

3'-O-Piv-dATP

Ribose-modified dATP analogue

3'-O-Piv-dATP is an analogue of dATP with blocked 3'-hydroxy function, potentially useful as a transient, pH-sensitive chain terminator during DNA polymerization. The pivaloyl group of 3'-O-Piv-dATP is sensitive towards alkaline conditions and can be split off to allow further subsequent polymerization steps.

BLG-D165T-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

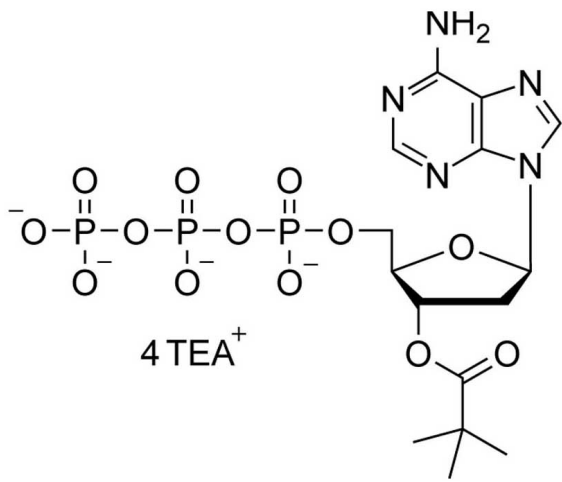
Ordering Information

[Order Online »](#)

BLG-D165T-01	1µmol
BLG-D165T-05	5x1µmol

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 2'-Deoxy-3'-O-pivaloyladenine-5'-O-triphosphate

Couple Type Modified nucleotides

Formula $C_{15}H_{24}N_5O_{13}P_3$

MW 575.3

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.6 mg

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-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