

Chloroquine Chloroquine diphosphate salt

| Cat.-No. | Amount |
|----------|--------|
| CPP-A03 | 1g |

For *in vitro* use only
Quality guaranteed for 12 months
Store at 4°C

EU Harmful

Harmful if swallowed. Do not breathe dust. Avoid contact with skin and eyes. Target organ(s): Eyes. Heart. Light sensitive.

Product Information

CAS Number: 50-63-5

Molecular Formula

$C_{18}H_{26}ClN_3 \cdot 2H_3PO_4$

Molecular Weight

515.9 Da

Melting point

193 °C

pKa

8.4

Solubility

soluble in water (50 mg/ml)

Description

During internalization some mechanisms lead to formation of intracellular vesicles. For release of cargo from these vesicles they must be destabilized. For this purpose chloroquine have been proven effective when added to the medium. Add chloroquine to the serum containing medium to achieve a concentration of 150 µM.

For more information see **general manual (4.2.2)**.

Selected References

1. *Clarke's Isolation and Identification of Drugs*, Ed. by Moffat, A. C. *et al.*, The Pharmaceutical Press, London (1986)
2. *Krajewski, W.A.* (1999) Effect of In Vivo Histone Hyperacetylation on the State of Chromatin Fibers. *J. Biomol. Struct. Dyn.* **16**:1097