# **Reagents for Determining Cellular Membrane Potentials**

| Ordering Information   | Storage Conditions                                   |
|--|--|
| Product Numbers: 21410 (25 mg), 21411 (25 mg), 21412 (5 mg), 21414 (25 mg) | Store at -20 °C, desiccated and protected from light |

# **Introduction**

DiBAC4(3), DiBAC4(5) and DiSBAC2(3) are a set of sensitive slow-response membrane potential probes that are widely used for measuring membrane potentials of many biological systems. In general, slow-response probes exhibit potential-dependent changes in their transmembrane distribution that are accompanied by a large fluorescence change. The magnitude of their optical responses is much greater than that of fast-response probes (typically a 1% fluorescence change per mV). Slow-response probes, which include cationic carbocyanines, rhodamines and anionic oxonols, are suitable for detecting changes in average membrane potentials of nonexcitable cells caused by respiratory activity, ion-channel permeability, drug binding and other factors.

# **Chemical and Physical Properties**

| Membrane Potential Indicators                                      | Catalog<br>Numbers | Excitation | Emission | Molecular<br>Weight | Solvent |
|--|--------------------|------------|----------|---------------------|---------|
| DiBAC4(3) [Bis-(1,3-dibutylbarbituric acid)trimethine oxonol]      | 21411              | 493 nm     | 516 nm   | 516.64              | DMSO    |
| DiBAC4(5) [Bis-(1,3-dibutylbarbituric acid)pentamethine oxonol]    | 21410              | 590 nm     | 616 nm   | 542.67              | DMSO    |
| DiSBAC30111  | 21412              | 535 nm     | 560 nm   | 366.42              | DMSO    |
| DiSBAC2(3) [Bis-(1,3-diethylthiobarbituric acid)trimethine oxonol] | 21414              | 535 nm     | 560 nm   | 436.55              | DMSO    |

# Assay Protocol with Membrane Potential (MP) Indicators (Example of DiSBAC2(3))

## **Brief Summary**

Prepare cells in growth medium  $\rightarrow$  Add dye-loading solution (100 µL/well for a 96-well plate or 25 µL/well for a 384-well plate)  $\rightarrow$  Incubate at RT or 37 °C for 30 min to 1 hour  $\rightarrow$  Read fluorescence intensity at Ex/Em = 490/525 nm (Cat. # 21411), 590/625 nm (Cat. # 21410) or 540/590 nm (Cat. # 21414)

Note: Following is our recommended protocol for live cells. It only provides a guideline, and should be modified according to your specific needs.

#### 1. Prepare cells:

- 1.1 For adherent cells: Plate cells overnight in growth medium at 40,000 to 80,000 cells/well/100 μL for 96well plates or 10,000 to 20,000 cells/well/25 μL for 384-well plates.
- 1.2 For non-adherent cells: Centrifuge the cells from the culture medium and then suspend the cell pellets in equal amount of HHBS and MP dye-loading solution (see Step 2.2 below) at 125,000 to 250,000 cells/well/100 μL for 96-well poly-D lysine plates or 30,000 to 60,000 cells/well/25 μL for 384-well poly-D lysine plates. Centrifuge the plates at 800 rpm for 2 minutes with brake off before the experiments. Note: Each cell line should be evaluated on an individual basis to determine the optimal cell density for the intracellular calcium mobilization.

#### 2. Prepare DiSBAC2(3) dye-loading solution (for 1 plate):

2.1 Prepare a 10 to 30 mM stock solution of DiSBAC2(3) in high-quality, anhydrous DMSO. The stock solution should be used promptly; any remaining solution need be aliquoted and frozen at ≤-20 °C. Note: Avoid repeated freeze-thaw cycles, and protect from light.

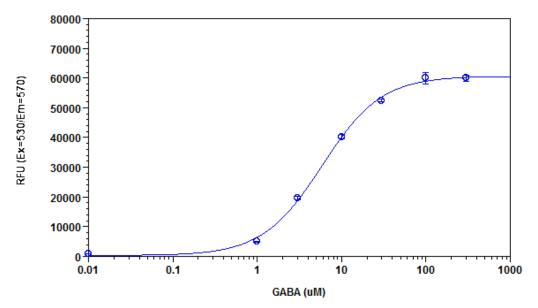
2.2 Prepare a 2X DiSBAC2(3) dye-loading solution: On the day of the experiment, either dissolve DiSBAC2(3) solid in DMSO or thaw an aliquot of the DiSBAC2(3) stock solution to room temperature. Prepare a 2X working solution of 20 to 40 μM in Hanks and 20 mM Hepes buffer (HHBS) or buffer of your choice, pH 7 with 0.04% to 0.08% Pluronic® F-127 (Cat. # 20053) and 2 mM Trypan Red Plus<sup>TM</sup> (Cat. # 2456). Mix them well by votexing. This working solution is stable for at least 2 hours at room temperature.

#### 3. Run Membrane Potential Assay:

- 3.1 Add 100 μL/well (96-well plate) or 25 μL/well (384-well plate) DiSBAC2(3) dye-loading solution (from Step 2.2) into the cell plate. *Note1: If your screen compounds interfere with growth medium and serum factors, replace the growth medium with equal volume of HHBS buffer before adding the DiSBAC2(3) dye-loading solution. Alternatively, cells can be grown in serum-free conditions. Note 2: Do NOT wash the cells after dye loading.*
- 3.2 Incubate the dye-loading plate in a cell incubator for 30 to 60 minutes. Note: In some cases, incubation at room temperature for 30 to 60 min may work better.
- 3.3 Prepare the compound plates by using HHBS or your desired buffer.
- 3.4 Run the membrane potential assay by monitoring the fluorescence intensity at Ex/Em = 540/590 nm (Cat. # 21414).

Note: It is important to run the signal test before your experiment. Different instruments have their own intensity range. Adjust the signal test intensity to the level of 10% to 15% of the maximum instrument intensity counts. For example, the maximum fluorescence intensity count for FLIPR-384 is 65,000, so the instrument settings should be adjusted to have its signal test intensity around 7,000 to 10,000.

### **Data Analysis**



**Figure 1.** GABA Dose Response was measured with DiSBAC2(3) in WSS-1 cells. WSS-1 cells were seeded overnight at 50,000 cells/100  $\mu$ L/well in a 96-well black wall/clear bottom costar plate. The cells were incubated with 100  $\mu$ L of DiSBAC2(3) dye loading solution for 30 minutes at room temperature. GABA (50  $\mu$ L/well) was added by FlexStation (Molecular Devices) to achieve the final indicated concentrations.

**Disclaimer:** This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.

©2008 by AAT Bioquest®, Inc., 520 Mercury Drive, Sunnyvale, CA 94085. Tel: 408-733-1055 Ordering: <u>sales@aatbio.com</u>; Tel: 800-990-8053 or 408-733-1055; Fax: 408-733-1304 Technical Support: <u>support@aatbio.com</u>; Tel: 408-733-1055