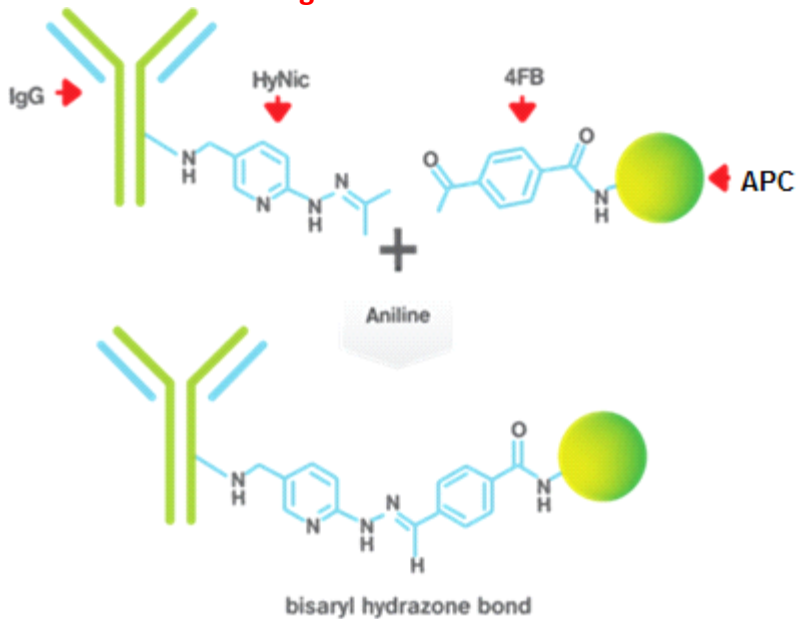


Product Data Sheet/Certificate of Analysis

APC-Antibody Conjugation Kit

Storage: **Store at 4°C. Protect 4FB-APC from light!**



Catalog Number: P-9903-001

Lot Number:

Components:	Component #:	Units:
S-HyNic	P-9903-001-01	1.0 mg
4FB-modified APC	P-9903-001-02	1.0 mg
10x Modification Buffer	P-9903-001-03	1.5 mL
10x Conjugation Buffer	P-9903-001-04	1.5 mL
0.5 mL Zeba spin Columns	P-9903-001-06	4
Anhydrous DMF	P-9903-001-07	1.0 mL
2-Sulfobenzaldehyde	P-9903-001-08	100 mg
10X MES Buffer	P-9903-001-09	1.5 mL
1.5 mL Collection Tubes	P-9903-001-10	12

Test	Specification	Result
Gel Analysis	Conjugation to Bovine IgG	Passed

QC Release Date:

Expiration Date:

Released By:



Product Description

4FB-modified APC is a 4-formylbenzamide-modified phycobiliprotein. This APC is covalently pre-activated with a precise number of aromatic aldehyde functional groups. 4FB-modified APC rapidly conjugates to HyNic-modified biomolecules to generate a highly fluorescent conjugate. Conjugates made with 4FB-APC retain their bright fluorescence.

Application

This kit is used for the conjugation of primary antibodies to APC for direct analysis in Flow, ELISA, immunofluorescence and Immunohistochemistry applications.

Comments

Store at 4°C. Protect 4FB-APC from light! Not for internal or external use in humans.

