

# Biotin-14-dATP, S pack

## Ordering Information

[Order Online »](#)

JBS-NU-835-BIO14-S	200µl
--------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for up to 1 year after delivery if stored correctly.
<b>Short Term Storage</b>	Ambient
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Biotin-14-N6-(6-Amino)hexyl-dATP, Triethylammonium salt, Bio-14-dATP
<b>Appearance</b>	colorless to slightly yellow solution
<b>Application Notes</b>	Incorporation into DNA/cDNA by – Nick Translation with DNase I/ DNA Polymerase I Primer Extension with Klenow fragment
<b>Couple Type</b>	Modified nucleotides
<b>Formula</b>	$C_{32}H_{54}N_9O_{15}P_3S$ (free acid)
<b>Formulation</b>	sterile solution in 10 mM Tris-HCl
<b>MW</b>	929.81 (free acid)
<b>Purity</b>	≥ 95 % (HPLC)
<b>Technical Info / Product Notes</b>	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)