
Product Information Sheet

Ordering Information

Product Number:	2503
Product Name:	ReadiUse™ CI-APC [Cross linked-AlloPhycocyanin] *Ammonium Sulfate-Free*
Unit Size:	1 mg
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	N/A
Soluble In:	Water
Excitation Wavelength:	651
Emission Wavelength:	662

Application Notes

Cross linked-AlloPhycocyanin (CI-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. CI-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.