

2'-Fluoro-dNTP Bundle

Ordering Information

[Order Online »](#)

JBS-NU-10405-20	4x50µl
-----------------	--------

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 2'-Fluoro-2'-deoxynucleoside-5'-triphosphates, Sodium salts, 2'-Deoxy-2'-fluoronucleoside-5'-triphosphates, 2'-F-2'-dNTP, 2'F-dNTP Bundle

Appearance colorless to slightly yellow solution

Couple Type Modified nucleotides

Formula
2'Fluoro-dATP: $C_{10}H_{15}FN_5O_{12}P_3$ (free acid)
2'Fluoro-dCTP: $C_9H_{15}FN_3O_{13}P_3$ (free acid)
2'Fluoro-dGTP: $C_{10}H_{15}FN_5O_{13}P_3$ (free acid)
2'Fluoro-dUTP: $C_9H_{14}FN_2O_{14}P_3$ (free acid)

Formulation solution in water

MW 2'Fluoro-dATP: 509.17 (free acid) 2'Fluoro-dCTP: 485.15 (free acid) 2'Fluoro-dGTP: 525.17 (free acid) 2'Fluoro-dUTP: 486.13 (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

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-fr@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