

Nile Red

Nile Red is a red fluorescent neutral lipid stain.



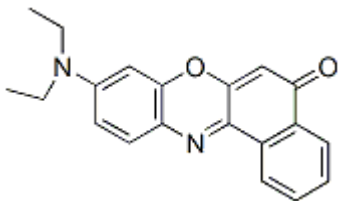
Product attributes

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
- $\lambda_{Abs}/\lambda_{Em}$ (MeOH) = 552/636 nm
- Store at 4 °C and protect from light
- C₂₀H₁₈N₂O₂M
- MWt: 318.37
- [7385-67-3]



This datasheet was generated on June 19, 2026 at 12:02:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/nile-red/>