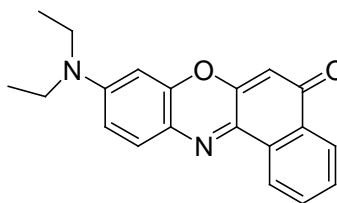


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Nile Red**CATALOG #** 60029**MOLECULAR INFORMATION:** C₂₀H₁₈N₂O₂
Mwt: 318.37**PROPERTIES:**

Color Dark red
Form Solid
Purity ≥ 95% by HPLC
Solubility Soluble in DMSO or DMF
Abs/Em: 552nm /636nm (methanol)

STORAGE AND HANDLING:

Store the solid at 4 °C and protect from light.

APPLICATION:

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.

Ref.: 1) *J Cell Biol* (1985), 100, 965

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.