

Product Information

Sambucus Nigra Lectin (SNA, EBL)

See [product page](#) for a full list of product names and catalog numbers.

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.

Conjugate Composition

These conjugates are provided at a concentration of 1 mg/mL. The storage buffer contains 10 mM HEPES, 0.15 M NaCl, 0.08% sodium azide, 0.1 mM CaCl₂ (pH 7.5).

Product Technical Information

See [product page](#) for spectral properties and other dye-specific technical information. See our [Spectra Viewer](#) to view and download the dye excitation and emission spectra.

Product Description

Sambucus nigra lectin, also known as elderberry bark lectin (EBL) and *Sambucus nigra* agglutinin (SNA), belongs to the group of type 2 ribosome-inactivating proteins. It is composed of an A-chain with enzymatic activity and a B-chain with carbohydrate-binding activity, and has a combined molecular weight of ~140,000 Da. SNA binds preferentially to sialic acid attached to terminal galactose in α -2,6 and to a lesser degree, α -2,3 linkage. SNA can induce apoptosis in insect cells and cultured mammalian cell cancer cells. This effect depends on the carbohydrate binding of the B-chain.

Sambucus nigra lectin is available conjugated to biotin and a selection of bright and photostable CF® Dyes. CF® Dyes are Biotium's line of next-generation fluorescent dyes with advantages in brightness, photostability, and conjugate specificity compared to other fluorescent dyes.

Experimental Protocols

Sambucus nigra lectin 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 ug/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)
29060-29063	CF® Dye PNA Lectin (<i>Arachis hypogaea</i>)
29096-29101	<i>Datura Stramonium</i> Lectin (DSL)
29102-29107	<i>Lycopersicon Esculentum</i> (Tomato) Lectin (LEL, TL)
29021-29076	Wheat Germ Agglutinin (WGA) Conjugates
29108-29113	<i>Ulex Europaeus</i> Agglutinin I (UEA I)
29114-29119	<i>Phaseolus Vulgaris</i> Leucoagglutinin (PHA-L)
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
40083...41040	NucSpot® Nuclear Stains
29030-29086	Streptavidin Conjugates
29067	Calcofluor White
40043	DAPI, 10 mg/mL in Water
40046	Hoechst 33342, 10 mg/mL in Water
30131-30135	CytoLiner™ Fixed Cell Membrane Stains
92170-96066	CF® Dye Tyramides
22027	Ready-to-Use Tyramide Amplification Buffer

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.