# **DATA SHEET**





# ■ CuSO₄ - click chemistry grade

(Copper(II)-Sulphate)

Cat. No.	Amount
CLK-MI004-50	5 x 10 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: CuSO<sub>4</sub>
Molecular Weight: 159.61 g/mol

CAS#: 7758-98-7

**EC number:** 231-847-6

Form: solid
Color: blue
Solubility: water

#### **Description:**

CuSO<sub>4</sub> can be used as a copper source for Cu(I)-catalyzed Alkyne-Azide click chemistry reactions (CuAAC).

Catalytically reactive Cu(I) ions are released by reduction reagents such as Na-Ascorbate.

A stock solution can be prepared in  $ddH_2O$  and subsequently be stored at  $4^{\circ}C.$ 

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:**

Sodium Ascorbate (Na-Ascorbate), #CLK-MI005 THPTA, #CLK-1010 BTTAA, #CLK-067

## **Selected References:**

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