

Biotin-16-dCTP, S pack

May require a license for import, please [contact us](#) for more information.

Ordering Information

[Order Online »](#)

JBS-NU-944-5	5mg
--------------	-----

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 m7GpF, 7-methyl-guanosine-5'-(α -fluoro)-monophosphate, 7-methyl-guanosine-5'-(1-fluoro)-monophosphate

Appearance white to off-white solid

CAS 1698011-01-6 (inner salt)

Couple Type Modified nucleotides

Formula $C_{11}H_{15}FN_5O_7P$

MW 379.24

Purity $\geq 95\%$ (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-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