



# Geenius Solution

A Game Changer



# Geenius HIV 1/2 Confirmatory

The Bio-Rad Geenius HIV 1/2 Confirmatory Assay is a single-use immunochromatographic test for the confirmation and differentiation of individual antibodies to Human Immunodeficiency Virus Types 1 and 2 (HIV-1 and HIV-2) in fingerstick whole blood, venous whole blood, serum or plasma samples.



HIV-1 & HIV-2 confirmation and differentiation



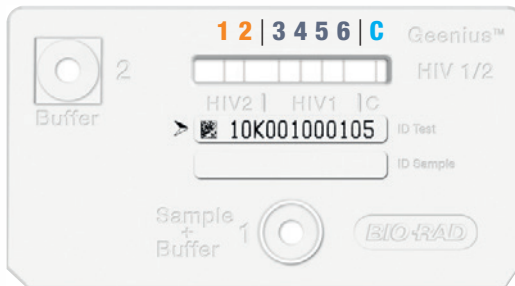
< 30 minutes



Room temperature storage



Automatic reading



- 1 Band 1** - gp36 (HIV-2, Envelope peptide) HIV-2 ENV
- 2 Band 2** - gp140 (HIV-2, Envelope peptide) HIV-2 ENV
- 3 Band 3** - p31 (HIV-1, Polymerase peptide) HIV-1 POL
- 4 Band 4** - gp160 (HIV-1, Envelope recombinant protein) HIV-1 ENV
- 5 Band 5** - p24 (HIV-1, Core recombinant protein) HIV-1 GAG
- 6 Band 6** - gp41 (Group M and O) HIV-1 ENV (HIV-1, Envelope peptide)
- C Control Band** - Protein A



- Listed on the following regulatory certifications: CE, FDA, TGA, ANVISA, Health Canada, CDSCO India, Thai FDA, Taiwan FDA
- Listed on the WHO prequalified IVD products list



# Geenius HCV Supplemental

Geenius HCV Supplemental Assay is a qualitative single-use immunochromatographic test for the confirmation of the presence of antibodies associated with infection caused by the Hepatitis C Virus (HCV) in human venous whole blood, serum or plasma specimens from blood donors or hospitalized patients.



Unitary HCV Supplemental Assay



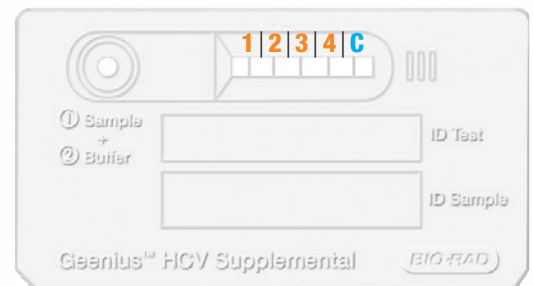
20 minutes



Room temperature storage



Automatic reading



- 1** Band 1 - NS4 antigen peptides
- 2** Band 2 - NS5 recombinant protein
- 3** Band 3 - Core antigen peptides
- 4** Band 4 - NS3 recombinant protein
- C** Control Band - Protein A



■ CE-marked

# Geenius Solution

A Game Changer

## One Patient. One Cartridge. One Result.



### Accurate Results

- Very high level of sensitivity and specificity
- Simple positivity criteria for easy overall interpretation
- Excellent differentiation rate



### Automatic Reading & Interpretation

- A tiny system with mighty possibilities, that can find its place easily in every laboratory
- Automated interpretation of each individual band
- Automatic conclusion of patient status using expert software and proprietary algorithms
- Intuitive and user friendly interface
- Minimal training required



### Full Traceability & Safety

- Unique datamatrix barcode on each cassette
- Automatic reader calibration
- Automatic checking of lot validation
- Positive identification of controls and patients
- Detailed report including cassette picture
- Secure data archiving
- Bidirectional interfacing with LIS



Geenius Controls



Geenius Assays

Visit [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) for more information.







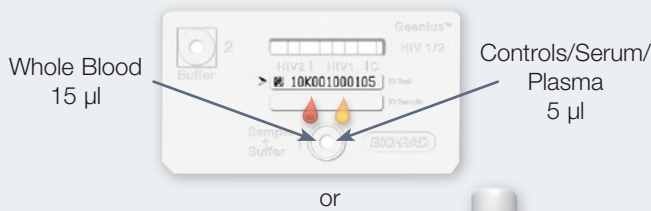
# Key Bio-Rad Milestones for Innovation in HIV and HCV Detection



Advance Discovery. Improve Lives.

# Protocol Geenius HIV 1/2 Confirmatory

**1** Dispense Controls or Samples into well 1



**2** Buffer  
2 Drops into well 1



Wait 5-7 minutes



**3** Buffer  
5 Drops into well 2



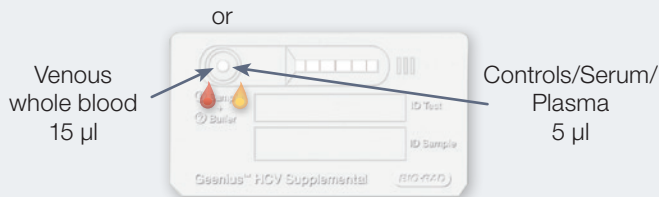
Wait 20 minutes

Read, interpret and report results



# Protocol Geenius HCV Supplemental

**1** Dispense Controls or Sample in "Sample + Buffer" Well



**2** Buffer  
4 Drops into "Sample + Buffer" Well



20 minutes

Read, interpret and report results



## Assays Ordering Information

Catalog #	Description
72460	Geenius HIV 1/2 Confirmatory Assay, 20 tests
72329	Geenius HIV 1/2 Confirmatory Controls (1 Positive + 1 Negative), 2 x 120 µl
92501	Geenius HCV Supplemental Assay, 20 tests
92502	Geenius HCV Supplemental Controls (1 Positive + 1 Negative), 2 x 150 µl

## System Ordering Information

Catalog #	Description
92465	Geenius Reader, 1 unit
12017081	Geenius PC + SW OUS (Dell 3510 Win10), 1 unit
92461	Geenius Software, 1 unit
94197	Anti-theft System, 1 unit

BIO-RAD and BIOPLEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.  
 GEENIUS, MONOLISA and DECISCAN are trademarks of Bio-Rad Europe, GmbH in certain jurisdictions.  
 All trademarks used herein are the property of their respective owner.

**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](http://www.bio-rad.com/diagnostics)*

Clinical  
Diagnostics Group

Website [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) U.S. 1-800-224-6723 Australia +61(2)9914-2800 Austria +43-1-877-8901 Belgium +32(3)710-53-00 Brazil +55(31)3689-6600  
 Canada 1-514-334-4372 China +86 21 6169 8500 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