

ReadiUse™ CL-APC [Cross linked-Allophycocyanin] *Ammonium Sulfate-Free*

Catalog Number: 2503, 2504

Unit Size: 1 mg, 10 mg

Product Details

Storage Conditions Refrigerated (2-8 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Powder blue

Molecular Weight ~105000

Soluble In Water (Add 100 µL of H₂O to Cat# 2503 or 1 mL to Cat# 2504 to reconstitute to 10 mg/mL in PBS)

Spectral Properties

Excitation Wavelength 651 nm

Emission Wavelength 660 nm

Applications

Cross linked-Allophycocyanin (CL-APC) is a phycobiliprotein isolated from *Spirulina* sp., a blue-green alga. Like other phycobiliproteins, APC is strongly fluorescent, with an extremely high absorptivity and a high quantum efficiency. It is a protein which can be easily linked to antibodies and other proteins by conventional protein cross-linking techniques without altering its spectral characteristics. APC is the least stable among the major phycobiliproteins, susceptible to dissociation at low concentrations including concentrations at which some assays are performed. CL-APC is chemically cross-linked between α and β subunits, and is much more stable than APC. The crosslinked APC has improved stability in aqueous solution. However, all the commercial APC materials are sold in concentrated ammonium sulfate buffers. The commercial APC materials from other vendors require the tedious dialysis or other purifications performed before it can be used for labeling purposes. AAT Bioquest offers this ReadiUse™ APC that can be readily used for any labelings without any purifications required. Our highly purified ReadiUse™ APC facilitates the rapid APC conjugations to antibodies and other proteins such as streptavidin and other secondary reagents.