

# CYTAG<sup>®</sup> CGH

## Labeling kit

Superior labeling efficiency and better dye incorporation results in less failed runs

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, the Enzo Life Sciences CGH Labeling Kits for oligo arrays produce high quality data using as little as 0.25 µg of genomic DNA, without a need for pre-amplification. Bulk quantity kit, 2 x 1000 reactions, is available upon special request. For the purification of labeled DNA, use PCR & Gel Clean-up columns, Prod No. [ENZ-GEN100](#).

Citations: 65

[View Online »](#)

### Ordering Information

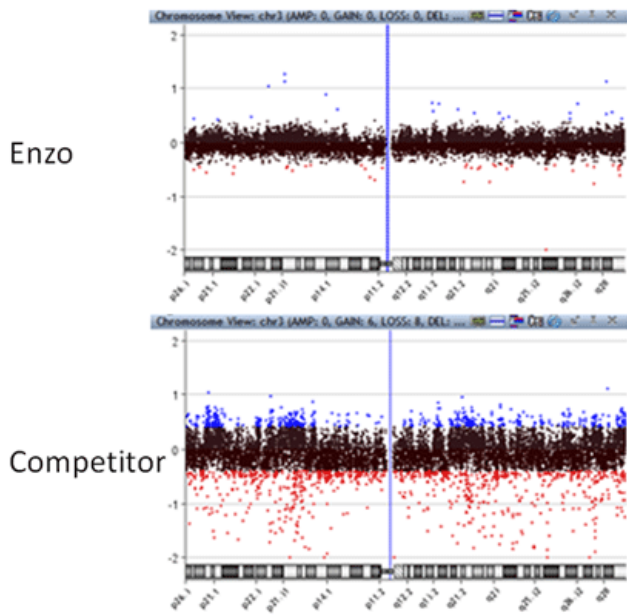
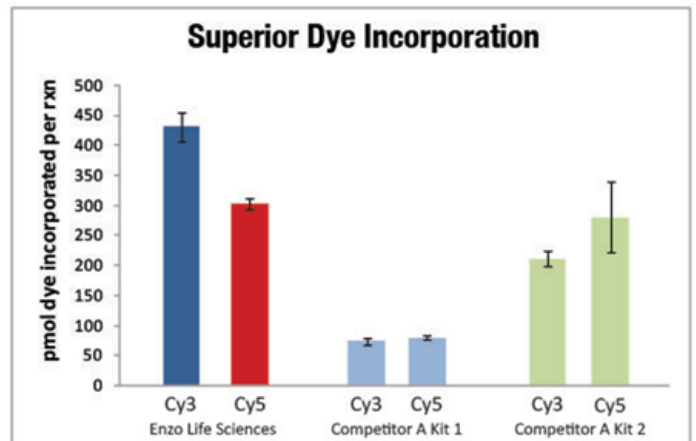
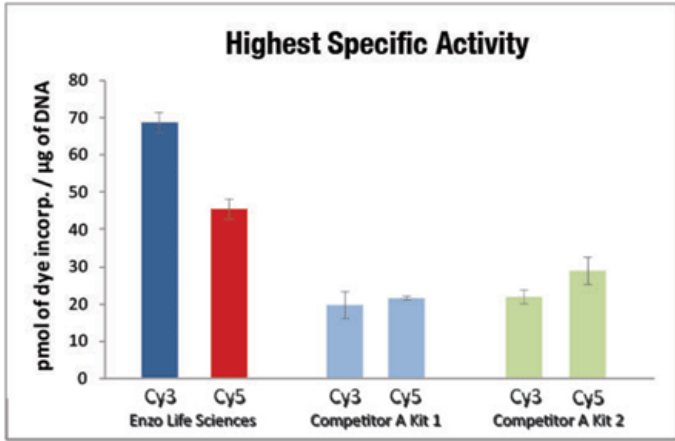
[Order Online »](#)

ENZ-42671-K010	2x10Reactions
ENZ-42671-K100	2x100Reactions

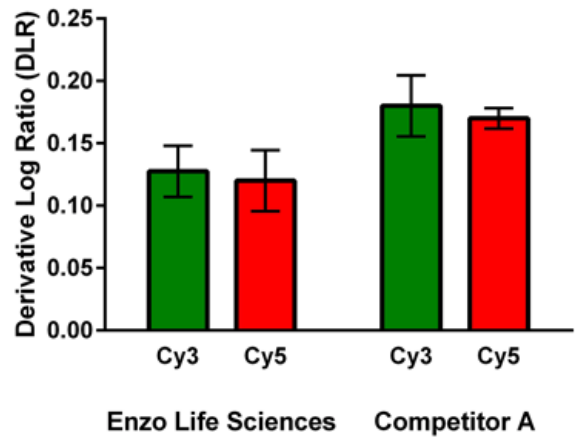
### Manuals, SDS & CofA

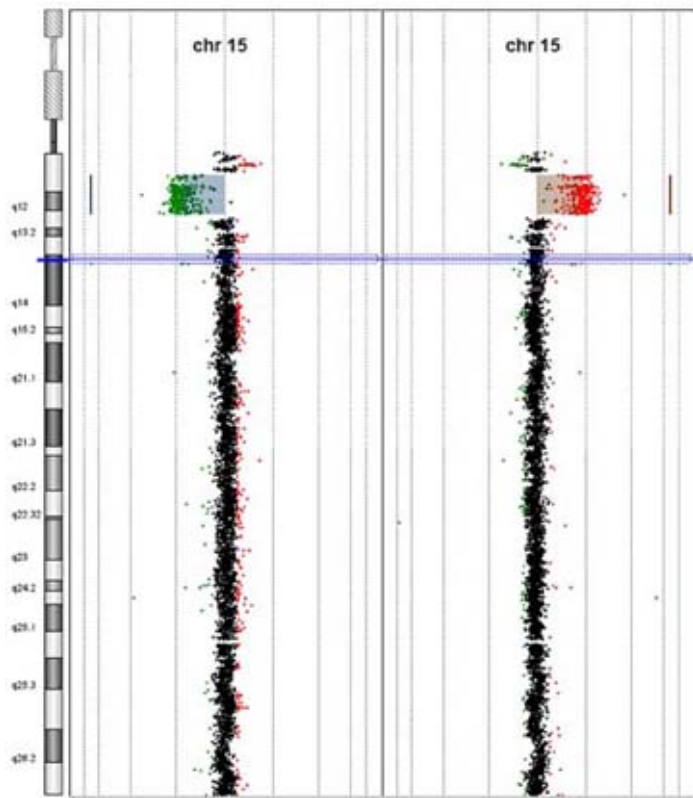
[View Online »](#)

- 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



### Lowest DLR scores result in less failed runs





Analysis of syndromic DNA using an oligonucleotide microarray (Agilent 4 x 180K) demonstrated the characteristic deletion in 15q11.2-q13 (chromosome 15) found in patients with Prader-Willi syndrome.



## Handling & Storage

Use/Stability	Stable until expiration date 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
Application Notes	Validated on Agilent SurePrint, OGT CytoSure™, and Illumina 24Sure Microarrays.
Contents	<p><b>-K010 size:</b> 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 DNA Polymerase, 1 x 20 µL; Stop Buffer, 1 x 100 µL; Nuclease-free Water, 1 x 1 mL</p> <p><b>-K100 size:</b> 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 DNA Polymerase, 1 x 0.2 mL; Stop Buffer, 1 x 1 mL; Nuclease-free Water, 1 x 10 mL</p>

## Technical Info / Product Notes

### Application Notes:

[Compatibility of CYTAG® CGH Labeling Kit from Enzo Life Sciences with SurePrint® G3 Human CGH 1x1M Microarrays from Agilent](#)

[Compatibility of CYTAG® CGH Labeling Kit with Illumina 24sure® Microarrays](#)

[CGH labeling protocol for small volumes of DNA sample](#)

[High Resolution Microarray Scanner used to Compare CGH Labeling Methods](#)

### Cited samples:

[For an overview of Cited Samples by Sample Type, please click here.](#)

[For an overview of Cited Samples by Array Type, please click here.](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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