

# CGH labeling kit for BAC arrays

For preparation of Cyanine 3- and Cyanine 5-labeled DNA for hybridization to BAC arrays. The CGH Labeling Kit for BAC Arrays has been optimized to label DNA for dual sample comparative genomic hybridization (CGH) assays on BAC arrays. All required reagents are provided including fully optimized nucleotide mixes that contain either cyanine 3-dUTP or cyanine 5-dUTP. The CGH Labeling Kit for BAC Arrays has been specifically tailored to meet the varying requirements of individual laboratories by providing sufficient amount of reagents for two different protocols. For the average user, the Standard Protocol calls for a 4-hour labeling in a 50µl reaction volume. For users with dilute DNA ( $\leq 20$  ng/µl), the Dilute DNA Protocol calls for a 4-16 hour labeling in a total volume of 100µl.

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-42670	20Reactions
-----------	-------------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	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	Primers/Reaction Buffer, 2 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
Quantity	Contains reagents sufficient for 20 reactions or 10 reactions in duplicate with a dye swap.

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)