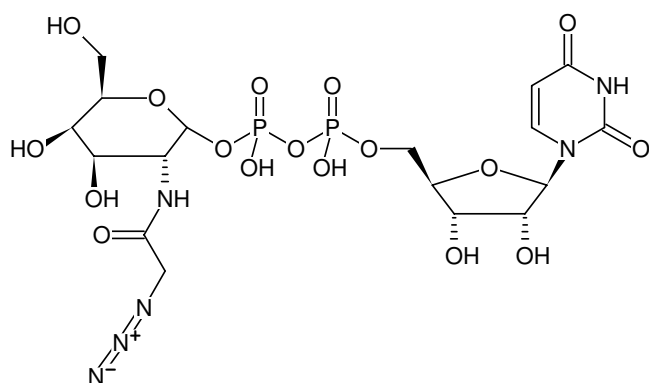


**UDP-GalNAz**

UDP-N-azidoacetylgalactosamine
 UDP-N-azidoacetylgalactosamine, Sodium salt

Cat. No.	Amount
CLK-077	0,5 mg



Structural formula of UDP-GalNAz

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₇H₂₆N₆O₁₇P₂ (free acid)

Molecular Weight: 648.37 g/mol (free acid)

Exact Mass: 648.08 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: colorless to white

Solubility: water, aqueous buffer pH 7-7.5

Spectroscopic Properties: λ_{max} 262 nm, ε 10.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Transfer of GalNAz moiety to GalNAc residues of proteins and monoclonal antibodies with β1,4-galactosyltransferase mutant Y289L-Gal-T1 [1-2].

The azide-functionalized target can subsequently be detected via Cu(I)-catalyzed or Cu(I)-free Click Chemistry that offers the choice to introduce

- a Biotin group (via Azides of Biotin or DBCO-containing Biotin, respectively) for subsequent purification tasks or
- a fluorophore (via Azides of fluorescent dyes or DBCO-containing fluorescent dyes, respectively) for subsequent microscopic imaging

Related Products:

UDP-6-azide-glucose, #CLK-076

Ac4GalNAz, #CLK-1086

Selected References:

[1] Hang *et al.* (2004) Probing glycosyltransferase activities with the Staudinger ligation. *J. Am. Chem. Soc.* **126**(1):6.

[2] Qasba *et al.* (2008) Site-specific linking of biomolecules via glycan residues using glycosyltransferases. *Biotechnol. Prog.* **24**(3):520.