

Triphosphate analogue with artificial base for directed mutagenesis experiments.

BLG-D114-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

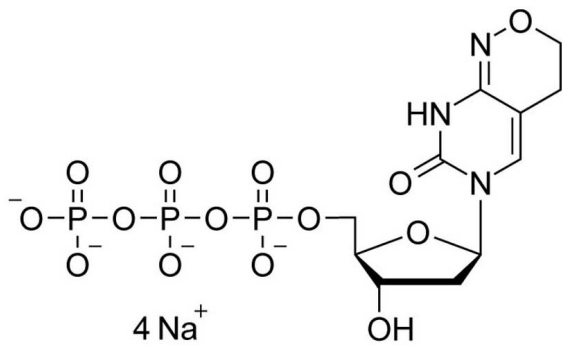
Ordering Information

[Order Online »](#)

BLG-D114-05	5 μ mol
BLG-D114-25	5x5 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-(β-D-2-Deoxyribofuranosyl)-3, 4-dihydro-8H-pyrimido-[4, 5-c]-[1, 2]-oxazin-7-one-5'-O-triphosphate, 2'-Deoxy-P-nucleoside-5'-O-triphosphate . tetrasodium salt

Appearance 10 mM aqueous solution

CAS 173964-83-5

Couple Type Modified nucleotides

Formula $C_{11}H_{18}N_3O_{14}P_3$ (free acid)

MW 509.2

Purity > 95% HPLC

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com