DiD

DiD (DiIC18(5); 1,1'-dioctadecyl-3,3,3',3'-tetramethylindodicarbocyanine, 4-chlorobenzenesulfonate salt) is a lipophilic carbocyanine dye similar to DiI, but with longer absorption and emission wavelengths.

ntötum 8

Product Description

DiD (DilC18(5); 1,1'-dioctadecyl-3,3,3',3'- tetramethylindodicarbocyanine, 4-chlorobenzenesulfonate salt) is a lipophilic carbocyanine dye similar to Dil, but with longer absorption and emission wavelengths. DiR, DiD, Dil and Neuro-DiO can be used in combination for multicolor imaging. DiD is also available in CellBriteTM Red, a ready-to-use cell labeling solution.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 644/663 nm
- ε= 193,000
- Dark blue solid soluble in ethanol, DMF or DMSO
- Store at 4°C and protect from light
- C₆₇H₁₀₃CIN₂O₃S
- MW: 1052.1
- [127274-91-3]

$$H_3C$$
 CH_3 H_3C CH_3 CH_3 CH_3 CH_3 CH_4 CH_5 CH_7 CH_8 CH_8

References

- 1. International Journal of Pharmaceutics (2014), doi: 10.1016/j.ijpharm.2014.11.012
- 2. Artificial Cells, Nanomedicine, and Biotechnology(2018), doi: 10.1080/21691401.2018.1445093

This datasheet was generated on November 15, 2025 at 03:19:00 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/did-diic185-or-11-dioctadecyl-3333-tetramethylindodicarbocyanine-4-chlorobenzenesulfonate-salt/

Product attributes

CAS number	127274-91-3
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
For live or fixed cells	For fixed cells, For live/intact cells
Assay type/options	Co-cultures, Extended staining (several days to weeks)
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
Colors	Far-red
Excitation/Emission	644/663 nm

Call us: 800-304-5357 Email: btinfo@biotium.com