



HUMAN HEALTH

ENVIRONMENTAL HEALTH

PERKINELMER ARRAY SOFTWARE SOLUTIONS

GENOGLYPHIX SOFTWARE

PerkinElmer Array Software Solutions

Not for distribution in the USA



Genoglyphix software is for Research Use Only.
Not for use in diagnostic procedures.

GENOGLYPHIX: TURNING 50,000 SAMPLES INTO MEANING

Genoglyphix® is a powerful data visualization software and database based on the analysis of over 50 000 validated cytogenetic samples.

Developed at Signature Genomics by Cytogeneticists for Cytogeneticists, Genoglyphix offers a proven analysis solution with a complete workflow

including sample tracking, aberration categorization, final report creation functionality, and data sharing options with other centres. It also offers the possibility to create user-specific database tracks. Genoglyphix provides intuitive data visualization and annotation features for streamlined, rapid and clear interpretation of CGX data.

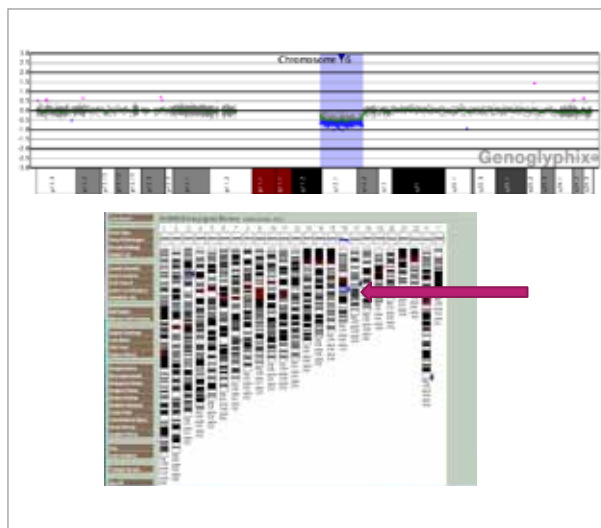
Genoglyphix software – bringing multiple genetic databases into one tool

- **Access to Signature's database** containing over 12,000 validated genetic alterations identified in over 50,000 samples
- **Direct links to other relevant resources** such as the DGV, OMIM, PubMed, UCSC and Ensembl databases for easy data interpretation
- **Secure web based (128-bit encryption)** access to Genoglyphix database and software

Genoglyphix guides cytogeneticists through the analysis with intuitive and easy-to use tools

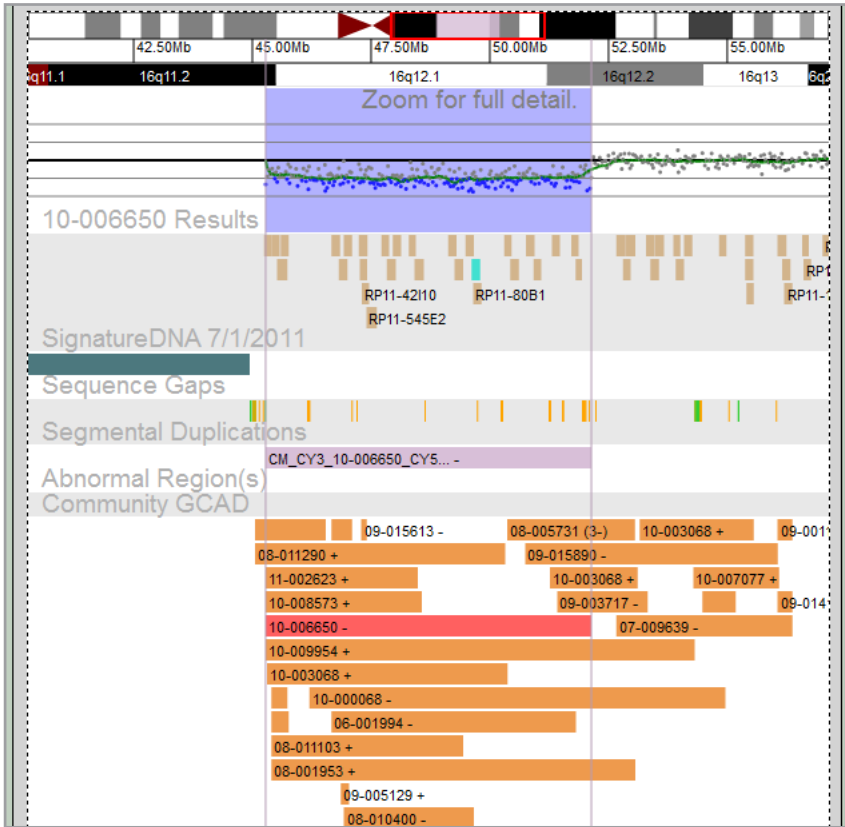
- Generation of user-defined databases and custom tracks displaying abnormal results, copy number variants and analysis notes
- Report creation functionality
- Optional sharing of data with other Genoglyphix users
- Availability of FISH probes for result confirmation

Figure 1: Case Review - Karyogram View



Data visualization in Genoglyphix in Chromosome and Karyogram View

Figure 2: Case Review - Genoglyphix Genome Browser

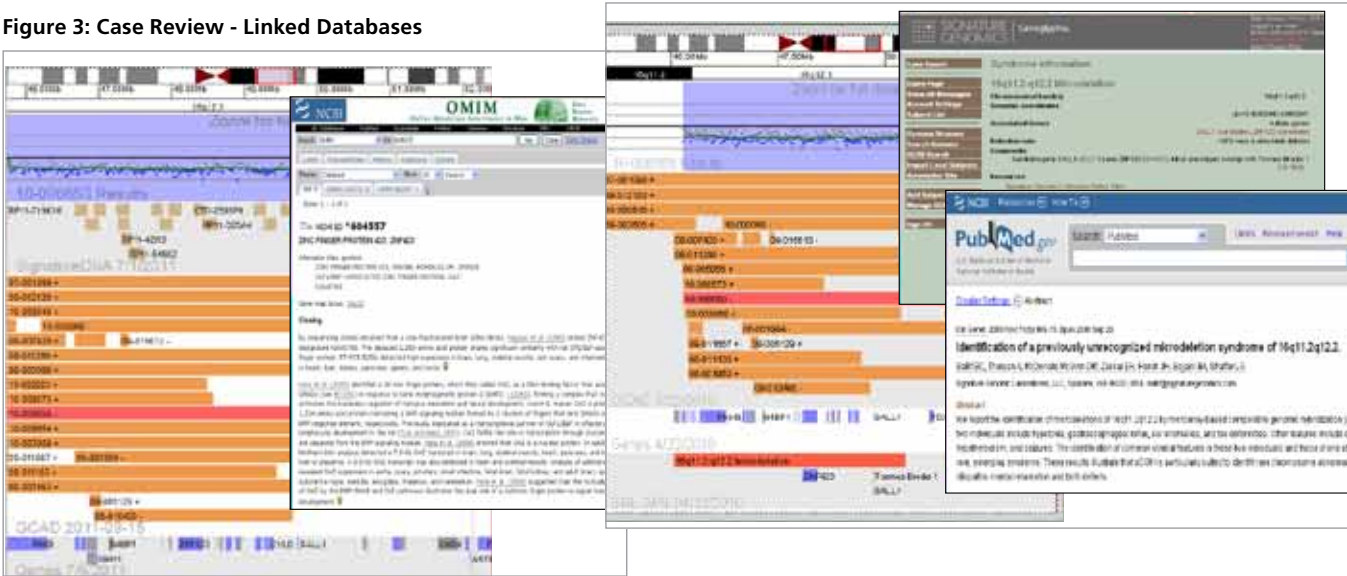


< Graphical representation of aberrations in the Genoglyphix Genome Browser

< Identify Signature BAC clones available for FISH confirmation

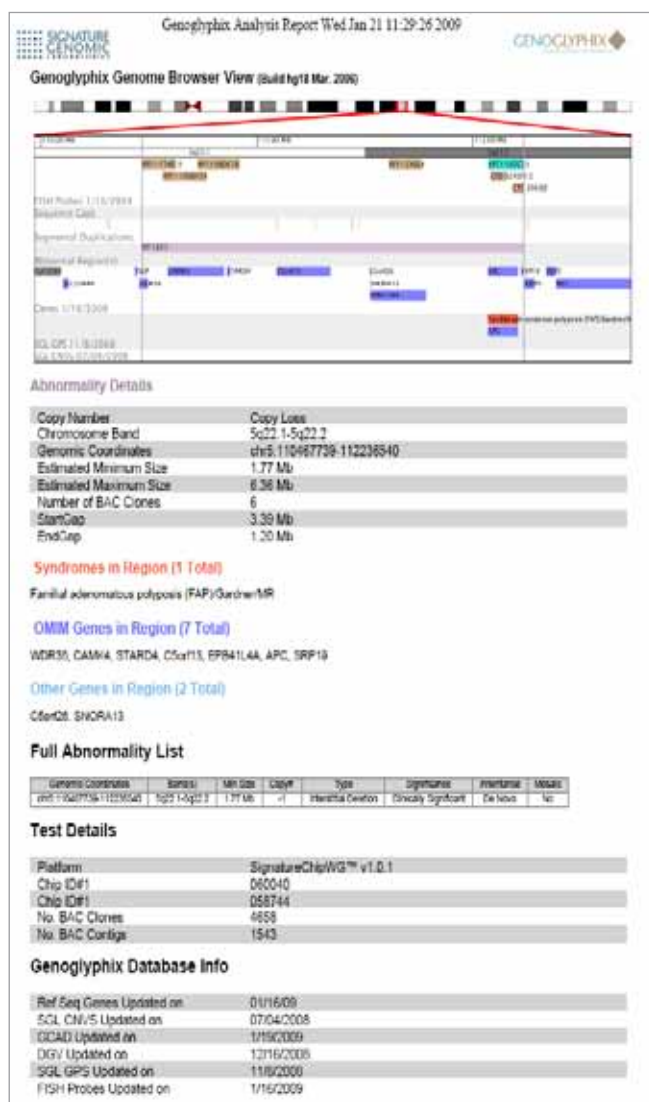
< Identify similar aberrations in Signature Genomics database of over 12,000 aberrations

Figure 3: Case Review - Linked Databases



Selection of the appropriate entries will automatically take you into databases such as OMIM, PubMed, and Genoglyphix Syndrome Detail Page

Figure 4: Final Report



Genoglyphix Workflow

1. Once CGX array results are uploaded, the data can be displayed either within whole-genome or chromosome specific graphs, or alternatively, by use of the Karyogram view (Figure 1).
2. Selecting aberrations of interest will automatically upload the Genoglyphix Genome browser (Figure 2). From here, different data bases such as OMIM, RefSeq, PubMed, DGV, as well as Signature's Genoglyphix Chromosome Aberration Database (GCAD) can be directly accessed (Figure 3). This database contains over 12,000 abnormal cases that can be viewed in context of the selected sample.
3. Data can be shared between users and can be made available as a community resource. The information will be displayed as a separate track within the Genoglyphix Genome Browser.
4. Interpretation of each abnormality within a DNA sample can be carried out within the Notation Review by classifying for example the aberration type, its clinical importance, and inheritance.
5. The Final Review and Report option offers the possibility to summarize all classified aberrations and report those within their genomic context. FISH information can be uploaded to be displayed within the same report (Figure 4).

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

PerkinElmer, Inc.
Wallac Oy
PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357

Signature Genomics/
PerkinElmer
2820 N. Astor Street
Spokane, WA 99207 USA
Phone: (877) 744-2447

ISO 13485
ISO 9001
CMDCAS
ISO 14001
OHSAS 18001



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

©2012 PerkinElmer, Inc. All rights reserved. PerkinElmer is a registered trademark of PerkinElmer, Inc. All other trademarks depicted are the property of their respective holders or owners. PerkinElmer reserves the right to change this document at any time and disclaims liability for editorial, pictorial or typographical errors.

1244-9985-01, March 2012