DilC1(5) (1,1'-dimethyl-3,3,3',3'-tetramethylindodicariodide)

rl stictum 3

The carbocyanine dye $DilC_1(5)$ has been used to measure mitochondrial membrane potential in apoptotic cells. Loss of membrane potential in apoptotic cells is reflected by a decrease of fluorescence signal in the infrared channel.

Product Description

The carbocyanine dye DilC₁(5) has been used to measure mitochondrial membrane potential in apoptotic cells. Loss of membrane potential in apoptotic cells is reflected by a decrease of fluorescence signal in the infrared channel.

- λ_{Ex}/λ_{Em} (MeOH) = 638/ 658nm
- Dark green solid soluble in DMSO or DMF
- Store at 4°C and protect from light
- C₂₇H₃₁IN₂
- MW: 510.46

References

1. Cytometry 33, 333 (1998).

This datasheet was generated on November 12, 2025 at 12:26:28 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/diic15-11-dimethyl-3333-tetramethylindodicarbocyanine-iodide/

Product attributes

Probe cellular localization	Mitochondria
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Cell permeability	Membrane permeant
Apoptosis/viability marker	Mitochondrial potential
Potential dependence	Mitochondrial potential-dependent
Colors	Far-red
Excitation/Emission	638/658 nm

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