CYTAG® SuperCGH+SNP Labeling Kit

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

Gorngarative genominate bridgettos (and all 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 »

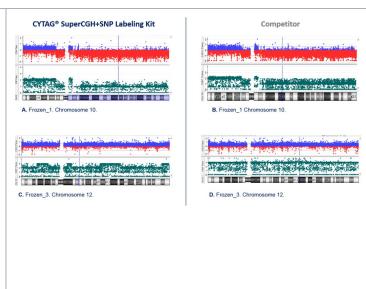
ENZ-GEN122-0010	2x10Reactions
ENZ-GEN122-0100	2x100Reactions

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





Handling & Storage

Use/Stability Stable for at least one year after reciept 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



info-

eu@enzolifesciences.com