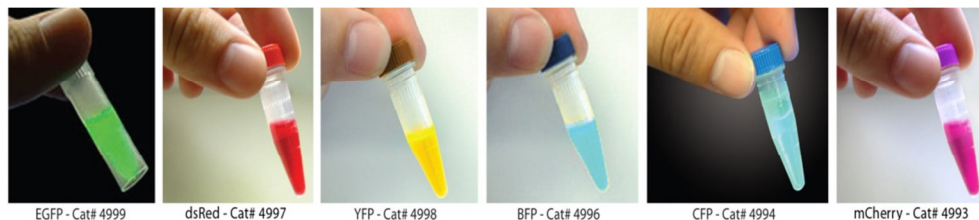


Fluorescent Protein Set

CATALOG #: K816-6-100
AMOUNT: 6 x 100 µg
LOT #: _____
SOURCE: *E. coli*
PURITY: ≥97% by SDS-PAGE and HPLC
ENDOTOXIN: ≤0.1 ng/ µg
FORM: Freeze Dried
RECONSTITUTION: Reconstitute with dH2O to 1 mg/ml
STORAGE: -80°C for long-term storage

CONTENTS:

| Protein | Color | #aa | Ex/Em (nm) | MW (kDa) | His-tag | ϵ (M ⁻¹ cm ⁻¹) at 280 nm | Part# |
|---------|--------|-----|-----------------|----------|----------|--|----------|
| EGFP | Blue | 293 | 488/507 | 32.7 | N/C-term | 21890 | 4999-100 |
| BFP | Orange | 259 | 308-383/440-447 | 29 | N-term | 19035 | 4994-100 |
| CFP | Purple | 284 | 458/480 | 31.3 | N-term | 26000 | 4996-100 |
| dsRed | Red | 245 | 557/585 | 27.6 | N-term | 34380 | 4997-100 |
| mCherry | Red | 256 | 587/610 | 28.8 | N-term | 34380 | 4993-100 |
| YFP | Yellow | 238 | 525/538 | 26.4 | None | 27390 | 4998-100 |



EGFP - Cat# 4999 dsRed - Cat# 4997 YFP - Cat# 4998 BFP - Cat# 4996 CFP - Cat# 4994 mCherry - Cat# 4993

PRODUCT DESCRIPTION:

100 µg of each of the six fluorescent proteins including the Enhanced Green fluorescent protein (EGFP), dsRed and mCherry Red fluorescent proteins, Blue fluorescent protein (BFP), Yellow fluorescent protein (YFP) and Cyan fluorescent protein (CFP). Recombinant fluorescent proteins were expressed and purified from transformed *E. coli* using a method that ensures highest purity and maximal fluorescence intensity. Fluorescent proteins can be used in a variety of applications. Fluorescent proteins antibodies and set are also available separately.

APPLICATIONS:

These proteins are suitable for a variety of applications including as control reagents for fluorescent protein expression studies, as standards for SDS-PAGE and Western blot analysis, as well as use for calibration of fluorometers and flow cytometers, fluorescence microscopy, and microinjection of fluorescent proteins into cells and tissues, etc. The recombinant fluorescent proteins can also be conjugated to other proteins. The presence of His-tag on most of these proteins allows their detection by both His-tag antibody and specific fluorescent protein antibodies on Western Blot analysis

RELATED PRODUCTS:

- Fluorescent Proteins
- Fluorescent Protein Conjugates (Biotin & FITC)
- Fluorescent Protein Antibodies
- His-tag Antibody
- GFP Quantification Kit
- pVision-GFP-C Vector
- pVision-GFP-N Vector
- Apoptosis Assay
- Cell Proliferation & Cytotoxicity Assays
- Fractionation Kits
- Metabolism Assays
- Cell Proliferation & Senescence
- Cell Damage & Oxidative Stress Assays
- Signal Transduction Kits & Reagents
- Adipocyte & Lipid Transfer
- Molecular Biology Kits & Reagents
- Reporter Assays
- Growth Factors and Cytokines
- Polyclonal and Monoclonal Antibodies

FOR RESEARCH USE ONLY! Not to be used on humans.