

CYTAG[®] TotalCGH

Labeling kit

Complete kit to digest and label DNA for
CGH+SNP arrays

Comparative genomic hybridization (aCGH) is a powerful diagnostic tool for detecting DNA copy number gains and losses associated with chromosome abnormalities. aCGH provides an understanding of genetic disorders, cancers and other genomic aberrations.

With optimized, proprietary reagents, Enzo's CYTAG[®] CGH Labeling Kits produce high quality data using as little as 0.25µg of genomic DNA, without a need for pre-amplification.

Bulk quantity kit, 2 x 1000 reactions, is available upon special request.

Ordering Information

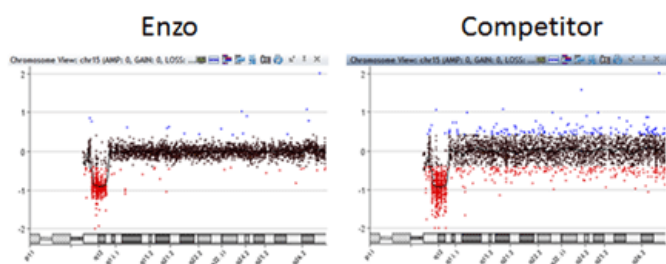
[Order Online »](#)

ENZ-42674-0100	2x100Reactions
ENZ-42674-0010	2x10Reactions

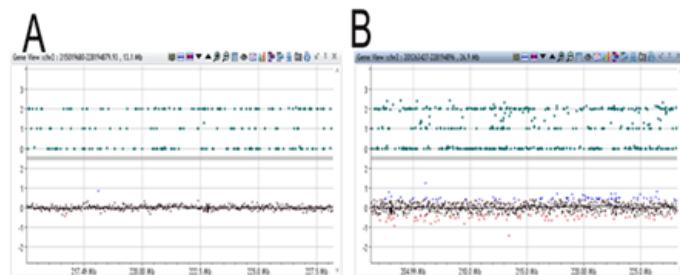
Manuals, SDS & CofA

[View Online »](#)

- Contains CYTAG[®] CGH Labeling kit, restriction enzymes *AluI* and *RsaI*, and PCR & Gel Clean-up Columns
- Compatible with CGH and CGH+SNP arrays
- Ability to perform total genomic DNA analysis without amplification or complexity reduction
- Provides excellent dynamic analytical range for challenging and complex, heterogeneous samples
- QC benchmarked and validated using high resolution arrays
- Superior DLR scores (0.09-0.12), exceeding industry standards



Comparative analysis of chromosome 15. Upon scanning, the quality of the labeling in DNA samples was visually inspected to demonstrate detection of known deletion in the Prader-Willi DNA. This demonstrates the superior labeling obtained with the CGH labeling kit from Enzo Life Sciences when compared with a competitor's kit.



SNP analysis of a segment of chromosome 2. DNA was labeled with the Enzo kit (left) or a competitor's kit (right). DNA was hybridized to Agilent's CGH+SNP microarray. The Enzo kit clearly identifies the sequence at SNPs while quantification of the SNP analysis of DNA labeled with a competitor's labeling system is not as clean.



Handling & Storage

Use/Stability	Store PCR & Gel Clean-up Columns at room temperature. All other components should be stored at -20°C. Stable for at least one year after receipt when stored as recommended.
Handling	Protect Cyanine 3- and 5-dUTPs from light at all times. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application CGH

Contents

-0010 size:

[ENZ-42671-K010](#): Primers/Reaction Buffer, 1 x 400 µl; Cyanine 3-dUTP Nucleotide Mix, 2 x 50 µl; Cyanine 5-dUTP Nucleotide Mix, 2 x 50 µl; Klenow Exo-DNA Polymerase, 1 x 20 µl; Stop Buffer, 1 x 100 µl; Nuclease-free Water, 1 x 1 ml

[ENZ-GEN100-0020](#): PCR & Gel Clean-up columns, 20 tests

[ENZ-GEN108-1000](#): AluI, 1000 units, 10X Cutting Buffer

[ENZ-GEN109-1000](#): RsaI, 1000 units, 10X Cutting Buffer

-0100 size:

[ENZ-42671-K100](#): Primers/Reaction Buffer, 1 x 4 ml; Cyanine 3-dUTP Nucleotide Mix, 1 x 1 ml; Cyanine 5-dUTP Nucleotide Mix, 1 x 1 ml; Klenow Exo-DNA Polymerase, 1 x 0.2 ml; Stop Buffer, 1 x 1 ml; Nuclease-free Water, 1 x 10 ml

[ENZ-GEN100-0200](#): PCR & Gel Clean-up columns, 200 tests

[ENZ-GEN108-1000](#): AluI, 1000 units, 10X Cutting Buffer

[ENZ-GEN109-1000](#): RsaI, 1000 units, 10X Cutting Buffer

Technical Info / Product Notes

Application Note:

[Using CYTAG[®] TotalCGH Labeling Kit for CGH + SNP Array](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: May 29, 2024