



Whole Blood Immunosuppressant Controls



Liquichek™ Whole Blood Immunosuppressant Control

A **liquid**, human blood based control to monitor immunosuppressant drugs including Everolimus. Suitable for immunoassay and chromatographic methods.

- Four individual levels to provide labs with choice and flexibility
- 3 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2°C – 8°C for most analytes

Lyphochek® Whole Blood Immunosuppressant Control

A unique **lyophilized** control with the flexibility of five levels to monitor three widely prescribed immunosuppressant drugs across a wide range of popular test procedures.

- Five individual levels that range from low to very high concentrations
- 3 year shelf life at 2°C – 8°C
- 14 day open-vial stability at 2°C – 8°C
- 30 day reconstituted, frozen stability at -20°C to -70°C

Analytes

Cyclosporine
Everolimus

Sirolimus
Tacrolimus

Analytes

Cyclosporine
Sirolimus

Tacrolimus

Catalog Number	Liquichek™ Whole Blood Immunosuppressant Control				Lyphochek® Whole Blood Immunosuppressant Control				
	Level 1	Level 2	Level 3	Level 4	Level 1	Level 2	Level 3	Level 4	Level 5
	12000404	12000405	12000406	12000407	274	275	276	277	278
Everolimus by Chromatology		●	●	●					
Typical range methods for Cyclosporine, Sirolimus and Tacrolimus	●	●	●		●	●	●		
Extended range methods for Tacrolimus and Cyclosporine, including C2 monitoring				●				●	●

Refer to the package insert for instrument-specific ranges for each level.

Ordering Information

Liquichek™ Whole Blood Immunosuppressant Control

Cat #	Quantity
12000404 Level 1	6 x 2 mL
12000405 Level 2	6 x 2 mL
12000406 Level 3	6 x 2 mL
12000407 Level 4	6 x 2 mL
12000408 4 Level MiniPak	4 x 2 mL

Lyphochek® Whole Blood Immunosuppressant Control

Cat #	Quantity
274 Level 1	6 x 2 mL
275 Level 2	6 x 2 mL
276 Level 3	6 x 2 mL
277 Level 4	6 x 2 mL
278 Level 5	6 x 2 mL
279X 5 Level MiniPak	5 x 2 mL

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



Bio-Rad
Laboratories

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

Clinical
Diagnostics Group

Website www.bio-rad.com/qualitycontrol Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil +55 (31)3689-6600 Canada 1-514-334-4372 China 86-21-61698500 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 1800-180-1224 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 +7-495-721-1404 Singapore 65-6415-3170 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden +46-8-555-127-00 Switzerland +41 (0)26-674-55-05/06 Taiwan 886-2-2578-7189 Thailand 662-651-8311 United Kingdom +44 (0)20-8328-2000