



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Allophycocyanin (APC), Cross-linked Lyophilized Solid, Ammonium-Free

High-quality allophycocyanin (APC) in a lyophilized format. Provided in a ammonium-free and ready-to-use format for conjugation reactions.



Product attributes

Product Description

Allophycocyanin (APC) 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 APC are commonly used to generate fluorescent probes through covalent conjugation to proteins and antibodies.

- Lyophilized, high-quality APC
- MW: ~105,000
- Absorption/Emission: 651/660 nm
- Convenient ammonium-free format, ready-to-use for conjugation

Ammonium sulfate is typically added as a stabilizer in most commercial APC products. However, under conditions commonly used for bioconjugation, ammonium will be converted to ammonia that can interfere with the labeling. Biotium's APC is provided in a convenient lyophilized and ammonium-free format that is read-to-use for conjugation reactions.

Biotium also offers [ammonium-free R-Phycoerythrin \(R-PE\)](#), as well as [reactive formats](#) of bright and photostable [CF® Dyes](#). For streamlined and optimized antibody labeling, learn about our [Mix-n-Stain™ Antibody Labeling Kits](#).

This datasheet was generated on December 22, 2025 at 12:16:48 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/allophycocyanin-apc-cross-linked-lyophilized-solid-ammonium-free/>