

Bio-Rad Laboratories
The Clear Choice in Autoimmune Testing

Anti-rTPO

Autoimmune EIA

Simple, Objective, Sensitive, Specific

Bio-Rad offers the most comprehensive menu of EIA assays available today, with a constantly expanding menu that includes new clinically relevant assays.

Using recombinant human TPO antigen, the Bio-Rad Kallestad™ Anti-Thyroid Peroxidase (anti-TPO) test is a quantitative/qualitative enzyme-linked immunosorbent assay (ELISA) for the detection of the IgG class of autoantibodies specific for thyroid peroxidase in human serum or EDTA, heparin or citrated plasma. It is intended to aid in the multicriterion diagnostic process of autoimmune thyroid disorders.

Product Highlights

- Common assay protocol and reagents
- Color-coded breakaway wells
- Quantitative and qualitative result interpretations
- Easy-to-automate

Clinical Utility

Autoimmune thyroid disorders encompass autoimmune destruction and stimulation; both states are associated with predominantly IgG local and circulating thyroid autoantibodies.

Trotter et al¹ and Roitt et al² subsequently showed that many patients with advanced thyroiditis had antibodies for a thyroid antigen distinct from Tg (Thyroglobulin). This was termed thyroid microsomal antigen (TMA). Considerable evidence indicates that TMA is antigenically related to thyroid peroxidase (TPO), a membrane-bound glycoprotein enzyme with an approximate mass of 101kD, whose in vivo function is the iodination of tyrosine in the synthesis of the thyroid hormones T3 and T4. Anti-TPO antibodies are found, often in conjunction with anti-thyroglobulin autoantibodies, in the majority of Hashimoto's Thyroiditis, Graves' disease and in cases of Primary Myxoedema. The relationship of autoimmune thyroid disease

BIO-RAD

Continued on the next page >



Anti-Thyroid Peroxidase (rTPO)

in pregnancy has been the subject of considerable interest, with the demonstration of TPO antibodies in most cases of post-partum thyroid disease syndrome³⁻⁶ and the association of thyroid autoantibodies with increased miscarriage risk.⁷ Anti-TPO antibodies are found in other non-thyroid conditions, e.g. pernicious anemia,^{8,9} diabetes mellitus,^{10,11} rheumatoid arthritis,¹² Addison's disease¹¹ and Sjögren's Syndrome.⁹ In addition, anti-TPO antibodies are detectable at low levels in 2-8% of apparently healthy individuals, particularly in the elderly and more often in women than in men, although the clinical significance of this is unclear.

Bio-Rad's new recombinant kit format maintains the same Bio-Rad Kallestad™ EIA methodology, calibration to NIBSC 66/387 thyroid microsomal antibody reference preparation, and kit assay protocol (60'/30'/30' incubation) as the rest of the Bio-Rad Kallestad™ autoimmune EIA product range.

The driving force behind this improvement is to make an assay that provides reproducible results through the use of DNA recombinant technology. This technology minimizes the potential of antisera to bind both to specific and unspecific thyroid cell components. The new and improved Bio-Rad Kallestad™ Anti-Thyroid Peroxidase kit is made with human recombinant TPO antigen, which has been demonstrated to be highly specific for TPO autoantibodies. Additional features include: a lower clinical cut-off reflecting increased sensitivity and specificity and a narrower measuring range ensuring greater accuracy and precision.

Result Interpretation	
Qualitative Absorbance Ratio	Result Interpretation
< 1.0	Negative
≥ 1.0	Positive
Quantitative	
< 10 IU/mL	Negative
≥ 10 IU/mL	Positive

REFERENCES

1. Trotter WR, et al. *Proc R Soc Med*, 50, 961, 1957.
2. Roitt IM, et al. *Lancet II*, 1027-1033, 1958.
3. Vargas VT, et al. *J Clin Endocrin Metab*, 67, 327-333, 1988.
4. Jansson R, et al. *J Clin Endocrin Metab*, 58, 681-687, 1984.
5. Stagnaro-Green A. *Thyroid Today*, 16, 1-11, 1993.
6. Lazurus JH. *Thyroid*, 9 No. 7, 685-689, 1999.
7. Stagnaro-Green A, et al. *Jama*, 264 No.11, 1422-1425, 1990.
8. Estienne V, et al. *European Journal of Endocrinol*, 141, 563-569, 1999.
9. Feldt-Rasmussen, et al. *Autoimmunity*, 9, 245-253, 1991.
10. Chang CC, et al. *European Journal of Endocrinol*, 139, 44-48, 1998.
11. Scherbaum WA. *Acta Endocrinol (Copenh) Suppl*, 281, 325-329, 1987.
12. Walker DJ, et al. *Ann Rheum Dis*, 45, 323-326, 1986.

Ordering Information

Catalog No.	Description
25087	Bio-Rad Kallestad™ Anti-Thyroid Peroxidase (anti-rTPO)96 tests



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com.

Clinical Diagnostics Group

Website www.bio-rad.com U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-2507-6191 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-331-99-99 Portugal 351-21-472-7700 Russia 7-095-721-14-00 Singapore 65-6415-3188 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Thailand 662-651-8311 United Kingdom 44-208-328-2000