

Technical Data Sheet

Codex Non-Wash Ratiometric Calcium Dye Kit

Product Information

Catalog Number:	CB-80500-421
Size:	Reagents for 10 plates
Components:	Ratiometric Calcium Dye~50 µg, 10 vials, lyophilized (Part No: 80500-121) 10X Calcium Dye Signal Enhancer, 10 ml (Part No: 80500-112)

Description

The Codex Non-Wash Ratiometric Calcium Dye Kit allows homogeneous measurement of intracellular calcium changes caused by activation of G-protein coupled receptors or calcium channels. The ratio of 340/380 nm excitation allows accurate measurements of the intracellular Ca²⁺ concentration. Measuring by ratio considerably reduces the effects of uneven dye loading, leakage of dye, and photobleaching, as well as problems associated with measuring Ca²⁺ in cells of unequal thickness. The assay involves only one step of dye addition and does not require any washing steps. It is user friendly and cost effective. The assay can be easily implemented in a high throughput environment.

Storage

Ratiometric Calcium Dye~50 µg	-20°C (protected from light)
10X Calcium Dye Signal Enhancer	Room Temp.

Materials not included

DMSO	Sigma D4540
------	-------------

ASSAY PROTOCOL**Prepare the cell plate:**

1. Seed 80 µl of cell suspension into each well of a 96-well plate or 20 µl of cell suspension into each well of a 384-well plate.
2. Grow the cells overnight in a CO₂ incubator

Prepare the buffers (for one plate):

On the 2nd day:

1. Prepare Buffer A (1X HBSS with 20 mM HEPES):
10 ml of 1M HEPES, pH 7.3 + 490 ml of 1X HBSS
2. Prepare 1 ml of 500 mM Probenecid (optional).
Dissolve 142 mg of Probenecid in 1 ml of 1N NaOH
3. Prepare stock solution of calcium dye
Add 10 µl of DMSO into each vial containing calcium dye for 1 plate
4. Prepare **2X Dye Loading Buffer** (1 plates).

Add 1 ml of Codex 10X Calcium Dye Signal Enhancer into 9 ml of Buffer A.

Add 100 μ l of 500 mM Probenecid (optional).

Add 10 μ l of calcium dye stock solution. Mix well by vortexing.

Assay:

1. Take the cell plate out from the incubator.
2. Add same volume of **2X Dye Loading Buffer** into each well, 80 μ l to a 96-well plate or 20 μ l to a 384-well plate.
3. Incubate at 37 °C incubator for 1 hr.
4. Take the cells out of the incubator and leave at room temp (in the dark) for 30 min (optional).
5. Put the plate into the instrument for assay

To perform the assays, use the following wavelength parameters. Excitation: 340 nm; Emission: 510 nm and Excitation: 380 nm; Emission: 510 nm

Note. Dispense speed and height for compound additions need to be optimized for each instrument

Appendix

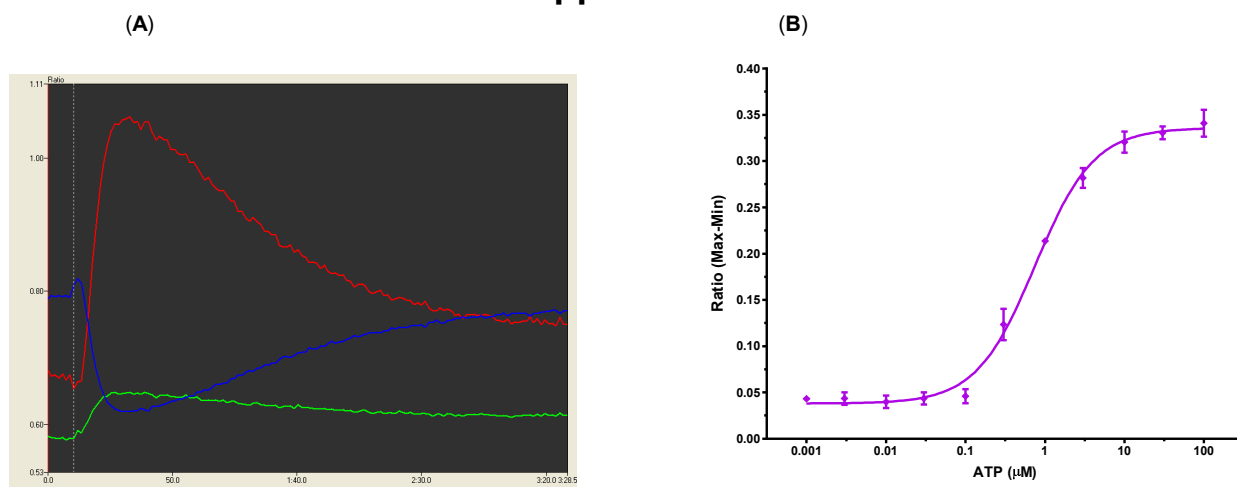


Figure 1. Response of endogenous P2Y receptors to ATP. HEK293 cells were plated overnight in 20 μ l culture medium on a 384 well black/clear plate. The next day, the cells were dye-loaded by adding 20 μ l of 2X Dye Loading Buffer and incubating for 1 hour at 37°C. ATP solution was added (10 μ l/well) by a FDSS 7000 (Hamamatsu), and the data was recorded simultaneously. **A.** Kinetic curve of calcium response to 100 μ M ATP (Green: emission at 340 nm; Blue: emission at 380 nm; Red: 340/380 ratio). **B.** ATP dose response curve (n = 4). EC50 = 0.75 μ M.

Limited Label License

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product for research use only by the buyer, where such research does not involve testing, analysis or screening services for any third party in return for compensation on a per test basis. The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data to third parties; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For more information, contact Business Development, Codex BioSolutions, Inc. 401 Professional Drive, Suite 160, Gaithersburg, MD 20878, USA. Tel: (240)-632-8810. Fax: (240)-632-8820

Codex Biosolutions, Inc. 401 Professional Drive, Suite 160, Gaithersburg, MD 20879
Tel: (240)-632-8810; Fax: (240)-632-8820 E-Mail: orders@codexbiosolutions.com
<http://www.codexbiosolutions.com/>