

8-Bu(3-yne)T-ATP

Alkyne-containing analogue of ATP

BLG-B302-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

Ordering Information

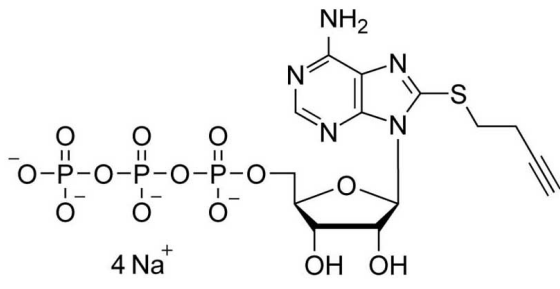
[Order Online »](#)

BLG-B302-01	1 μ mol
BLG-B302-05	5x1 μ mol

Manuals, SDS & CofA

[View Online »](#)

The alkyne-containing 8-Bu(3-yne)T-ATP is an analogue of ATP and a suitable molecular tool for the copper-catalysed Huisgen azide-alkyne [3+2] cycloaddition (CuAAC), which has attracted particular interest for the introduction of reporter groups and for the labelling of various biomolecules like ATP-binding proteins by the so-called click-chemistry technique.



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-(3-Butynylthio)adenosine-5'-O-triphosphate . sodium salt

Formula $C_{14}H_{20}N_5O_{13}P_3S$ (free acid)

MW 591.3 (free acid)

Purity > 95% HPLC

Quantity 1 μ mol = ~0.6 mg

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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