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
BPD-DIA-HE1A-04	400µg

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 NotesCan 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 *in*

vitro from a monoclonal hybridoma.

Last modified: May 29, 2024

