

Product datasheet

Cellular Membrane Potential Assay Kit (Fluorometric - Orange) ab176764

[2 Images](#)

Overview

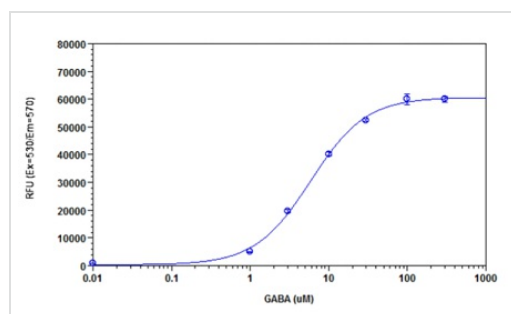
Product name	Cellular Membrane Potential Assay Kit (Fluorometric - Orange)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Membrane potential is the difference in voltage between the interior and exterior of a cell. The membrane potential allows a cell to function as a battery, providing power to operate a variety of "molecular devices" embedded in the membrane. In electrically excitable cells such as neurons, membrane potential is used for transmitting signals between different parts of a cell. Opening or closing of ion channels at one point in the membrane produces a local change in the membrane potential, which causes electric current to flow rapidly to other points in the membrane. Ion channels have been identified as important drug discovery targets.</p> <p>Abcam's Cellular Membrane Potential Assay Kit (Fluorometric - Orange) (ab176764) is a homogeneous assay with fast read time. It uses a proprietary long wavelength membrane potential indicator to detect the membrane potential change that is caused by the opening and closing of the ion channels. The red fluorescence of the membrane potential indicator used in the kit has enhanced fluorescence upon entering cells and minimizes the interferences resulted from the screening compounds and/or cellular autofluorescence.</p>
Platform	Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

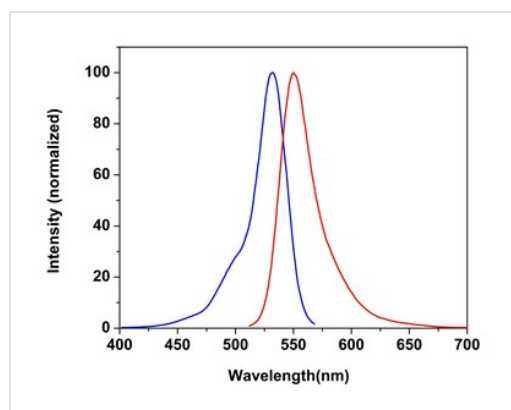
Components	1 x 96 tests	10 x 96 tests	100 x 96 tests
10X Assay buffer	1 x 1ml	1 x 10ml	1 x 100ml
HHBS	1 x 9ml	1 x 100ml	0 x 0ml
MP Sensor	1 x 15µl	1 x 150µl	10 x 150µl

Images



GABA Dose Response in WSS-1 cells measured with Cellular Membrane Potential Assay Kit (Fluorometric-Orange) (ab176764)

WSS-1 cells were seeded overnight in 50,000 cells per 100 μ L per well in a 96-well black wall/clear bottom plate. The cells were incubated with 100 μ l of [ab176765](#) for 30 minutes at room temperature. GABA (50 μ l/well) was added to achieve the final indicated concentrations.



Excitation and Emission Spectra for Cellular Membrane Potential Assay Kit (Fluorometric - Orange) (ab176764)

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors