

UDP-GalNAc

Nucleotide sugar and donor substrate for N-acetylgalactosaminyltransferases (GalNAcTs)

BLG-U027-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

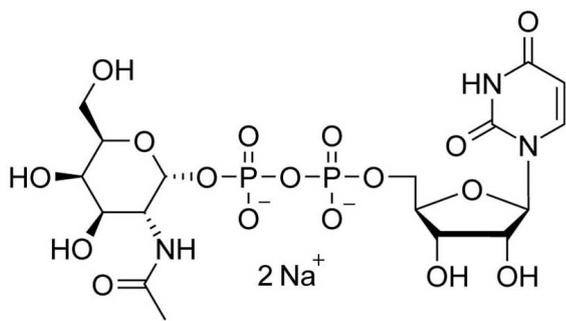
[Order Online »](#)

BLG-U027-10	10 μ mol
BLG-U027-50	5x10 μ mol

Manuals, SDS & CofA

[View Online »](#)

UDP-N-acetyl- α -d-galactosamine (UDP-GalNAc) disodium salt is a nucleotide sugar and the donor substrate for eukaryotic and prokaryotic N-acetylgalactosaminyltransferases (GalNAcTs) involved in mucin, glycosphingolipid, and glycosaminoglycan synthesis.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Uridine-5'-diphospho-N-acetylgalactosamine . sodium salt

CAS 108320-87-2

Formula $C_{17}H_{27}N_3O_{17}P_2$ (free acid)

MW 607.4 (free acid)

Purity > 98% HPLC

Quantity 10 µmol are equivalent to ~ 6.1 mg



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