

DiOC5(3)

DiOC5(3) and DiOC6(3) are two of the most widely used carbocyanine dyes used for membrane potential measurements.



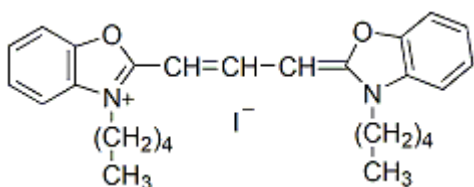
Product attributes

CAS number	53213-81-3
For live or fixed cells	For live/intact cells
Potential dependence	Slow response (translational) membrane potential dye
Colors	Green
Excitation/Emission	482/497 nm

Product Description

DiOC5(3) and DiOC6(3) are two of the most widely used carbocyanine dyes used for membrane potential measurements.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 482/497 nm
- Orange solid soluble in DMSO or DMF
- Store at 4 °C and protect from light
- $C_{27}H_{33}IN_2O_2$
- MW: 544.47
- [53213-81-3]



References

1. J Biol Chem 270, 3788 (1995).

This datasheet was generated on June 17, 2026 at 01:03:41 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/dioc53-33-dipentylloxycarbocyanine-iodide/>