



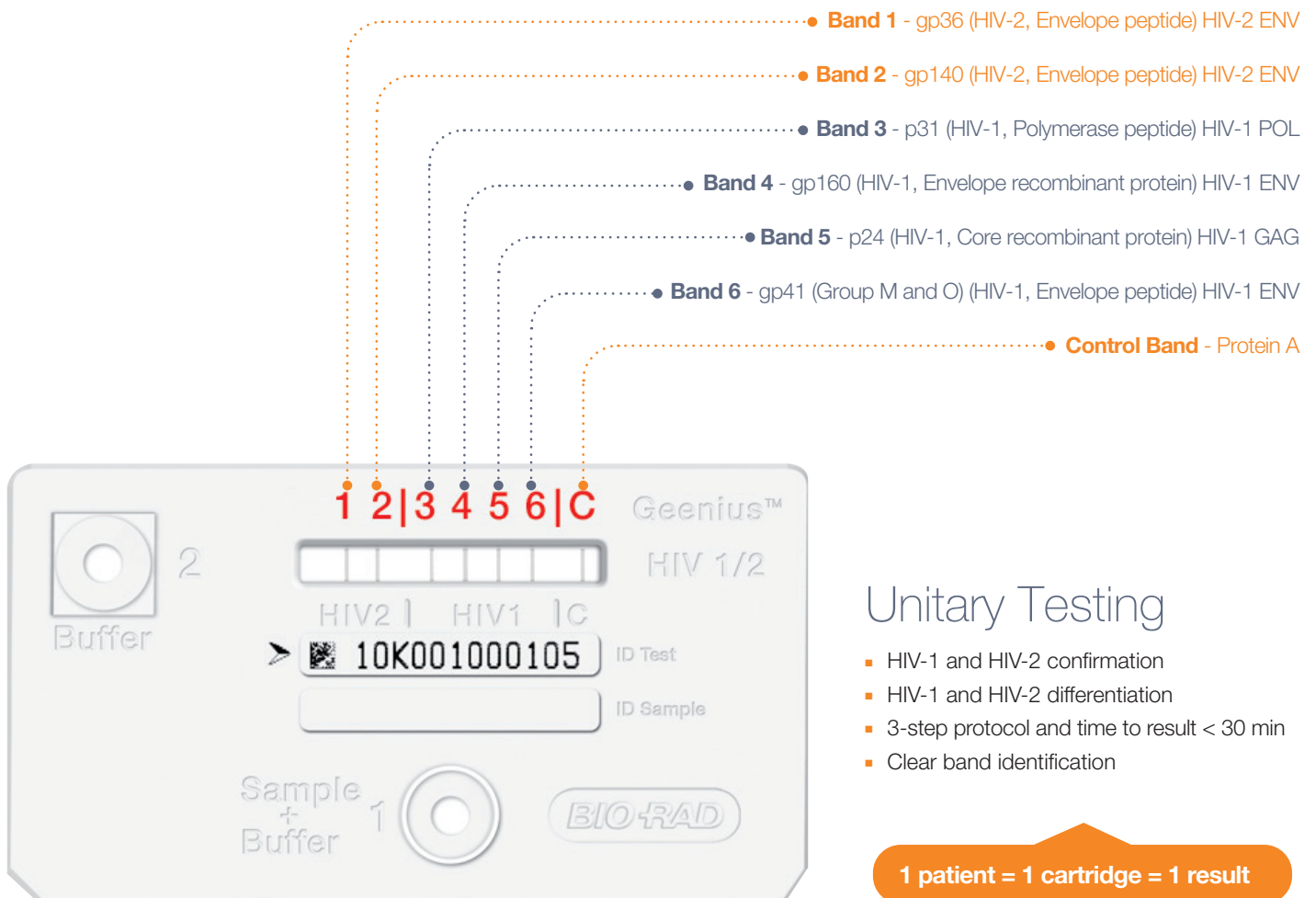
Geenius HIV 1/2 Confirmatory System

A Game Changer



Geenius HIV 1/2 Confirmatory System

A Game Changer



Accurate Results

- High specificity: very low rate of indeterminate results
- High sensitivity performance
- Simple positivity criteria for easy overall
- Interpretation
- Excellent differentiation rate



Geenius Reader



Notebook
(validated for Geenius Software)



Geenius Software

Geenius Report

Automatic Reading and Interpretation

- Objective reading
- Automated interpretation of each individual band
- Qualitative interpretation of individual HIV-1 and HIV-2 assays
- Automatic conclusion using expert software and proprietary algorithms
- Intuitive and user friendly interface
- Minimal training required interpretation

Full Traceability and Safety

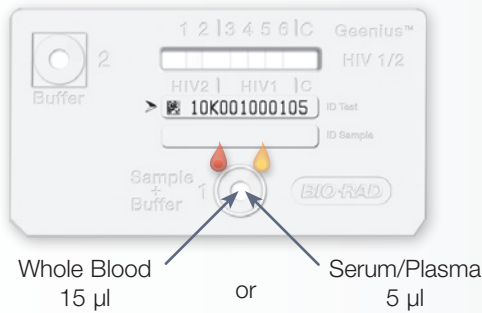
- Unique Datamatrix barcode on each cassette
- Automatic reader calibration
- Automatic checking of lot validation
- Positive identification of controls and patients
- Detailed report including cassette picture
- Secure Data archiving
- Bidirectional interfacing with LIS



Geenius HIV 1/2 Confirmatory Assay

Protocol

- 1 Dispense 15 µl of whole blood or 5 µl of serum/plasma into well 1



- 2 Buffer
2 Drops into well 1



Wait 5-7 minutes
All the blue colored test lines should have disappeared from the **Test** and **Control** window



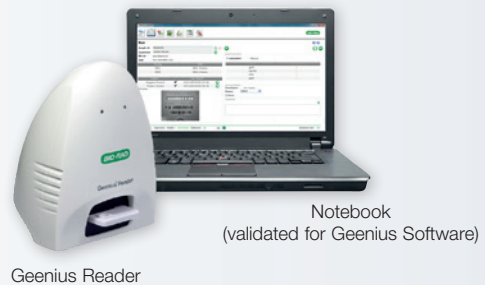
- 3 Buffer
5 Drops into well 2



Wait 20 minutes



- 4 Read, interpret and report results
Do not read results later than 30 minutes after the addition of the buffer into well 2



Ordering Information

Catalog #	Description
72460	Genius HIV 1/2 Confirmatory Assay, 20 tests
72329	Genius HIV 1/2 Confirmatory Controls (1 Positive + 1 Negative), 2 x 120 µl
92465	Genius Reader, 1 unit
12017081	Genius PC + SW OUS (Dell 3510 Win10), 1 unit
92461	Genius Software, 1 unit
94197	Anti-theft System, 1 unit



Visit www.bio-rad.com/diagnostics for more information.



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



**Bio-Rad
Laboratories**

Clinical
Diagnostics Group

Website 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 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