

Isolectin B₄ (Bandeiraea simplicifolia), (FITC conjugate)

BS-I is a tetrameric lectin consisting of two types of subunits designated A and B. There are five BS-I isolectins with different subunit composition: BSI-B₄, BSI-AB₃, BSI-A₂B₂, BSI-A₃B and BSI-A₄. BSI-B₄ is blood group B specific and has an exclusive affinity for terminal α-D-galactosyl residues, whereas BSI-A₄ has blood group A specificity and has a major affinity for terminal N-acetyl-α-D-galactosaminy residues.

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-650-001F-MC05	FITC - 0.5mg
-------------------	--------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Prepare fresh dilutions prior to each application.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Activity	Agglutinates blood group B glutaraldehyde-fixed human red blood cells at a concentration of ~2µg/ml. No agglutination of blood group A, glutaraldehyde-fixed human red blood cells occurred at concentrations up to ~100µg/ml.
Alternative Name	Isolectin B4 (<i>Griffonia simplicifolia</i>), BSI-B4
Concentration	1mg/ml
Formulation	Liquid. In 10mM HEPES, pH 7.5, containing 0.15M sodium chloride, 0.08% sodium azide and 0.1mM Ca ²⁺ .
Quality Control	Free Fluorescein F/P (molar) Active Conjugate
Source	Isolated from <i>Griffonia (Bandeiraea) simplicifolia</i> seeds.
Specificity	Binds to α-(1,3) galactosyl galactoside epitopes.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com