

JC-1 (chloride salt)

A fluorescent mitochondrial dye that can be used for ratiometric detection of mitochondrial potential changes.

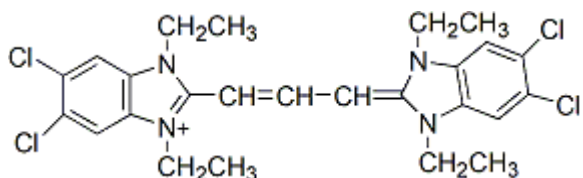


Product Description

JC-1 (5,5',6,6'-Tetrachloro-1,1',3,3'-tetraethylbenzimidazolylcarbocyanine chloride) is a mitochondrial dye that stains mitochondria in living cells in a membrane potential-dependent fashion. JC-1 monomer is in equilibrium with so-called J-aggregates, which are favored at higher dye concentration or higher mitochondrial membrane potential. The monomer JC-1 has green fluorescence ($\lambda_{Em} = 527 \text{ nm}$), while the J-aggregates have red fluorescence ($\lambda_{Em} = 590 \text{ nm}$). Therefore, it is possible to use the fluorescence ratio technique to study mitochondrial membrane potentials. JC-1 is particularly useful for apoptosis studies. In apoptotic cells, the dye stays in the cytoplasm and fluoresces green. It has also been used in high throughput drug screening applications.

While the original salt form of the dye is JC-1 iodide salt ([70011](#)), Biotium developed JC-1 chloride salt for applications where iodide ions are not desired. Biotium also offers [Aquaphile™ JC-1](#), an improved and more soluble formulation of the JC-1 dye which minimizes false positive J-aggregate signal outside of cells.

- $\lambda_{Ex}/\lambda_{Em}$ 510/527 nm (monomer); 585/590 nm (aggregate)
- ϵ (MeOH) = 190,000
- Red solid soluble in DMSO
- Store at 4 °C and protect from light, especially in solution
- $C_{25}H_{27}Cl_5N_4$
- MW: 561.5



References

1. PNAS 88, 3671 (1991).
2. Biochemistry 30, 4480 (1991).

This datasheet was generated on January 20, 2026 at 04:10:15 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/jc-1-chloride-salt-5566-tetrachloro-1133-tetraethylbenzimidazolylcarbocyanine-chloride/>

Product attributes

Probe cellular localization	Mitochondria
For live or fixed cells	For live/intact cells
Assay type/options	No-wash staining, Real-time imaging
Cell permeability	Membrane permeant
Apoptosis/viability marker	Mitochondrial potential
Potential dependence	Mitochondrial potential-dependent
Indicator type	Ratiometric
Colors	Green/Red
Excitation/Emission	510/527 nm (monomer); 585/590 nm (aggregate)