

Technical Data Sheet

Codex ACTOne™ Non-Wash Calcium Dye Kit

Product Information

Catalog Number:	CB-80500-311
Size:	Reagents for 100 plates
Components:	Codex Calcium Dye, 100 vials, lyophilized (Part No: 80500-110) 10X Calcium Dye Signal Enhancer, 100 ml (Part No: 80500-112)

Description

The Codex ACTOne™ Non-Wash Calcium Dye Kit allows homogeneous measurement for identifying cardiotoxic compounds using iPSC-CMs. It is a fast, simple, and reliable fluorescence-based method. The dye showed very little effect on beating rate and peak amplitude over a very long period. The dye can be used to quantify and report drug-induced early-after depolarization (EAD)-like waveforms, cardiomyocyte ectopic beats, changes in beating rate, and peak amplitude from a variety of agents. It is user friendly and cost effective. The assay can be easily implemented in a high throughput environment.

Storage

Codex Calcium Dye	-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. Human iPSC-CMs from Codex BioSolutions were plated in 96-well plates that had been pre-coated with fibronectin overnight at 37 °C. The cell density is 60,000 cells/well.
2. Cells were cultured for 24 h in plating media and, after that, switched to the maintenance media. The maintenance media was changed every other day and 24 h prior to the day of the experiment.
3. Cells were tested 2 weeks after plating at which time they formed a spontaneously beating monolayer.

Prepare the Dye Loading Solution:

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 stock solution of calcium dye
Add 8µl of DMSO into each well containing 50 µg of calcium dye
3. Prepare **2X Dye Loading Solution** (1 plates).
Add 0.8 ml of Codex 10X Calcium Dye Signal Enhancer into 7.2 ml of Buffer A.
Add 8 µl 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 Solution** into each well, 80 μ l to a 96-well plate
3. Incubate at 37 °C incubator for 1 hr.
3. Place the cell plates on a FDSS or FLIPR plate form and Ca²⁺ simultaneous transient signals were collected from 96-well plates. The standard filters for calcium assays (same as Fluo-4) is used

Appendix

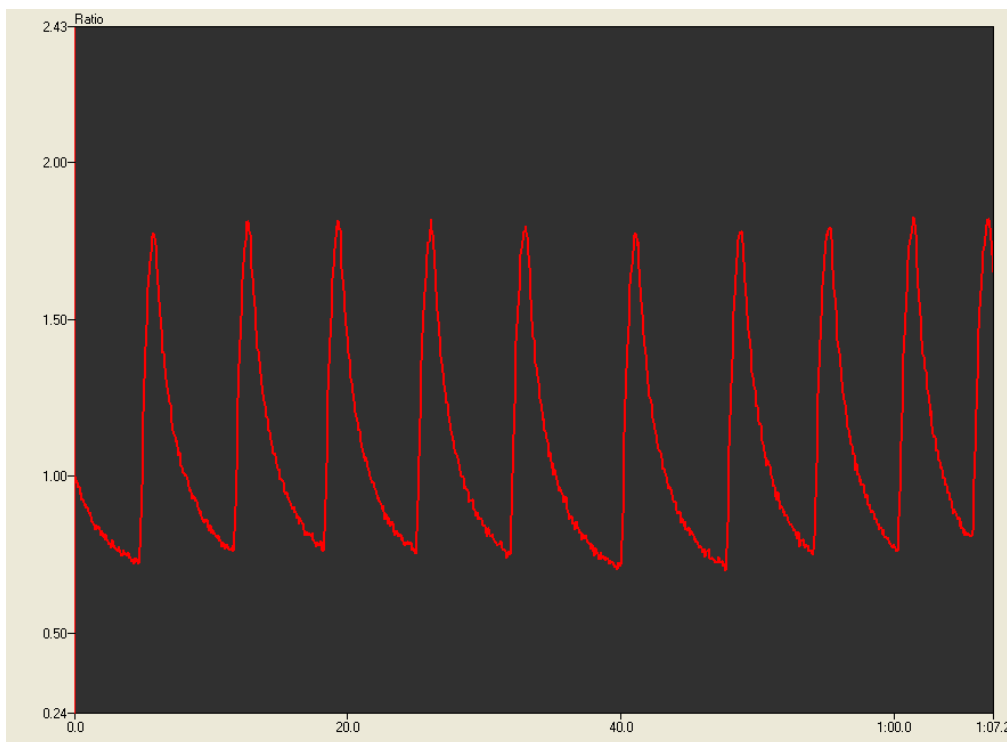


Figure 1. Human iPSC-CMs from Codex BioSolutions were plated in 96-well plates that had been pre-coated with fibronectin overnight at 37 °C. The cell density is 60,000 cells/well. Cells were cultured for 24 h in plating media and, after that, switched to the maintenance media. The maintenance media was changed every other day and 24 h prior to the day of the experiment. Cells were tested 2 weeks after plating at which time they formed a spontaneously beating monolayer. On experiment day, cells were incubated with Codex ACTOne® dye for 1 h at 37 °C, 5% CO₂. FDSS7000 imaging platform (Hamamatsu Ltd., Hamamatsu, Japan) was used to simultaneously collected Ca²⁺ transient signals from 96-well plates, at a sampling rate of 20 Hz.

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