

Rp-UTP- α -S

Useful for inhibition or activation of UTP-responsive receptors and for determination of their stereospecificity. Increased metabolic stability compared to UTP. Not accepted by RNA polymerases.

BLG-U002-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

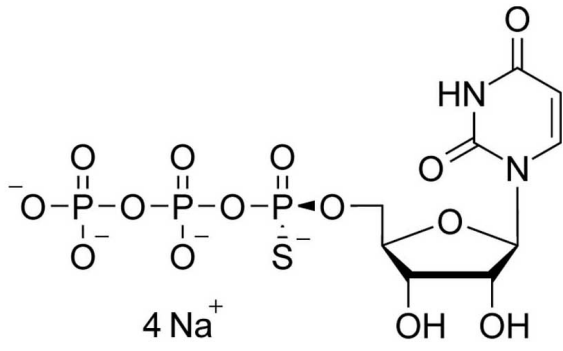
Ordering Information

[Order Online »](#)

BLG-U002-05	5 μ mol
BLG-U002-25	5x5 μ mol

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 Uridine-5'-O-(1-thiotriphosphate) . tetrasodium salt, Rp-isomer

Appearance 10 mM aqueous solution

CAS 71214-30-7

Couple Type Modified nucleotides

Formula $C_9H_{15}N_2O_{14}P_3S$ (free acid)

MW 500.2 (free acid)

Purity > 95% HPLC

Quantity 5µmol≈2.5mg

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