

7-N₃Ac-AParg-7-C-dATP

Azide-containing analogue of dATP

The azide-containing 7-N₃Ac-AParg-7-C-dATP is an analogue of dATP 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 dATP-binding proteins by the so-called click-chemistry technique. Azide-functionalised 7-deaza purine nucleotides can replace their natural counterpart dATP in polymerase reactions, resulting in functionalized DNA that can subsequently be further reacted and modified using click chemistry methods.

BLG-A383-0125 (5 x 0.25 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

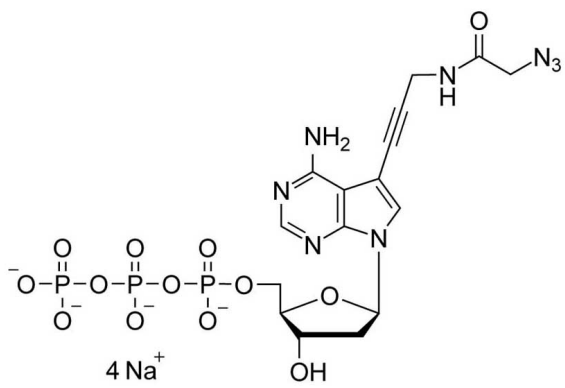
Ordering Information

[Order Online »](#)

BLG-A383-0125	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 7-(3-(2-Azidoacetyl)aminopropargyl)-7-deaza-2'-deoxyadenosine-5'-O-triphosphate . sodium salt

CAS 2570307-48-9

Couple Type Modified nucleotides

Formula $C_{16}H_{21}N_8O_{13}P_3$ (free acid)

MW 626.3 (free acid)

Purity > 95% HPLC

Quantity 0.25µmol = ~157µg



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