, Protein

Synthetic construct expressed in *E. coli* cells

Catalog #	B620-30H
Lot #	H2620-8P3
Purity	>90%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°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.