

# CYTAG® SuperCGH Labeling Kit

Superior labeling efficiency for low input samples, with greater sensitivity and a simpler workflow. **Note: purification columns are not included; use the Columns recommended in the kit manual or your preferred purification protocol.**

With optimized, proprietary reagents, Enzo's CYTAG® SuperCGH Labeling Kit for oligo arrays produces high quality data using as little as 50 ng of genomic DNA, without a need for pre-amplification for precious low input samples. This kit does not include purification columns; use the Columns recommended in the kit manual or your preferred purification protocol.

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.

## Ordering Information

[Order Online »](#)

ENZ-GEN121-0010	10Reactions
ENZ-GEN121-0100	100Reactions

## Manuals, SDS & CofA

[View Online »](#)

- Suitable for previously limiting low input samples, including precious biopsy samples
- Validated for blood, amniotic fluid, postnatal, and oncology samples
- Provides excellent dynamic analytical range down to 50 ng
- Shorter workflow — samples ready for hybridization in <4 hours
- Superior DLR scores (0.09–0.12) exceeding industry standards
- Ability to perform total genomic DNA analysis without amplification or complexity reduction
- Purification columns not included — use the Columns recommended in the kit manual or your preferred purification protocol

# Handling & Storage

Use/Stability	Refer to label for expiration date. Store as recommended.
Handling	Protect Super Cyanine 3- and 5-dUTPs from light at all times. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Dry Ice

**Regulatory Status** RUO - Research Use Only

# Product Details

**Application** CGH

**Contents**

**-0010 size (10 reactions):**  
Primers/Reaction Buffer, 1 x 400 µL; Super Cyanine 3-dUTP® Nucleotide Mix, 1 x 100 µL; Super Cyanine 5-dUTP® Nucleotide Mix, 1 x 100 µL; Klenow DNA Polymerase, 1 x 40 µL; Stop Buffer, 1 x 100 µL; Nuclease-free Water, 1 x 1 mL

**-0100 size (100 reactions):**  
Primers/Reaction Buffer, 1 x 4 mL; Super Cyanine 3-dUTP® Nucleotide Mix, 1 x 1 mL; Super Cyanine 5-dUTP® Nucleotide Mix, 1 x 1 mL; Klenow DNA Polymerase, 1 x 0.4 mL; Stop Buffer, 1 x 1 mL; Nuclease-free Water, 1 x 10 mL

**Quality Control** Expiration as stated on the label when stored as recommended.

**Specificity** Validated on Agilent SurePrint Microarrays

**Technical Info / Product Notes**

**Application Notes:**  
[Using CYTAG® SuperCGH Labeling Kit for the Analysis of FFPE Oncological Samples by CGH.](#)  
[Automated Liquid Handling for High Throughput DNA Labeling with CYTAG® SuperCGH Labeling Kit.](#)  
[Sensitivity of Comparative Genomic Hybridization Using Artificial Mosaicism and Small DNA Quantities.](#)



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)