



PRODUCT INFORMATION SHEET

<u>Product Number:</u>	AD0201
<u>Compound:</u>	Streptavidin conjugated to SureLight®-Allophycocyanin for protein binding assays
<u>Amount:</u>	1 mg total protein
<u>Lot Number:</u>	APC1P051
<u>Storage:</u>	4°C in the dark. After product has been reconstituted, DO NOT FREEZE.
<u>Handling:</u>	Before use, reconstitute to 1.0 ml with distilled deionized water and allow to sit on ice for 20-30 minutes. Gently mix by vortexing or tapping tube before use.
<u>Buffer:</u>	Supplied lyophilized. Upon reconstitution with 1.0 mL distilled deionized water, the product is in 10 mM Tris (pH 8.2), 150 mM sodium chloride, 50 mM sucrose, 0.1% BSA and 1 µg/mL pentachlorophenol.
<u>Stability:</u>	Upon reconstitution, product should be stored at 4°C and be used within 8 months.
<u>Date of Manufacture:</u>	September 12, 2024
<u>Shelf Life:</u>	5 years from test date
<u>Test Date:</u>	9/16/24

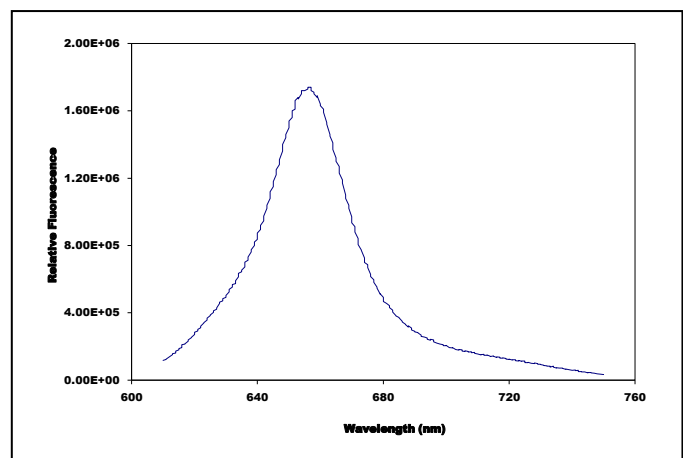
Intended use: The Streptavidin:SureLight®-Allophycocyanin conjugate can be used as an acceptor for the energy emitted by Eu³⁺-chelates in the LANCE™ homogeneous assay format. Product is for research purposes only, the product is not intended for any other use (such as cosmetic, agricultural, drug, food or household uses). Assay optimization should be based on the Streptavidin concentration. One can assume that 1 Streptavidin binds two biotinylated peptides or one biotinylated protein.

Lot specific information:
(upon reconstitution)

Concentration: 1.0 mg/ml total protein

SureLight®-APC molarity = 8.9 µM

Streptavidin molarity = 2.9 µM



SureLight® is a registered trademark of Columbia Biosciences Corporation
LANCE™ is a trademark of Revvity Health Sciences, Inc