



## Geenius HCV Supplemental Assay

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

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



# Protocol

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

or

15  $\mu$ l venous whole blood

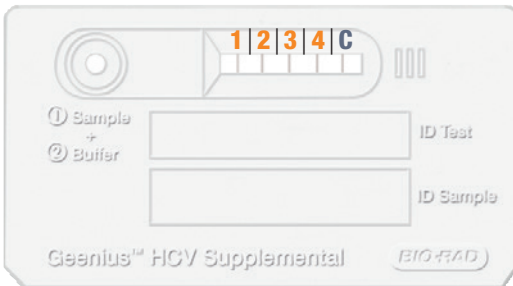
5  $\mu$ l control/serum/plasma

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

20 minutes

Read, interpret and report results

# Results Interpretation



- 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

## Ordering Information

Catalog #	Description		
92501	<b>Geenius HCV Supplemental Assay</b> , 20 tests	92465	<b>Geenius Reader</b> , 1 unit
92502	<b>Geenius HCV Supplemental Controls</b> (1 Positive + 1 Negative), 2 x 150 $\mu$ l	12017081	<b>Geenius PC + SW OUS (Dell 3510 Win10)</b> , 1 unit
		92461	<b>Geenius Software</b> , 1 unit
		94197	<b>Anti-theft System</b> , 1 unit

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.  
 GEENIUS is a trademark of Bio-Rad Europe, GmbH in certain jurisdictions.



Clinical  
Diagnostics Group

**Website** [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) **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** +86-21-61698500 **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** +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 26 674 55 05/06 **Taiwan** +886-2-2578-7189 **Thailand** +662-651-8311 **United Kingdom** +44(0)20-8328-2000

