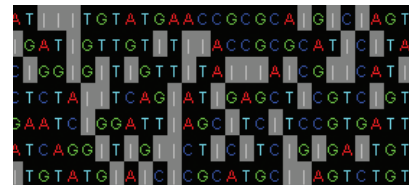


**FPQuest™ and InfoQuest™ FP Bioinformatics Software**



**Comprehensive Solutions for the Study of Biological Relationships**



## Bio-Rad Bioinformatics Software

FPQuest and InfoQuestFP modular software packages offer customizable applications to meet a variety of laboratory informatics requirements. FPQuest software offers advanced analysis and statistical tools for analysis of multiple fingerprints or one-dimensional (1-D) gels. Powerful InfoQuestFP software combines the functionality of FPQuest with the ability to analyze many other data types, for more comprehensive studies of biological relationships. Use the selection guide (see last page) to customize the software for your needs. The modular software can accommodate changing needs for multiple types of sample data, analytical functionality, and display of analyses. For more information, go to [www.bio-rad.com/bioinformatics/](http://www.bio-rad.com/bioinformatics/)

## Integrate and Analyze Data From Multiple Bio-Rad Products\*

- BioLogic DuoFlow™ software
- iQ™5 real-time PCR detection system software
- Bio-Plex Manager™ 4.1 software
- Experion™ software

\* Contact Technical Support for more information.

# Solutions for Massive Amounts of Biological Information

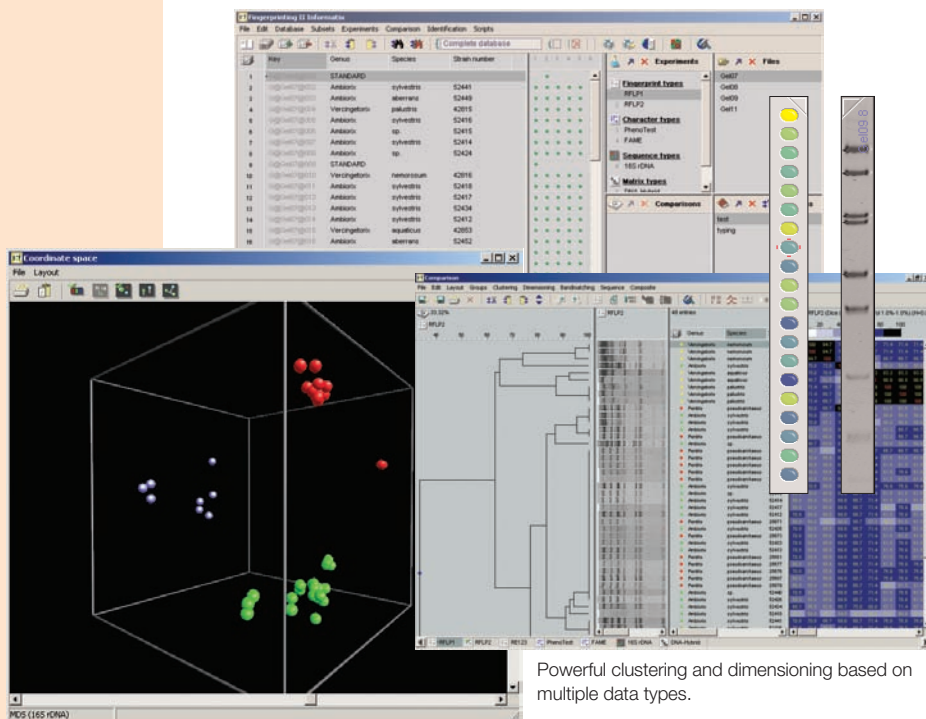
## Analysis of All Types of Data

InfoQuestFP and FPQuest bioinformatics software are universal tools for studying biological relationships. The programs include a relational database that combines all types of biological data, including phenotypic, genotypic, electrophoretic, array, and sequence information. Whatever your organism or sample — microbial strains, viruses, animals, plants, or tissues — you can collect, compare, and analyze data using the software's powerful clustering and statistical tools.

## Modularity

The modular software packages expand with your growing needs for multiple types of sample data, analytical functionality, and display options. Modules that accommodate various sample data types allow tailoring of InfoQuestFP for the sample types that best fit your experimental data. Optional analysis modules customize the software to your data analysis needs. Whether your bioinformatics needs are for strain typing, identification of unknown organisms, or comparisons of large numbers of genes, InfoQuestFP offers the algorithms and statistical features you need.

The software recognizes an extensive set of file formats, including text, EMBL, GenBank, FASTA, and Excel, and a variety of image formats.

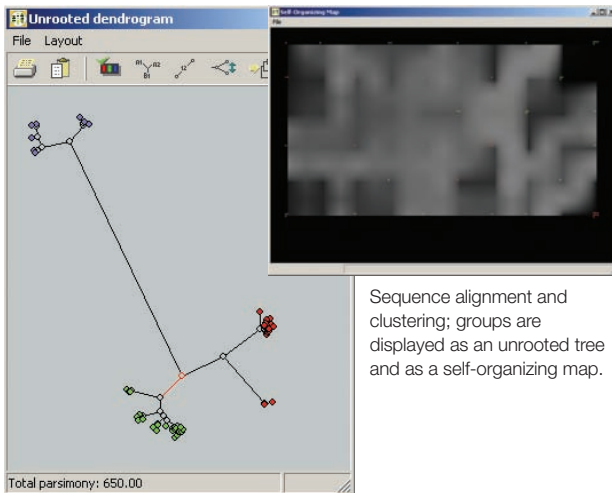
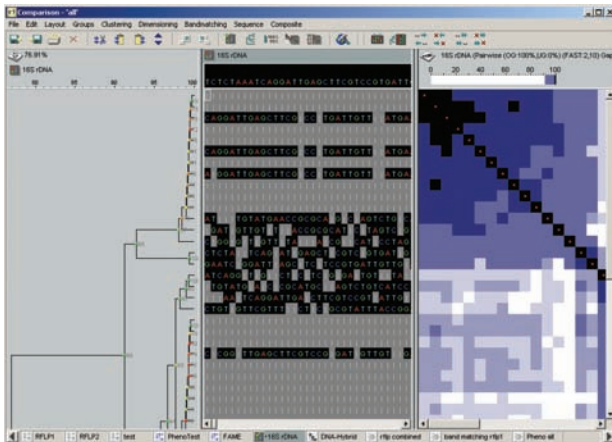


Powerful clustering and dimensioning based on multiple data types.

## Data Sharing and Export

InfoQuestFP has powerful data export and database sharing capabilities that simplify information exchange. The optional database sharing module is an excellent fit for laboratories with a large number of users on a network, for research collaborations, or when access to a central public or private database is needed. Users can query the database, exchange data, and create automated reports or alerts. InfoQuestFP allows:

- Printing of reports and analyses
- Customized reporting using scripts
- Peer-to-peer data exchange or client-server projects
- Storing of the database as a connected database linked to Oracle, SQL Server, or Access
- Multiple network licensing on a local area network



## Fingerprint Types

- Electrophoretic gels
- HPLC and other chromatographic patterns
- Spectrophotometric and densitometric curves

## Character Types

- Binary or continuous arrays
- Microarrays and gene chips
- Enzyme activity test panels
- Fatty acid profiles

## Sequence Types

- Nucleic acids
- Amino acids

## Matrix Types

- External similarity and distance matrices

## Cluster Analysis

- Dendrograms from individual experiments and composite data sets
- Phylogenetic tree construction
- Cluster and group significance
- Congruence of techniques

## Identification and Library Manager

- Identification of unknowns against a library of entries of the same taxon or group
- Quick similarity searches
- Construction of libraries

## Dimensioning Techniques

- Principal component analysis
- Multidimensional scaling
- Discriminant analysis and MANOVA
- Self-organizing maps

## Database Sharing Tools

- Peer-to-peer data exchange over the Internet
- External database access
- Client-server projects

**Bioinformatics Software Selection Guide**

Software Analysis Features	Catalog # for Modules	
	FPQuest	InfoQuestFP
<b>Application Modules (Type of Data Handled)</b>		
1-D gels, chromatograms, density curves	Basic software (170-9300)*	Basic fingerprint types (170-9310)*
Microarrays, profiles, test panels, dot blots, etc.	—	Character types (170-9311)*
Sequence data; direct import of EMBL, GenBank, Flat A, and FASTA formats	—	Sequence types (170-9312)*
Similarity and distance matrices; partial (DNA homology) matrices accepted	—	Matrix types (170-9313)*
<b>Analysis Tools (Type of Analysis Performed)</b>		
Dendrogram generation and cluster analysis	Cluster analysis (170-9301)	Cluster analysis (170-9314)
Database screening and library generation	Identification and library manager (170-9302)	Identification and library manager (170-9315)
Comparative quantitation of bands between groups of patterns	Comparative quantitation and polymorphism analysis (170-9303)	Basic fingerprint types (170-9310); cluster analysis (170-9314)
Principal component analysis and nonhierarchical grouping by MDS	Dimensioning techniques (170-9304)	Dimensioning techniques (170-9316)
Data sharing over the Internet	Database sharing tools (170-9305)	Database sharing tools (170-9317)

\* At least one is required.

**Ordering Information**

Catalog # Description

**FPQuest Software**

170-9300 FPQuest Software

**Optional FPQuest Analysis Tools**

170-9301 FPQuest Cluster Analysis  
 170-9302 FPQuest Identification and Library Manager  
 170-9303 FPQuest Comparative Quantitation and Polymorphism Analysis  
 170-9304 FPQuest Dimensioning Techniques  
 170-9305 FPQuest Database Sharing Tools

**InfoQuestFP Software\***

170-9310 InfoQuestFP Basic Fingerprint Types Module  
 170-9311 InfoQuestFP Character Types Module  
 170-9312 InfoQuestFP Sequence Types Module  
 170-9313 InfoQuestFP Matrix Types Module

**Optional InfoQuestFP Analysis Tools**

170-9314 InfoQuestFP Cluster Analysis  
 170-9315 InfoQuestFP Identification and Library Manager  
 170-9316 InfoQuestFP Dimensioning Techniques  
 170-9317 InfoQuestFP Database Sharing Tools

\* At least one module is required.

Access, Excel, and SQL Server are trademarks of Microsoft Corporation. GenBank is a trademark of the US Dept. of Health and Human Services. Oracle is a trademark of Oracle Corporation.

Bio-Rad's iQ5 real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400  
 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050  
 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280  
 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723  
 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000