

## DiO/DPA Membrane Potential Detection Kit



DiO/DPA is a highly sensitive FRET-based dye system that reports membrane potential change on submillisecond time scale.

### Product Description

DiO/DPA is a non-invasive optical membrane potential reporter system with low cytotoxicity, high speed, and high sensitivity.

The DiO/DPA system detects cytoplasmic membrane potential changes using the principle of fluorescence resonance energy transfer (FRET). The green fluorescent lipophilic dye DiO acts a stationary FRET donor in the plasma membrane. The anionic quencher DPA acts as a mobile FRET acceptor, rapidly partitioning in and out of the membrane in response to membrane potential, thus modulating DiO fluorescence. The DiO/DPA system has been reported to produce large fluorescence signal changes of >56% in HEK-293 cells and >25% in neuronal cultures and brain slices per 100 mV membrane potential change, and to report membrane potential *in situ* on submillisecond time scale (see [Ref. 1](#)).

The DiOC<sub>16</sub>(3) in this kit is an improved version over the DiOC<sub>16</sub>(3) from other suppliers; it is formulated with gluconate as the counter ion to increase dye solubility, for brighter and more uniform cell staining.

DPA is a hazardous material due to its explosive nature. To improve its safety during shipping and handling, it is provided as a dilute solution in the kit. We also offer DPA solid (catalog no. [60037](#)).

### References

1. J Neurosci (2009) 29(29), 9197. [DOI:10.1523/JNEUROSCI.1240-09.2009](https://doi.org/10.1523/JNEUROSCI.1240-09.2009)
2. Brit J Pharmacol (2011) 164, 667–680. [DOI:10.1111/j.1476-5381.2011.01396.x](https://doi.org/10.1111/j.1476-5381.2011.01396.x)

This datasheet was generated on June 19, 2026 at 01:37:23 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/diodpa-membrane-potential-detection-kit/>

### Product attributes

For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Potential dependence	Fast-response membrane potential dye
Colors	Green
Excitation/Emission	484/501 nm
CAS number	DiO: 34215-57-1, Dipicrylamine: 131-73-7
Storage Conditions	Store at room temperature, Protect from light