# **DATA SHEET**





# Na-Ascorbate - click chemistry grade

L-Ascorbic acid sodium salt

Cat. No.	Amount
CLK-MI005-1G	5 x 200 mg

For general laboratory use.

Shipping: shipped at ambient temperature Storage Conditions: store at ambient temperature Additional Storage Conditions: store dry Shelf Life: 12 months after date of delivery Molecular Formula: C<sub>6</sub>H<sub>7</sub>NaO<sub>6</sub> Molecular Weight: 198.11 g/mol CAS#: 134-03-2 Form: solid Color: white to off-white Solubility: water

## Description:

Na-Ascorbate can be used as a reduction reagent for Cu(I)-catalyzed Alkyne-Azide click chemistry reactions (CuAAC).

It catalyzes the reduction of Cu(II) sources such as  $CuSO_4$  thereby releasing catalytically reactive Cu(I) ions.

Ideally, solutions should be freshly prepared in  $ddH_2O$  shortly before use. Alternatively, a stock solution can be prepared, stored at -20°C and freshly by diluted shortly before use.

Please note: Do not use solutions that appear brown. Freshly prepared, fully functional Na-Ascorbate solutions are colorless and turn brown upon oxidization thereby loosing their reduction capability.

Presolski *et al.*<sup>[1]</sup> and Hong *et al.*<sup>[2]</sup> provide a general protocol for CuAAC reactions that may be used as a starting point for the set up and optimization of individual assays.

#### **Related Products:**

Copper (II)-Sulphate (CuSO<sub>4</sub>), #CLK-MI004 THPTA, #CLK-1010 BTTAA, #CLK-067

### Selected References:

 Presolski et al. (2011) Copper-Catalyzed Azide-Alkyne Click Chemistry for Bioconjugation. *Current Protocols in Chemical Biology* 3:153.
Hong et al. (2011) Analysis and Optimization of Copper-Catalyzed Azide-Alkyne Cycloaddition for Bioconjugation. *Angew. Chem. Int. Ed.* 48:9879.

