

3'-O-Ac-dGTP

Ribose-modified dGTP analogue

3'-O-Ac-dGTP is an analogue of dGTP with blocked 3'-hydroxy function, which can be used as a transient, pH-sensitive chain terminator during DNA polymerization. The acetyl group of 3'-O-Ac-dGTP is sensitive towards alkaline conditions and can be easily split off to allow further subsequent polymerization steps.

BLG-A278-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

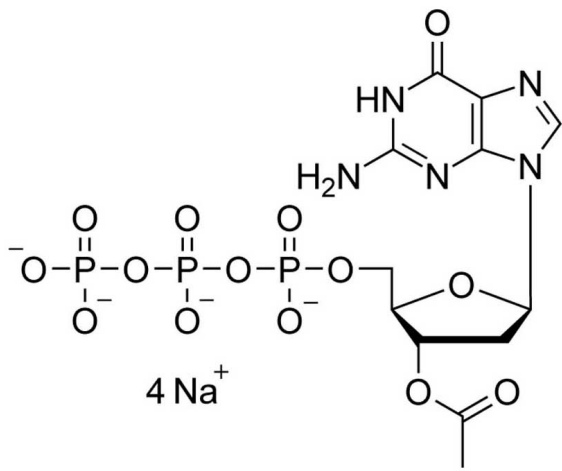
Ordering Information

[Order Online »](#)

BLG-A278-05	5 μ mol
BLG-A278-25	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 3'-O-Acetyl-2'-deoxyguanosine-5'-O-triphosphate . sodium salt

CAS 90290-77-0

Formula $C_{12}H_{18}N_5O_{14}P_3 \cdot 4 Na$

Formulation 10 mM aqueous solution

MW 549.2 (free acid)

Purity > 95% HPLC



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-be@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