

BioPlex 2200 Syphilis Total & RPR Assay

Simultaneous detection and differentiation of Treponemal and non-Treponemal antibodies

The BioPlex Syphilis Total & RPR assay is the only dual treponemal/non-treponemal immunoassay test that offers laboratories a global one-step testing solution for Syphilis screening. The BioPlex Syphilis Total & RPR kit is intended for the qualitative detection of Total (IgG/IgM) antibodies to *Treponema pallidum* and the qualitative detection and/or titer determination of non-treponemal reagin antibodies in human serum or plasma.

- Syphilis Total (IgG/IgM)
- RPR (Rapid Plasma Reagin)

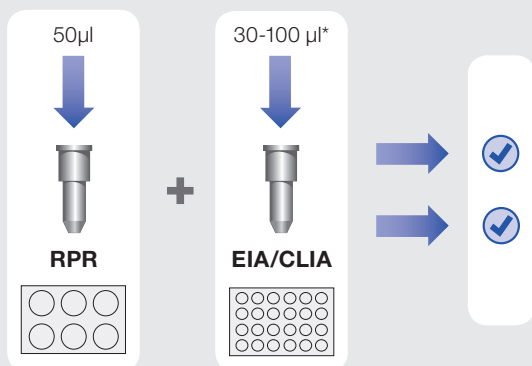


Simplify your Workflow

Traditional or Reverse Algorithm

Screen with a reagent test or Treponemal assay (EIA/CLIA). Confirmation by an orthogonal test.

- 2 sample aliquots
- 2 test results



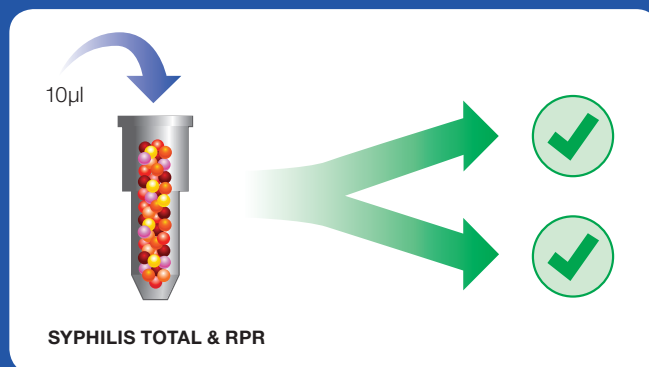
2 tests, 2 methods = 2 results

*method dependent

BioPlex Test

Tests for both Syphilis Total & RPR, with the flexibility to order only one and/or both analytes.

- 1 sample aliquot
- 2 test results/sample



SYPHILIS TOTAL & RPR

1 test = 2 results

Performance

- Excellent negative agreement with commercially available methods for both Syphilis Total (98.9%) and RPR (99.5%)
- Excellent positive agreement with commercially available methods for medically diagnosed samples for Syphilis Total (99.1%) and RPR (100%)
- Excellent total precision for Syphilis Total ($\leq 11.2\%$) and RPR ($\leq 13.4\%$)

Improved Laboratory Workflow

- Simultaneous detection of Syphilis Total & RPR for higher total throughput, time and labor savings and improved patient management
- Fully automated RPR assay reduces hands-on time with accurate and objective result reporting and LIS traceability
- On-board RPR titer/dilutions (1:4, 1:8, 1:16, 1:32, 1:64)
- Time to first result in 45 minutes
- Up to 200 results per hour; up to 2 different reportable test results from 1 patient sample
- Ready-to-use calibrators and controls
- Calibrator curve stability up to 30 days

Ordering Information

Catalog #	Description	
12000650	BioPlex 2200 Syphilis Total & RPR Reagent Pack, 100 samples.....	1 pack
12000651	BioPlex 2200 Syphilis Total & RPR Calibrator Set.....	1 set
12000653	BioPlex 2200 Syphilis Total & RPR Control Set	2 sets

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

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

Clinical
Diagnostics Group

Web site www.bio-rad.com/diagnostics USA 1 800 224 6723 Australia 61 2 9914 2800 Austria 43 1 877 8901 Belgium 32 03 710 53 00
 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