



Immunoglobulin E (IgE)

Human Myeloma

Replace the undependable supply of human sourced materials with the consistency of our cell culture alternative.

Bio-Rad Critical Raw Materials IgE antigen is readily available, consistent, and free from concern of viral contamination. It is from a human Myeloma cell line, ensuring an analogue to the more erratic and uncertain material purified from human fluids.

	Method Used	Acceptable Result
Protein Quantitation	cobas IgE Assay	> 5.0 mg/mL
Product Quality	Centaur IgE Assay	> 2,083,330 IU/mL
Total Protein	OD280	< 75 O. D. Units

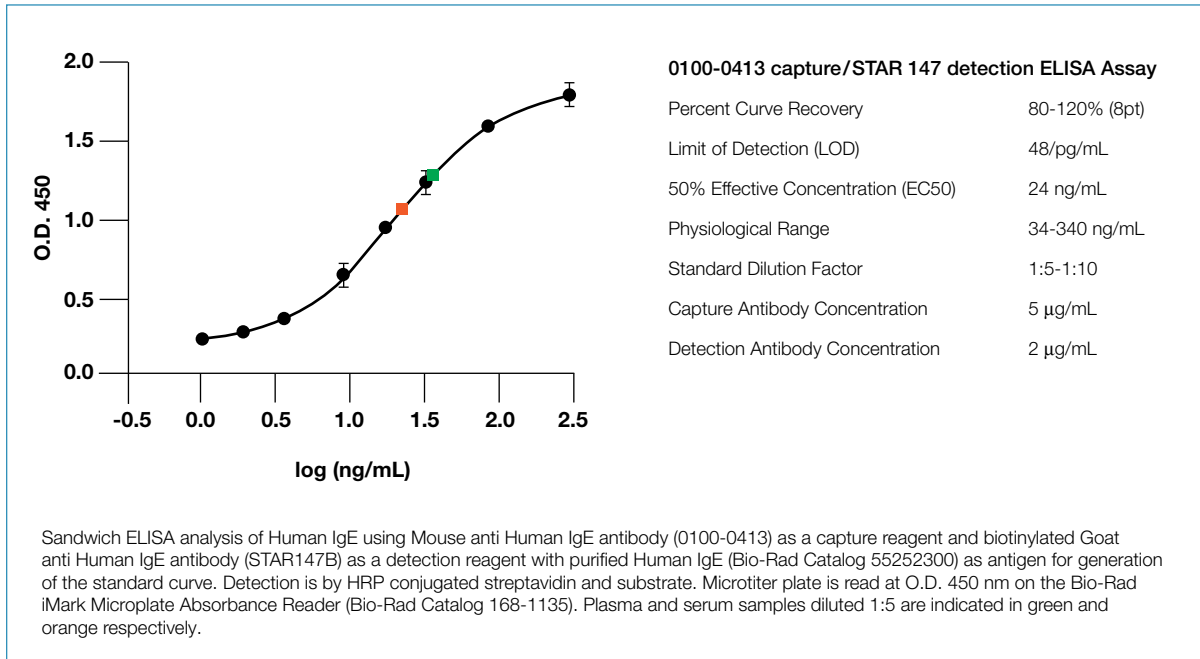
- Grade: Partially Purified
- Source: Human IgE Myeloma Cell Line
- Form: Liquid in 0.1M Tris/0.2M NaCl pH 7.5 ± 0.1
- Storage: -20°C
- Shelf Life: 4 years at -20°C

Immunoglobulin E (MW: 150Kd-900Kd) is a gamma-globulin that functions as part of the immune system in protecting the body against infections. They occur in humans as five different classes including IgG, IgM, IgA, IgE, and IgD and are synthesized by the B lymphocytes. Diseases resulting from deficiencies of the immunoglobulins typically result in recurrent infections.



This partially purified product has comparable commutability results to native IgE material on major diagnostic testing systems and is recommended for use in multi analyte controls, or in single analyte standards or controls.

The graph and information below demonstrate evidence of the performance of the IgE Ag., CC, PP 55252300 within a Sandwich ELISA.



Ordering Information

Cat #	Source	Grade
IgE Ag., CC PP		
55252300	Human Cell Culture	Partially Purified



**Bio-Rad
Laboratories**

For further information or to request a sample, please contact
CriticalRM_info@bio-rad.com or learn more online at www.bio-rad.com/crm

Clinical
Diagnostics Group

Website www.bio-rad.com/crm United States Bio-Rad Laboratories, 1045 Riverside St, Portland, Maine 04103 • Phone 207-615-0571 • Telefax 207-553-2523