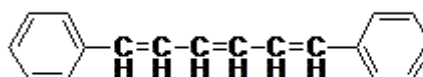


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 1,6-Diphenyl-1,3,5-hexatriene**CATALOG #** 60021**MOLECULAR INFORMATION:** C₁₈H₁₆
Mwt: 232.32**PROPERTIES:****Color** Yellow to tan**Form** Solid**Purity** ≥ 95% by HPLC**Solubility** Soluble in DMSO or DMF**STORAGE AND HANDLING:** Store the solid at -20 °C and protect from light.**APPLICATION:**

1,6-Diphenyl-1,3,5-hexatriene (DPH, excitation/emission=350 nm/420 nm) is almost non-fluorescent in water, but it exhibits a strong increase in fluorescence after intercalation into membranes and can be used as a probe for viscosity, polarity and lipid order. DPH has been used for the high-content imaging of neutral lipid droplets¹ as well as a reporter of inner spore membrane fluidity in bacterial spores².

Ref.: 1) *BioTechniques* (2011), doi 10.2144/000113702; 2) *J. Microbiol. Methods* (2013), <http://dx.doi.org/10.1016/j.mimet.2013.11.009>

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.