

**CuSO₄ - click chemistry grade**

(Copper(II)-Sulphate)

Cat. No.	Amount
CLK-MI004-50	5 x 10 mg

For research use only!**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:** CuSO₄**Molecular Weight:** 159.61 g/mol**CAS#:** 7758-98-7**EC number:** 231-847-6**Purity:** > 98 %, anhydrous**Form:** blue solid**Solubility:** water**Description:**

CuSO₄ 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 ddH₂O and subsequently be stored at 4°C.

Presolski *et al.*^[1] and Hong *et al.*^[2] 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

BTAA, #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.