Glowing products for science

## **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

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

## **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<sub>27</sub>H<sub>33</sub>IN<sub>2</sub>O<sub>2</sub>
- MW: 544.47
- [53213-81-3]

## References

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

This datasheet was generated on November 10, 2025 at 03:40:34 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/dioc53-33-dipentyloxacarbocyanine-iodide/">https://biotium.com/product/dioc53-33-dipentyloxacarbocyanine-iodide/</a>