

5-N₃Ac-AA-dCTP

Azide-containing analogue of dCTP

The azide-containing 5-N₃Ac-AA-dCTP is an analogue of dCTP and a suitable molecular tool for the copper(I)-catalysed Huisgen azide-alkyne [3+2] cycloaddition (CuAAC) or copper(I)-free strain-promoted Alkyne-Azide Click Chemistry (SPAAC) with strained alkynes, respectively. Both reactions have attracted particular interest for the introduction of reporter groups and for the labelling of various biomolecules like dCTP-binding proteins by the so-called click-chemistry technique. 5-Azide-functionalised pyrimidine nucleotides can replace their natural counterpart dCTP in polymerase reactions, resulting in functionalized DNA that can subsequently be further reacted and modified using click chemistry methods.

BLG-A385-0125 (5 x 0.25 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

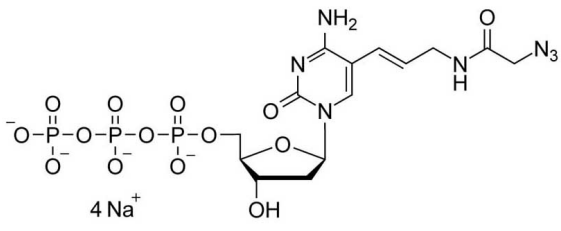
[Order Online »](#)

BLG-A385-0125

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-(2-Azidoacetyl)aminoallyl)-2'-deoxycytidine-5'-O-triphosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{14}H_{22}N_7O_{14}P_3$ (free acid)

MW 605.3 (free acid)

Purity > 95% HPLC

Quantity 0.25µmol = ~151µg



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