

The logo for CYTAG, featuring the word "CYTAG" in a bold, dark blue sans-serif font, followed by a registered trademark symbol (®).

CYTAG®

SuperCGH+SNP Labeling Kit

Optimized to label DNA for dual comparative genomic hybridization (CGH+SNP or CGH

only) for low DNA input samples. 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.

The CYTAG® SuperCGH+SNP Labeling Kit has been designed to provide higher labeled DNA yield, while still maintaining high specific activity for low-input DNA samples, as low as 100 ng for CGH+SNP and 50 ng for CGH only.

Ordering Information

[Order Online »](#)

ENZ-GEN122-0010	2x10Reactions
ENZ-GEN122-0100	2x100Reactions

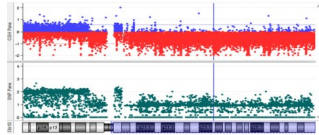
Manuals, SDS & CofA

[View Online »](#)

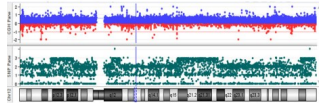
- Suitable for previously limiting low input samples, including precious biopsy samples
- Validated for blood, amniotic fluid, postnatal, and oncology samples
- Provides excellent dynamic analytical range down to 50 ng
- Shorter workflow has samples ready for hybridization in < 4 hours
- Superior DLR scores (<0.18)
- Ability to perform total genomic DNA analysis without amplification or complexity reduction



CYTAG® SuperCGH+SNP Labeling Kit

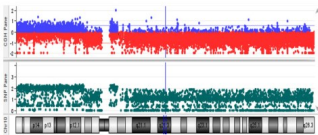


A. Frozen_1, Chromosome 10.

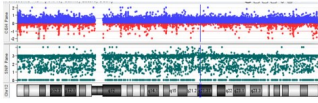


C. Frozen_3, Chromosome 12.

Competitor



B. Frozen_1 Chromosome 10.



D. Frozen_3, Chromosome 12.

Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect Super 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 Microarrays
Contents	Labeling Kit and Restriction Enzymes

Last modified: September 26, 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)