

IgE (non-immune, azide) (human)

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: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

BPD-DIA-HE1A-1	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°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 0.15M sodium chloride and 0.09% sodium azide.
Purity	≥98% (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-HE1A 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.

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-be@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