

PRODUCT INFORMATION SHEET

Streptavidin conjugated to SureLight®-Allophycocyanin for kinase assays

<u>Product Number:</u>	CR130-150
<u>Compound:</u>	Streptavidin conjugated to SureLight®-Allophycocyanin for kinase assays
<u>Amount:</u>	50.0 mg total protein
<u>Lot Number:</u>	APC1K112
<u>Storage:</u>	4°C in the dark. After product has been reconstituted, DO NOT FREEZE.
<u>Handling:</u>	Before use, reconstitute to 25.0 ml with distilled deionized water and allow vial 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 25.0 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 with water, product should be stored at 4°C and be used within 8 months.
<u>Expiration:</u>	9/2020 when stored lyophilized.

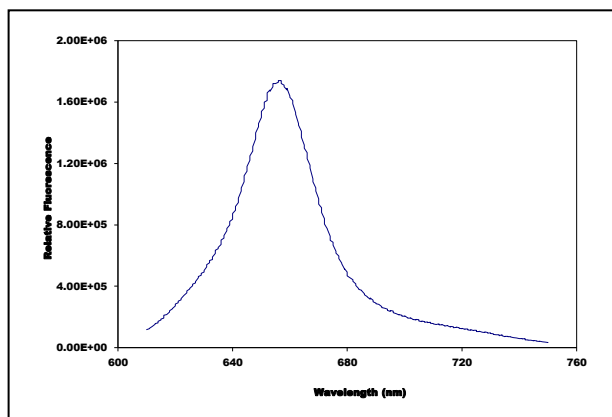
Intended use: The Streptavidin: SureLight®-Allophycocyanin conjugate can be used as an acceptor for the energy emitted by Eu³⁺-chelates in the LANCE™ homogeneous assay format. Product is for research purposes only, the product is not intended for any other use (such as cosmetic, agricultural, drug, food 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:
(after reconstitution)

Concentration: 2.07 mg/ml total protein


SureLight®-APC molarity = 12.4 µM

Streptavidin molarity = 13.9 µM



SureLight® is a registered trademark of Columbia Biosciences Corporation

Limited Warranty: This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.