

# Rhodamine-12-dUTP, S pack

## Ordering Information

[Order Online »](#)

JBS-NU-803-RHOX-S	10µl
-------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Rhodamine-X-5-Aminoallyl-dUTP, 5/6-Rhodamine-X-5-Aminoallyl-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

**Appearance** orange solution

**Application Notes** Incorporation into DNA/cDNA by – Nick Translation with DNase I/ DNA Polymerase I

**Couple Type** Modified nucleotides

**Emission Maximum** 530 nm

**Excitation Maximum** 505 nm

**Formula**  $C_{39}H_{43}N_6O_{19}P_3$  (free acid)

**Formulation** solution in water

**MW** 992.71 (free acid)

**Purity** ≥ 95 % (HPLC)

**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](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)