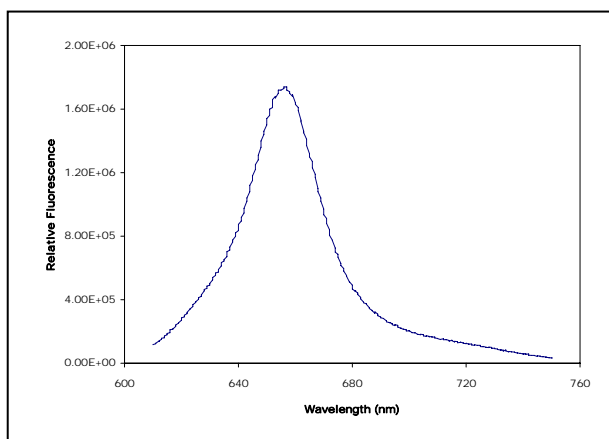


PRODUCT INFORMATION SHEET

<u>Product Number:</u>	CR130-100
<u>Compound:</u>	Streptavidin conjugated to SureLight®-Allophycocyanin for kinase assays
<u>Amount:</u>	1 mg total protein
<u>Lot Number:</u>	
<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 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 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 with water, product should be stored at 4°C and be used within 8 months.
<u>Expiration:</u>	3 years from date of manufacture when 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.



SureLight® is a registered trademark of Martek Biosciences Corporation
LANCE[™] is a trademark of Perkin Elmer Corporation