



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29385

NBP2-29385 Protocol

Product Handling Protocol (NBP2-29385)

Product Handling Protocol (NBP2-29385): https://www.novusbio.com/products/quinacrine-inhibitor_nbp2-29385

1. To make 100 mM stock solution, dissolve 200 mg quinacrine in 4.2 ml water by gentle vortex.
2. Filter sterilize through a 0.22 um filter.
3. Store at 4oC (Note: Quinacrine solution is light sensitive.).
4. For TLR signaling inhibition study, perform a pilot inhibitory test with the different concentrations of quinacrine ranging from 1 to 10 mM to optimize your experiments.