

3'-O-Ac-dCTP

Ribose-modified dCTP analogue

3'-O-Ac-dCTP is an analogue of dCTP 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-dCTP is sensitive towards alkaline conditions and can be easily split off at pH > 8 to allow further subsequent polymerization steps.

BLG-A277-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

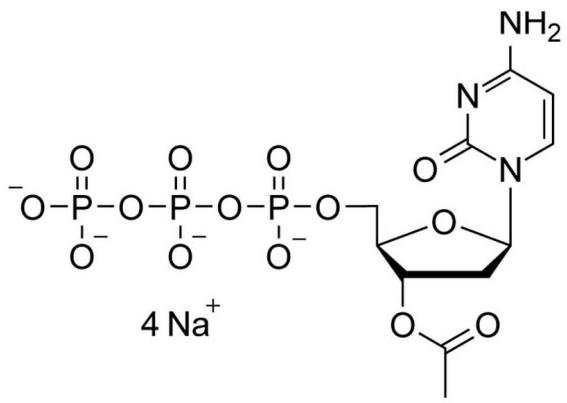
Ordering Information

[Order Online »](#)

BLG-A277-05	5µmol
BLG-A277-25	5x5µmol

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'-deoxycytidine-5'-O-triphosphate . sodium salt

CAS 90290-80-5

Formula $C_{11}H_{18}N_3O_{14}P_3 \cdot 4 Na$

Formulation 10 mM aqueous solution

MW 509.2 (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