



Bio-Rad Laboratories
The Clear Choice in Autoimmune Testing



Autoimmune

Education and Training Tools

From the Worldwide Leader in Autoimmune Testing

As the worldwide leader in autoimmune diagnostics, Bio-Rad offers premier customer support and technical service both in product offering and ongoing customer education and support. This leadership position is further enhanced with the acquisition of both Sanofi Diagnostics Pasteur and Helix™ Diagnostics along with the well-established product branding under the Kallestad™ name.

Bio-Rad offers the most comprehensive customer education program. From workstation aids to sophisticated web-based training software, Bio-Rad understands how critical it is to obtain and maintain proficiency in this complex area of testing, and is committed to facilitating access to continuing education programs and day-to-day technical support.

Choose and stay with Bio-Rad autoimmune products because of the strength of our product line in quality and breadth of menu and our superior ability to train and support your staff.

Educational Tools and Programs

- AutoimmuneWEB™
- AutoimmuneTUTOR™
- Autoimmune Atlas™
- Pattern Plus Auditor™
- Patient Sample Pattern Interpretation Services
- InServices
- Worldwide seminars and workshops
- Extensive technical support and sales network

A
U
T
O
I
M
M
U
N
E
T
E
S
T
I
N
G



BIO-RAD



Education and Training Tools

AutoimmuneWEB™

The only comprehensive web-based product for autoimmune testing. AutoimmuneWEB™ is the premier service for online, interactive autoimmune customer education that offers an innovative, state-of-the-art approach for training, optimizing expertise, and competency assessment.

- Interactive – AutoimmuneWEB™ uses video and digital animation to keep you engaged in the learning process.
- Dynamic – Updated competency exams, relevant links to autoimmune websites, and expansion of IFA pattern images and FAQs on a continual basis.
- Simple – Provides the flexibility of self-paced learning and reduces the time and costs associated with off-site training seminars.

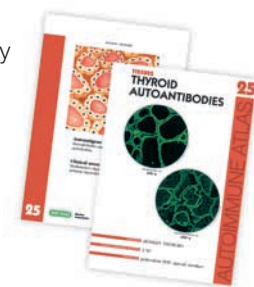
AutoimmuneTUTOR™

AutoimmuneTUTOR™ is an interactive computer program that enhances the understanding of autoantibody testing. It helps the laboratory comply with licensing requirements for initial instruction, continuing education (accredited for 5 contact hours with ASCLS P.A.C.E.® program) and competency assessment.



Autoimmune Atlas™*

Autoimmune Atlas™ is a unique collection of fluorescent antibody patterns and histologic illustrations. Fifty autoantibody panels covering typical and unusual HEp-2 staining patterns, twenty-five autoantibody panels covering tissue staining patterns on rodent and monkey substrates are included.



Pattern Plus Auditor™

Pattern Plus Auditor™ provides a quarterly mailing of four high quality prints of immunofluorescent images. Assemble these prints to create your own reference binder. Information about the image includes how to identify the pattern, the antibody and the clinical correlation.



Ordering Information

Catalog No. Description

AutoimmuneWEB™

465-0001	AutoimmuneWEB™, English Version	1 year subscription (per site)
465-0004	AutoimmuneWEB™, French Version	1 year subscription (per site)
465-0007	AutoimmuneWEB™, Spanish Version	1 year subscription (per site)

(2 and 3 year subscriptions of AutoimmuneWEB™ are also available)

AutoimmuneTUTOR™

425-1150	AutoimmuneTUTOR™ CD Ver. 1.0, English/Unlimited Version	1 unit
425-1151	AutoimmuneTUTOR™ CD Ver. 1.0, French/Spanish/Unlimited Version	1 unit
425-1156	AutoimmuneTUTOR™ CD Ver. 1.0, English/15-day Demo	1 unit

Autoimmune Atlas™*

17521	Autoimmune Atlas™*	1 unit
-------	--------------------------	--------

Pattern Plus Auditor™

102799	Pattern Plus Auditor™	4 patterns per quarter
--------	-----------------------------	------------------------

* Not available in the US.



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 5521-3461-5202 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