

1,6-Diphenyl-1,3,5-hexatriene

1,6-Diphenyl-1,3,5-hexatriene (DPH) 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.



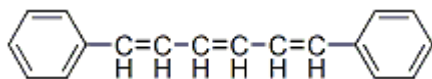
Product attributes

CAS number	1720-32-7
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Blue
Excitation/Emission	350/420 nm

Product Description

1,6-Diphenyl-1,3,5-hexatriene (DPH) 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.

- $\lambda_{Ex}/\lambda_{Em} = 350/420$ nm
- Yellow solid soluble in DMF and DMSO
- Store at -20 °C and protect from light
- C₁₈H₁₆
- MWt: 232.32
- [1720-32-7]



References

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