

Isolectin B₄ (Bandeiraea simplicifolia), (biotin conjugate)

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-650-001B-MC05	Biotin - 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
Formulation	Liquid. In 10mM HEPES, pH 7.5, containing 0.15M sodium chloride, 0.08% sodium azide and 0.1mM Ca ²⁺ .
Quality Control	Active Conjugate Lectin Biotinylated
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](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)