



Glowing products for science

## DiB

DiB is a lipophilic cationically charged blue fluorescent dye.



## Product Description

DiB is a lipophilic cationically charged blue fluorescent dye.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 353/442 nm
- Light yellow solid
- Soluble in DMSO, DMF, or EtOH
- MW: 1074

Note: For staining live or fixed cells, we recommend using [CellBrite™ Blue Cytoplasmic Membrane Staining Kit](#) (catalog #30024), which includes DiB solution and a cell loading buffer that is required for efficient cell staining.

## References

1. PloS One (2012) doi:10.1371/journal.pone.0041434

This datasheet was generated on December 6, 2025 at 09:40:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dib/>

Call us : [800-304-5357](tel:800-304-5357)

### Product attributes

<b>Probe cellular localization</b>	Membrane/cell surface, Membrane/vesicular
<b>For live or fixed cells</b>	For fixed cells, For live/intact cells
<b>Fixation options</b>	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
<b>Colors</b>	Blue
<b>Excitation/Emission</b>	353/442 nm