

# 4-Methyltetrazine-PO- (EO)<sub>4</sub>-P-AH-dCTP / 4- Methyltetrazine-PEG<sub>4</sub>- AH-dCTP

4-Methyltetrazine-PEG<sub>4</sub>-AH-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-M166-0125 (5 x 0.25 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

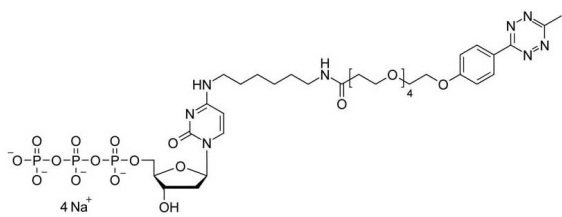
## Ordering Information

[Order Online »](#)

BLG-M166-0025	0.25µmol
BLG-M166-0125	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** N4-(6-(15-((6-Methyl-1,2,4,5-tetrazine-3-yl)phenoxy)-(tetraethoxy)-propionyl)aminoethyl)-2'-deoxyucytidine-5'-O-triphosphate . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{35}H_{55}N_8O_{19}P_3$

**Formulation** 1 mM aqueous solution

**MW** 984.8

**Purity** > 95% HPLC

**Quantity** 0.25µmol are equivalent to ~246 µg



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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