

UDP-6-azide-glucose

Ordering Information

[Order Online »](#)

JBS-CLK-076	0.5mg
-------------	-------

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 UDP-6-N3-Glc, 6-N3-UDPG

Appearance white to off-white solid

Application Notes Glycosylation of 5-hydroxymethylcytosine (5-hmC) with T4 β-Glycosyltransferase (T4 β-GT). The resulting azide-functionalized, glycosylated 5-hmC moiety can subsequently be detected via Cu(I)-free click chemistry that offers the choice to introduce a (Desthio)Biotin group for subsequent purification tasks (via DBCO-containing (Desthio)Biotin) to introduce fluorescent group for subsequent microscopic imaging (via DBCO-containing fluorescent dyes). 0.5 mg UDP-6-azide-glucose are sufficient to prepare 283 μl of a 3 mM solution.

Formula $C_{15}H_{23}N_5O_{16}P_2$

MW 591.32

Purity ≥ 98 % (HPLC)

Solubility water, aqueous buffer such as PBS

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