

## DiD

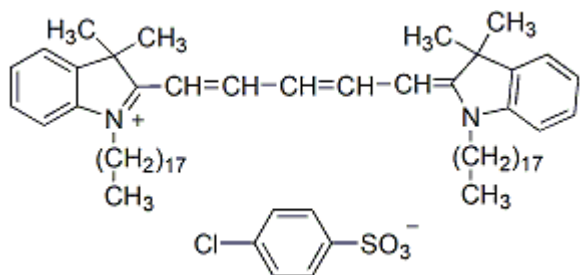
DiD (DiI18(5); 1,1'-dioctadecyl-3,3,3',3'-tetramethylindodicarbocyanine, 4-chlorobenzenesulfonate salt) is a lipophilic carbocyanine dye similar to DiI, but with longer absorption and emission wavelengths.



## Product Description

DiD (DiI18(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. DiR, DiD, DiI and Neuro-DiO can be used in combination for multicolor imaging. DiD is also available in [CellBrite™ Red](#), a ready-to-use cell labeling solution.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 644/663 nm
- $\epsilon$  = 193,000
- Dark blue solid soluble in ethanol, DMF or DMSO
- Store at 4 °C and protect from light
- $C_{67}H_{103}ClN_2O_3S$
- MW: 1052.1
- [127274-91-3]



## References

1. International Journal of Pharmaceutics (2014), doi: [10.1016/j.ijpharm.2014.11.012](https://doi.org/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 January 12, 2026 at 02:36:31 PM. 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