

Product Information

R-Phycoerythrin (R-PE), Lyophilized Solid, Ammonium-Free

Catalog Number: 29091-10mg, 29091-100mg, 29091-1000mg

Unit Size: 10 mg, 100 mg, 10 x 100 mg

Storage and Handling

Store at 4°C and protect from light. Lyophilized solids are stable for at least 12 months from date of receipt when stored as recommended.

Stock solutions may be prepared in water or PBS. Solutions can be stored protected from light at 4°C, for at least one month. For longer-term storage, single-use aliquots can be stored at -20°C for at least 3 months. Aliquots should not be refrozen after thawing.

Molecular Weight: ~240,000

Color and Form: Pink solid

Solubility: Water or aqueous buffer

Absorption/Emission: 496, 546, 565/576 nm

Product Description

R-Phycoerythrin (R-PE or PE) is a highly stable and soluble fluorescent phycobiliprotein from red algae. Phycobiliproteins have exceptional fluorescence characteristics due to their biological role as light harvesters in cyanobacteria and certain algae. For this reason, phycobiliproteins like R-PE are commonly used to generate fluorescent probes through covalent conjugation to proteins and antibodies.

Ammonium sulfate is typically added as a stabilizer in most commercial R-PE products. However, under conditions commonly used for bioconjugation, ammonium will be converted to ammonia that can interfere with the labeling. Thus, R-PE with ammonium sulfate needs to undergo buffer exchange to remove the ammonium, which adds an additional tedious step to the conjugation process. For greater convenience, Biotium's R-PE is lyophilized in a proprietary buffer that is both ammonium-free and protective of the spectral and chemical properties of the fluorescent protein. No buffer exchange is required, simply dissolve the lyophilized R-PE in your buffer of choice for conjugation.

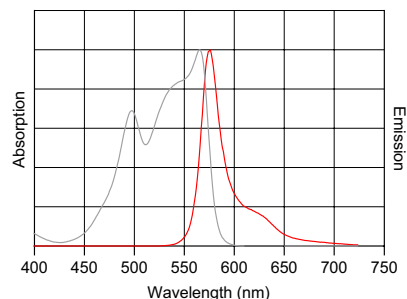


Figure 2. Absorbance (grey) and emission (red) spectra of R-PE.

Related Products

Cat. No.	Product
92452, 92298, 92299	Mix-n-Stain™ R-PE Antibody Labeling Kit
92442, 92443	Mix-n-Stain™ R-PE-CF@583R Antibody Labeling Kit
92340, 92341, 92346	Mix-n-Stain™ R-PE-CF@647T Antibody Labeling Kit
92453, 92306, 92307	Mix-n-Stain™ APC Antibody Labeling Kit
92230... 92463	Mix-n-Stain™ CF@ Dye Antibody Labeling Kits
92208... 92228	CF@ Dye & Biotin SE Protein Labeling Kits
29044	Streptavidin, R-PE Conjugate
29045	Annexin V, R-PE Conjugate
29092	Allophycocyanin (APC), Cross-linked Lyophilized Solid, Ammonium-Free
30071	AccuOrange™ Protein Quantitation Kit
22004	Ultrafiltration vial, 10K MWCO (pack of 5)
22018	Ultrafiltration vial, 3K MWCO (pack of 5)
41024-4L	Water, Ultrapure Molecular Biology Grade

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent CF@ Dye antibody conjugates, antibody labeling kits, reactive dyes, protein gel stains, and other products for life science research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.