

Cyanine 3-NHS ester pack

Enzo cyanine 3 NHS-ester is a reactive, water-soluble fluorescent dye that provides bright orange signal. The dye can be chemically linked to either nucleic acids or proteins and is suitable for a wide variety of applications including microarray analysis and protein tagging. In microarray experiments, DNA or RNA samples are labeled with the cyanine 3 and cyanine 5 dyes for differential display analysis. Since NHS esters react only with aliphatic amine groups, which nucleic acids lack, nucleotides must be modified first with aminoallyl groups thru incorporating aminoallyl-modified nucleotides during synthesis. Alternatively, labeling can be performed directly by nick translation using cyanine dUTPs (Gold 550 dUTP, Prod. No. ENZ-42521, and Red 580 dUTP, Prod. No. ENZ-42844). For protein labeling, it is critical to control the number of dye molecules affixed per protein to maintain maximal activity. For example, no more than 8 dye molecules should be bound per antibody molecule, while annexin V will only tolerate 1-2 dye molecules per protein without loss of activity. Wavelength Maxima: Excitation 553nm, Emission 570nm

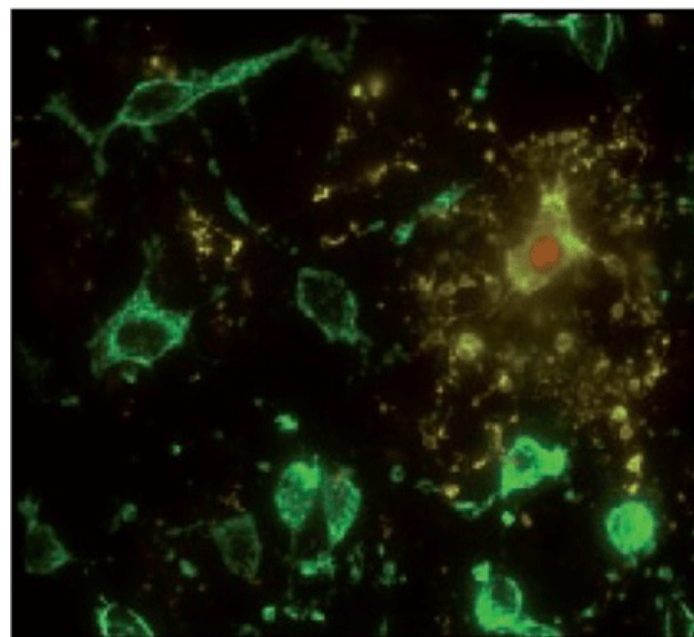
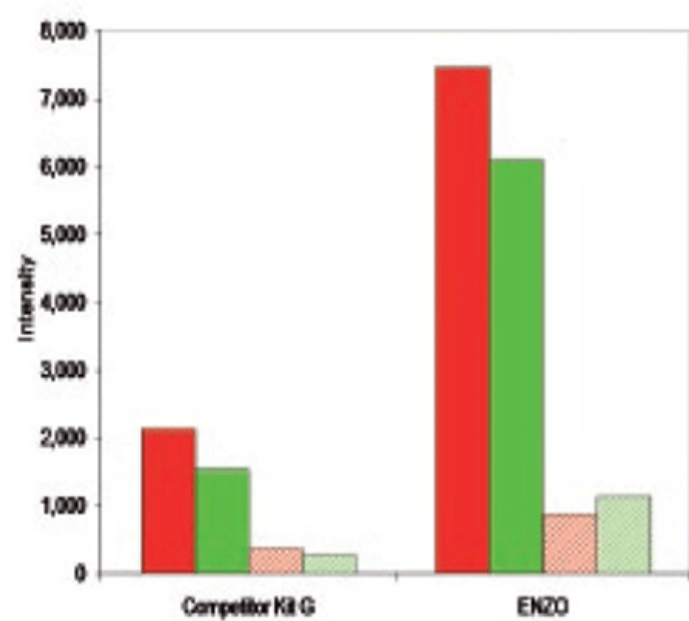
Ordering Information

[Order Online »](#)

ENZ-42541	12x50nmol
-----------	-----------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture. Protect from light.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Technical Info / Product Notes

This product is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.

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)