

Product Information

Phaseolus Vulgaris Leucoagglutinin (PHA-L)

Product List

Cat. No.	Conjugate	Ex/Em (nm)
29115	CF@488A	490/516
29116	CF@568	562/584
29117	CF@594	593/615
29118	CF@640R	642/663
29119	CF@680	681/698
29114	Biotin	N/A

Unit Size: 1 mL

Storage and Handling

Store at 4°C, protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

Product Description

Phaseolus vulgaris (red kidney bean) agglutinin (phytohemagglutinin; PHA) is a family of lectins, each consisting of four subunits. There are two different types of subunits. The "E" (erythroagglutinin) subunit causes red cell agglutination, and the "L" (lymphocyte) subunit can trigger lymphocyte agglutination and mitogenic activity. These subunits combine to produce five isolectins. *Phaseolus vulgaris* Leucoagglutinin (PHA-L) has four "L" type subunits and does not agglutinate red blood cells but is a potent mitogen and triggers lymphocyte agglutination. Note that CF@ Dye PHA-L conjugates are formulated with sodium azide and are not suitable for use *in vivo* or live cell culture.

Biotium's next-generation fluorescent CF@ Dyes offer advantages in brightness and photostability compared to other fluorescent dyes. We also offer a variety of lectin conjugates, including *Lycopersicon esculentum* (Tomato) Lectin (LEL, TL), *Datura stramonium* lectin (DSL), *Ulex europaeus* Agglutinin I (UEA I), and PNA Lectin (*Arachis hypogaea*) CF@ Dye conjugates.

Conjugate Composition

Supplied at 1 mg/mL in 10 mM HEPES pH 7.5, 0.15 M NaCl, 0.08% sodium azide, 0.1 mM CaCl₂.

Experimental Protocol

PHA-L conjugate applications are reported in published literature. We recommend consulting the literature to find a protocol that will be suitable for your application. For immunofluorescence of tissue sections, we recommend staining with 5 µg/mL lectin conjugate in PBS for 15 minutes, followed by washing in PBS as a starting point for optimization.

Related Products

Cat. No.	Product
00070-00079	Cholera Toxin Subunit B CF@ Dye Conjugates
29015...29080	CF@ Dye Concanavalin A (Con A)
29030-29086	Streptavidin Conjugates
29060-29063	CF@ Dye PNA Lectin (<i>Arachis Hypogaea</i>)
29102-29107	<i>Lycopersicon Esculentum</i> (Tomato) Lectin (LEL, TL)
29096-29101	<i>Datura Stramonium</i> Lectin (DSL)
29021-29076	Wheat Germ Agglutinin (WGA) Conjugates
29108-29113	<i>Ulex Europaeus</i> Agglutinin I (UEA I)
30021-30024	CellBrite® Cytoplasmic Membrane Dyes
30088-30090	CellBrite® Fix Membrane Stains
30105-30109	CellBrite® Steady Membrane Staining Kits
30092-30104	MemBrite® Fix Cell Surface Staining Kits
29067	Calcofluor White
40043	DAPI, 10 mg/mL in Water
40046	Hoechst 33342, 10 mg/mL in Water
30015	DAB Substrate Kit
92170-96066	CF@ Dye Tyramides
22027	Ready-to-Use Tyramide Amplification Buffer
22016	Permeabilization Buffer
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF@ Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.