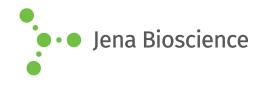
DATA SHEET





6-Azido-Trehalose

6-TreAz

| Cat. No. | Amount |
|----------|--------|
| CLK-078 | 1 mg |

Structural formula of 6-Azido-Trehalose

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: C₁₂H₂₁N₃O₁₀ Molecular Weight: 367.31 g/mol

Exact Mass: 367.12 g/mol

Purity: ≥ 90 % (TLC)

Form: solid

Color: white to off-white **Solubility:** water, DMSO

Applications:

Metabolic Lipopolysaccharide labeling of Myobacteria

Description:

Myobacteria are characterized by a trehalose-lipopolysaccharidecontaining outer cell membrane.

The Azide-functionalized 6-Azido-Trehalose provides an attractive approach for the detection of *Myobacteria* in living cultures: It is cell-permeable, intracellularly processed and incorporated instead of its natural counterpart.

The resulting Azide-functionalized lipopolysaccharides can subsequently be detected via Cu(I)-catalyzed or Cu(I)-free Click Chemistry that offers the choice to introduce a Biotin group (via Alkynes of Biotin or DBCO-containing Biotin, respectively) for subsequent purification tasks or to introduce fluorescent group (via Alkynes of fluorescent dyes or DBCO-containing fluorescent dyes, respectively) for subsequent microscopic imaging.

Myobacteria strains tested^[1]:

Mycobacterium smegmatis mc2155, Mycobacterium tuberculosis

H37Rv,Mycobacterium bovis BCG

Selected References:

[1] Swarts *et al.* (2012) Probing the Mycobacterial Trehalome with Bioorthogonal Chemistry. *J. Am. Chem. Soc.* **134 (39)**:16123.