



Liquichek™ Autoimmune Controls



Liquichek™ Autoimmune Controls

Individual controls for monitoring the performance of autoimmune assays using indirect immunofluorescence (IFA)

- Liquid, human serum based
- 2 year shelf life at 2–8°C
- 6 month open-vial stability at 2–8°C

Package Configuration

- Negative and weak positive controls packaged separately
- 0.5 mL per vial

Refer to the package insert of currently available lots for specific analyte and stability claims



Ordering Information

Cat #	Liquichek™ Autoimmune Controls	Quantity
107	Liquichek™ ANA Control, Homogeneous Pattern	.3 x 0.5 mL
108	Liquichek™ ANA Control, Speckled Pattern	.3 x 0.5 mL
111	Liquichek™ ANA Control, Nucleolar Pattern	.3 x 0.5 mL
112	Liquichek™ ANA Control, Centromere Pattern	.3 x 0.5 mL
118	Liquichek™ ANA Control, Mitotic Spindle Pattern	.2 x 0.5 mL
131	Liquichek™ ANA Controls Set, Positive (Homogeneous, Speckled, Centromere and Nucleolar Patterns)	.4 x 0.5 mL
113	Liquichek™ Anti-SS-B Control	.2 x 0.5 mL
114	Liquichek™ Anti-SS-A Control	.2 x 0.5 mL
115	Liquichek™ Anti-Sm Control	.2 x 0.5 mL
116	Liquichek™ Anti-RNP Control	.2 x 0.5 mL
117	Liquichek™ Anti-Scl-70 Control	.2 x 0.5 mL
119	Liquichek™ Anti-nDNA Control	.3 x 0.5 mL
127	Liquichek™ Anti-Mitochondrial Control	.3 x 0.5 mL
129	Liquichek™ Anti-Smooth Muscle Control	.3 x 0.5 mL
130	Liquichek™ Autoimmune Negative Control	.3 x 0.5 mL



These controls are featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

Discover the power of Unity™ Solutions at www.QCNet.com.



**Bio-Rad
Laboratories**

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

**Clinical
Diagnostics Group**

Website www.bio-rad.com/qualitycontrol **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400 **Canada** 1-514-334-4372 **China** 86-21-64260808 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00 **Germany** +49-(0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 1-800-180-1224 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** +52 (55) 5488-7670 **The Netherlands** +31-318-540666 **New Zealand** 64-9-415-2280 **Norway** 47-23-38-41-30 **Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** 7-495-721-14-04 **Singapore** 65-6415-3170 **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-(0)20-8328-2000