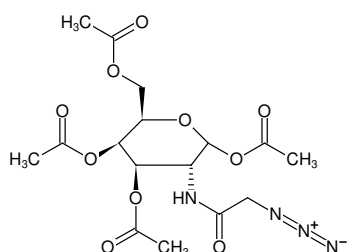


**Ac₄GalNAz**N-azidoacetylgalactosamine-tetraacylated (Ac₄GalNAz)

Cat. No.	Amount
CLK-1086-5	5 mg
CLK-1086-25	25 mg
CLK-1086-100	100 mg

Structural formula of Ac₄GalNAz**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₁₆H₂₂N₄O₁₀**Molecular Weight:** 430.37 g/mol**Exact Mass:** 430.13 g/mol**CAS#:** 653600-56-7**Purity:** mass identification (ESI-MS)**Form:** amorphous solid**Color:** off-white to grey**Solubility:** DMSO, DMF, MeOH**Applications:**

Glycoconjugate synthesis monitoring by metabolic labeling

Description:

The tetraacylated N-Azidoacetyl-galactosamine (Ac₄GalNAz) provides a non-radioactive alternative for glycoconjugate visualization. It is cell-permeable, intracellularly processed and incorporated instead of its natural monosaccharide counterpart N-Acetylgalactosamine (GalNAc).

The resulting Azide-functionalized glycoconjugates 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 to introduce fluorescent group (via Azides of fluorescent dyes or DBCO-containing fluorescent dyes, respectively) for subsequent microscopic imaging.

Recommended concentration for metabolic labeling: 25-75 μM. This concentration range may serve as a starting point for an individual assay set-up.