



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NBP2-26243](http://www.novusbio.com/NBP2-26243)

## NBP2-26243 Protocol

# Product Handling Protocol (NBP2-26243)

---

Product Handling Protocol (NBP2-26243): [https://www.novusbio.com/products/curcumin\\_nbp2-26243](https://www.novusbio.com/products/curcumin_nbp2-26243)

1. Add DMSO to bring curcumin to desired concentration; Solubility is at 11mg/ml
2. Dissolve curcumin in DMSO completely by gentle vortex.
3. Divide into useable aliquots and store them at -80C (Stock solutions are stable for up to 3 months at -80C).
4. Thaw stock solution briefly in a 37C water bath just prior to use.
5. Perform a pilot inhibitory assay with different concentrations of curcumin ranging from 5 to 100 M to optimize your experiments.