

# Isolectin B<sub>4</sub> (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-B<sub>4</sub>, BSI-AB<sub>3</sub>, BSI-A<sub>2</sub>B<sub>2</sub>, BSI-A<sub>3</sub>B and BSI-A<sub>4</sub>. BSI-B<sub>4</sub> is blood group B specific and has an exclusive affinity for terminal α-D-galactosyl residues, whereas BSI-A<sub>4</sub> 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 (Griffonia simplicifolia), 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 <sup>2+</sup> .
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.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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