

Nile Red

Nile Red is a red fluorescent neutral lipid stain.



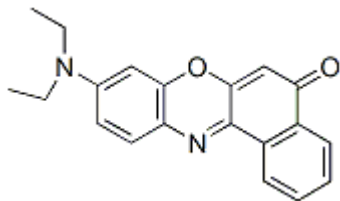
Product attributes

| | |
|-----------------------------|----------------|
| CAS number | 7385-67-3 |
| Probe cellular localization | Lipid droplets |
| Colors | Red |
| Excitation/Emission | 552/636 nm |

Product Description

Nile Red is used to localize and quantitate lipids, particularly neutral lipid droplets within cells. Nile Red is almost nonfluorescent in water and other polar solvents but undergoes fluorescence enhancement and large absorption and emission blue shifts in nonpolar environments. For lipid droplet dyes with improved specificity and more wavelength options compared to Nile Red, see our [LipidSpot™ Lipid Droplet Stains](#)

- Dark red solid soluble in DMF and DMSO
- $\lambda_{\text{Abs}}/\lambda_{\text{Em}}$ (MeOH) = 552/636 nm
- Store at 4 °C and protect from light
- $\text{C}_{20}\text{H}_{18}\text{N}_2\text{O}_2$
- MWt: 318.37
- [7385-67-3]



This datasheet was generated on January 17, 2026 at 09:40:00 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/nile-red/>