

# IgE (non-immune) (human), (biotin conjugate)

IgE comes from a monoclonal cell line. This antibody is well-suited as a standard in IgE-quantifying assays due to its very low batch-to-batch variation. Other applications include immunochemistry and cellular immunology research.

Citations: 13

[View Online »](#)

## Ordering Information

[Order Online »](#)

BPD-DIA-HE1B-015	150µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Do not freeze.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Immunoglobulin E
Application Notes	Can be used as standard in IgE quantifying assays. Functional Application: cell stimulation.
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.
Purity	≥90% (SDS-PAGE)
Purity Detail	Protein L purified. The original antibody-producing B cells were obtained from a healthy donor tested negative by US-FDA approved tests against HIV, HCV and Hepatitis B. Although the antibody comprising BPD-DIA-HE1B is purified from a monoclonal hybridoma line derived from the antibody-producing donor cell, it must be handled as potentially infectious as all human material. As the IgE comes from a monoclonal cell line, there is no contamination of antibodies of other isotypes. The remaining contaminants are mainly components from fetal bovine serum.
Source	Purified human IgE with kappa light chains produced <i>in vitro</i> from a monoclonal hybridoma.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.943.0420  
[usa@enzolifesciences.com](mailto: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-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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