



License String Document -DEMO

Fingerprinting[®] II and Fingerprinting II Informatix[®] Software

This software is for demonstration purposes only. To obtain a license for the software, please send e-mail to LSG.TechServ.US@bio-rad.com with the SPRO-number found on the dongle in this package. You will be granted a License String valid for 50 hours of operation of the software. The License String can be renewed at no cost when 50 hours have expired.

Fingerprinting II Software

- 170-9300 Basic Software
- 170-9301 Cluster Analysis
- 170-9302 Identification and Library Manager
- 170-9303 Comparative Quantification and Polymorphism Analysis
- 170-9304 Dimensioning Techniques
- 170-9305 Database Sharing Tools

Fingerprinting II Informatix Software

- 170-9310 Fingerprint Types
- 170-9311 Character Types
- 170-9312 Sequence Types
- 170-9313 Matrix Types
- 170-9314 Cluster Analysis
- 170-9315 Identification and Library Manager
- 170-9316 Dimensioning Techniques
- 170-9317 Database Sharing Tools

➤ **SPRO Number:**

➤ **License String:**

(will be provided in e-mail)

➤ **Order number:**

(internal use only)