

5-Methyltetrazine-PO- (EO)₄-P-AA-dCTP / 5- Methyltetrazine-PEG₄- AA-dCTP

5-Methyltetrazine-PEG₄-AA-dCTP is an analogue of dCTP and a suitable molecular tool for the so-called "Tetrazine-Alkene-Ligation", a copper(I)-free inverse electron demand Diels-Alder (IEDDA) reaction. The IEDDA reaction is a prime example of the click chemistry technique and has attracted particular interest for the introduction of reporter groups into oligonucleotides and for the labelling of various biomolecules.

BLG-M171-0125 (5 x 0.25 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

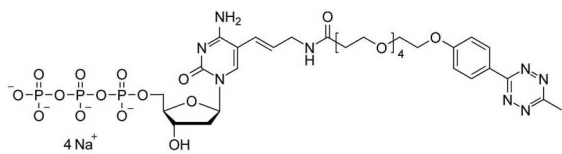
Ordering Information

[Order Online »](#)

BLG-M171-0025	0.25µmol
BLG-M171-0125	5x0.25µmol

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 5-(3-(15-((6-Methyl-1,2,4,5-tetrazine-3-yl)phenoxy)-(tetraethoxy)-propionyl)aminoallyl)-2'-deoxycytidine-5'-O-triphosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{32}H_{47}N_8O_{19}P_3$

Formulation 1 mM aqueous solution

MW 940.7

Purity > 95% HPLC

Quantity 0.25µmol are equivalent to ~235 µg



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