



PRODUCT INFORMATION SHEET

<u>Product Number:</u>	AD0059F
<u>Compound:</u>	Anti-FLAG® IgG conjugated to SureLight®-Allophycocyanin
<u>Amount:</u>	1 mg total protein
<u>Lot Number:</u>	APC1354
<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 with water, product should be stored at 4°C and be used within 3 months.
<u>Date of Manufacturing:</u>	October 9, 2024
<u>Shelf Life:</u>	5 years from test date
<u>Test Date:</u>	10/14/24

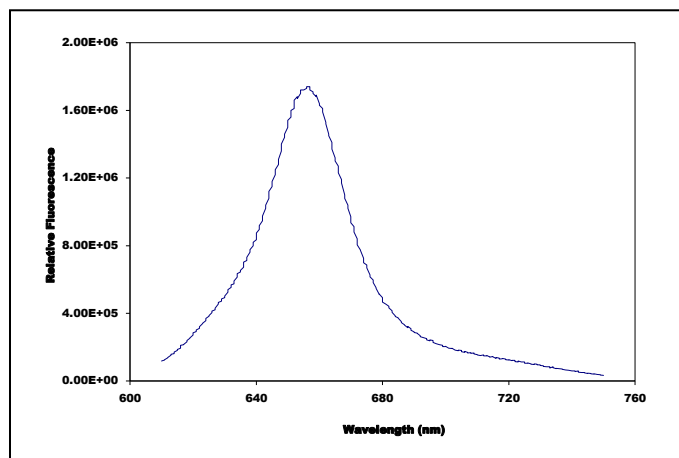
Intended use: The Anti-FLAG® IgG :SureLight®-Allophycocyanin conjugate can be used as an acceptor for the energy emitted by Eu³⁺-chelates in the LANCE™ homogeneous assay format. Product is for research purposes only, the product is not intended for any other use (such as cosmetic, agricultural, drug,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

Lot specific information:
(upon reconstitution)

Concentration: 1.09 mg/ml total protein

SureLight®-APC molarity = 4.15 µM

Anti-FLAG® IgG molarity = 4.10 µM



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