



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29386

NBP2-29386 Protocol

Product Handling Protocol (NBP2-29386)

Product Handling Protocol (NBP2-29386): https://www.novusbio.com/products/chloroquine-inhibitor_nbp2-29386

1. To make 100 mM stock solution, dissolve 200 mg chloroquine in 3.9 ml water by gentle vortex.
2. Filter sterilize through a 0.22 μ m filter.
3. Store at 4°C (Note: Chloroquine solution is light sensitive).
4. For TLR signaling inhibition study, perform a pilot inhibitory test with the different concentrations of chloroquine ranging from 1 to 10 mM to optimize your experiments.