Glowing products for science

DiO/DPA Membrane Potential Detection Kit

Mintum

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 $DiOC_{16}(3)$ in this kit is an improved version over the $DiOC_{16}(3)$ 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 2. Brit J Pharmacol (2011) 164, 667–680. DOI:10.1111/j.1476-5381.2011.01396.x

This datasheet was generated on November 3, 2025 at 12:25:22 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/diodpa-membrane-potential-detection-kit/

Product attributes

Call us: 800-304-5357

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

Email: btinfo@biotium.com