



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### **TRPM2-HEK293 Recombinant Cell line Cat #: 90331**

#### **Description**

Recombinant HEK293 cell line expressing tetracycline-inducible human TRPM2 (transient receptor potential cation channel, subfamily M, member 2, accession number NM\_003307).

#### **Host cell**

HEK293 cells, tetracycline-inducible

#### **Format**

Each vial contains  $1.5 \times 10^6$  cells in 1 ml of 10% DMSO.

#### **Storage**

Immediately upon receipt, store in liquid nitrogen.

#### **Mycoplasma testing**

The cell line has been screened using the PCR-based Venor<sup>®</sup>GeM Mycoplasma Detection kit (Sigma-Aldrich) to confirm the absence of Mycoplasma species.

#### **Introduction**

TRPM2 channel belongs to the transient receptor potential channel (TRP) superfamily which is divided into seven main subfamilies. TRPM2 channel is a nonselective, calcium-permeable cation channel highly expressed in the brain, vascular smooth muscle, endothelial cells and immune cells. It is activated by oxidative stressors such as hydrogen peroxide, cyclic ADP ribose, and NADP, leading to an increase in the intracellular free calcium concentration and cell death. TRPM2 is a potential drug target in a variety of human diseases, including cardiovascular and neurodegenerative diseases.

#### **Applications**

- Monitor TRPM2 calcium channel activity
- Screen for activators or inhibitors of TRPM2 calcium channel

#### **Functional validation**

N<sup>-</sup> terminal FLAG tagged human TRPM2 channel has been stably integrated into HEK293 cells and its expression can be induced by tetracycline (doxycycline). The tetracycline-inducible expression of TRPM2 was confirmed by Western blotting.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

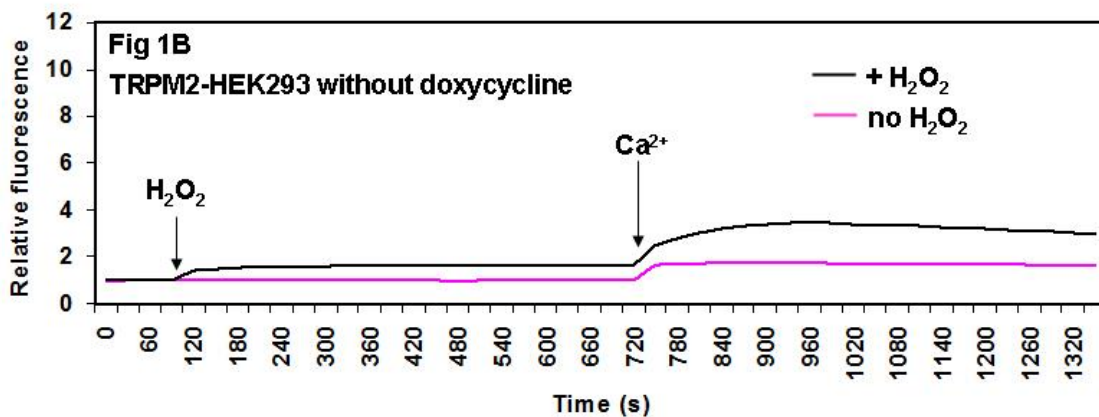
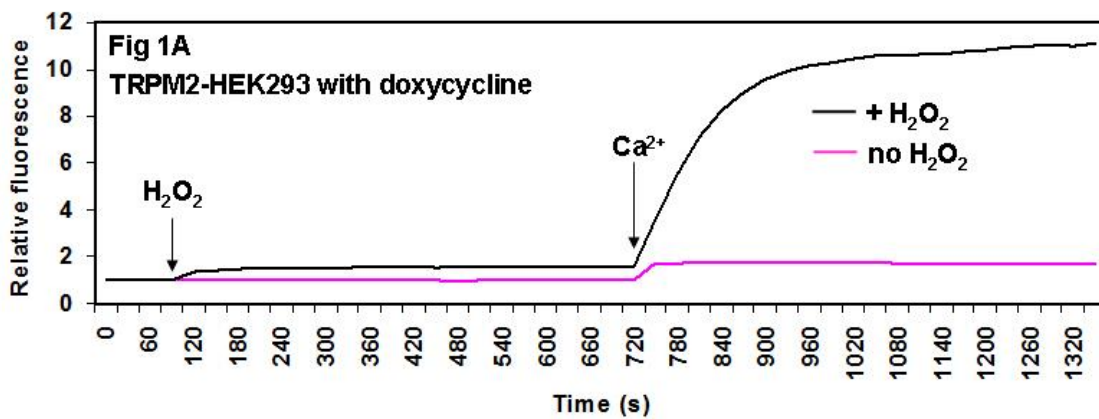
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

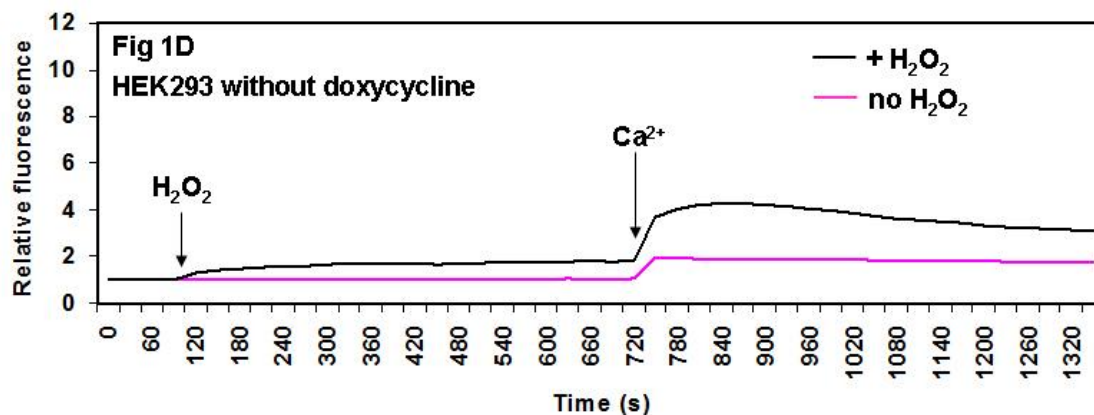
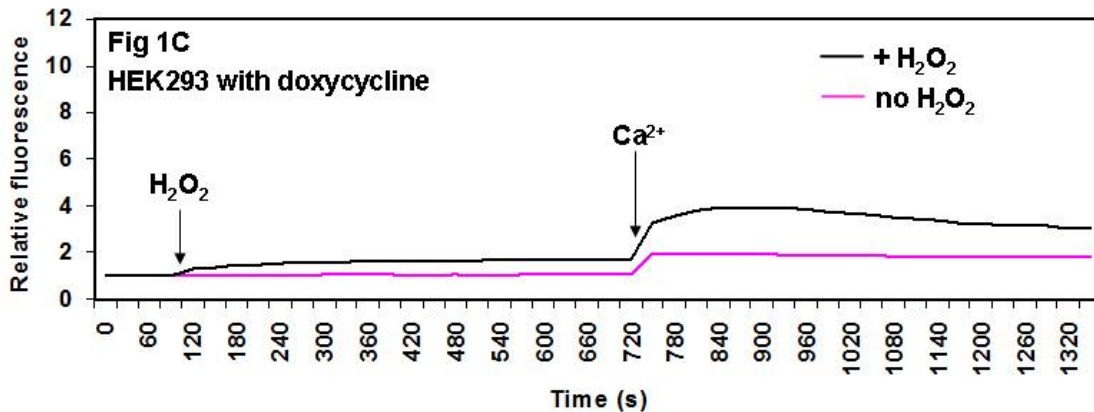
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

The function of TRPM2 was characterized by calcium assay. TRPM2-HEK293 cells were pre-treated with doxycycline to induce the expression of TRPM2, then stimulated with  $H_2O_2$ , and calcium influx was observed. The  $H_2O_2$ -induced calcium influx via TRPM2 could be blocked by N-(p-Amylcinnamoyl) anthranilic Acid (ACA), a TRP channel blocker.

**Figure 1. TRPM2 expression in HEK293 cells produced calcium influx after extracellular addition of  $H_2O_2$ .**



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.  
 To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



A) TRPM2-HEK293 cells with doxycycline treatment; B) TRPM2-HEK293 cells without doxycycline treatment; C) Control HEK293 cells (tetracycline-inducible) with doxycycline treatment; D) Control HEK293 cells (tetracycline-inducible) without doxycycline treatment.

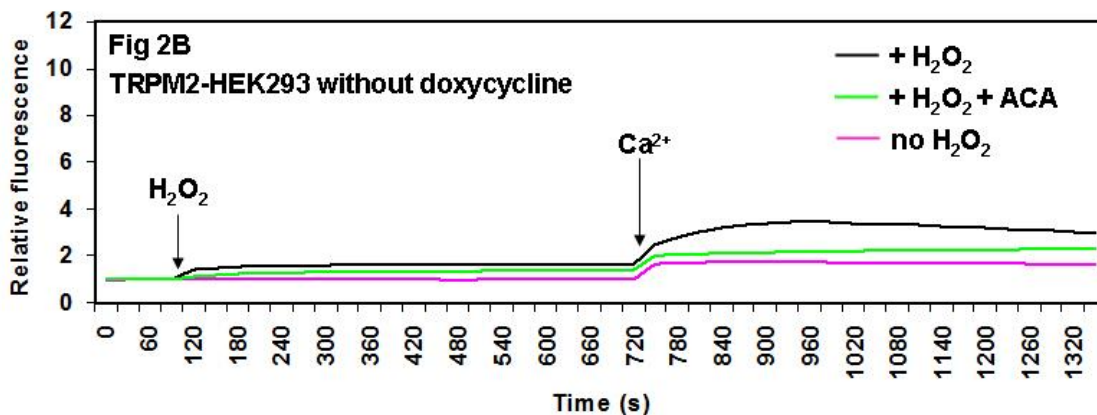
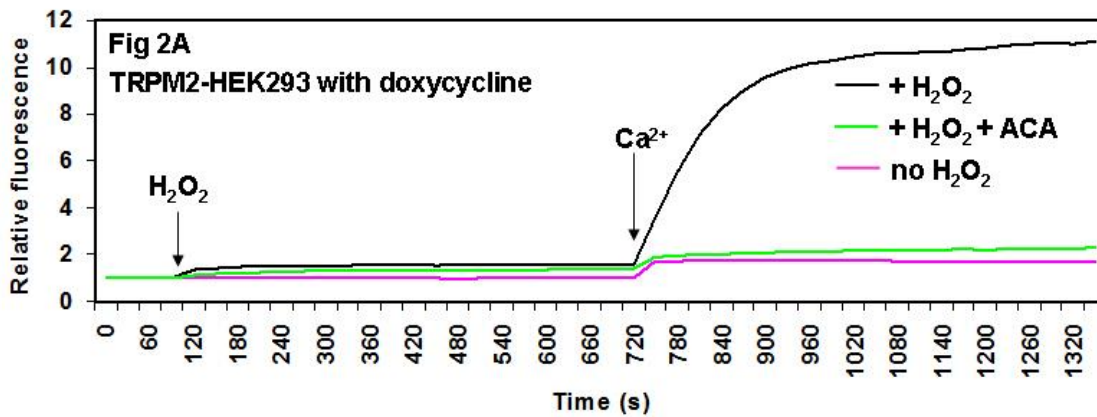
TRPM2-HEK293 or parental HEK293 cells (tetracycline-inducible) cells were pre-treated with doxycycline (A and C) or without doxycycline (B and D) for 24 hours. The next day cells were loaded with the fluorescent Ca<sup>2+</sup> indicator Fluo-8 and incubated in the absence of extracellular Ca<sup>2+</sup>, then treated with (black) or without (pink) H<sub>2</sub>O<sub>2</sub> (~ 6.7 mM). Following this, Ca<sup>2+</sup> (final 10 mM) was added to cells where indicated. The calcium influx was measured by Fluo-8 fluorescence (excitation 485±10nm and emission 528±10nm).

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.  
 To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Figure 2. H<sub>2</sub>O<sub>2</sub>-induced Ca<sup>2+</sup> influx in doxycycline-treated TRPM2-HEK293 cells was blocked by N-(p-Amylcinnamoyl) anthranilic Acid (ACA), a TRP channel blocker.**

A) TRPM2-HEK293 cells with doxycycline treatment; B) TRPM2-HEK293 cells without doxycycline treatment;

TRPM2-HEK293 cells were pre-treated with (A) or without (B) doxycycline for 24 hours. The next day cells were loaded with the fluorescent Ca<sup>2+</sup> indicator Fluo-8 and incubated with or without ACA (10 μM) in the absence of extracellular Ca<sup>2+</sup>, then treated with or without H<sub>2</sub>O<sub>2</sub> (~ 6.7 mM). Following this, Ca<sup>2+</sup> (final 10 mM) was added to cells where indicated. The calcium influx was measured by Fluo-8 fluorescence (excitation 485±10nm and emission 528±10nm).



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.  
 To place your order, please contact us by Phone 1.858.829.3082 Fax 1.858.481.8694  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Culture conditions

**Thaw Medium 1 (BPS Cat. #60187):** MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Invitrogen #26140-079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone SV30010.01)

**Complete Growth Medium:** Thaw Medium 1 (BPS Cat. #60187) plus 5 µg/ml Blastidicin (Life Technologies # R210-01), and 400 µg/ml of Geneticin (Life Technologies #11811031).

Cells should be grown at 37°C with 5% CO<sub>2</sub> using complete growth medium (Thaw Medium 1, Blastidicin, and Geneticin). It may be necessary to adjust the percentage of CO<sub>2</sub> in the incubator depending on the NaHCO<sub>3</sub> level in the basal medium. Cells should exhibit a typical cell division time of ~24 hours.

**To thaw the cells,** it is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, transfer to a tube containing 10 ml of Thaw Medium 1 (**no Blastidicin and Geneticin**), spin down the cells, resuspend cells in pre-warmed Thaw Medium 1 (**no Blastidicin and Geneticin**) in a T25 flask, and seed appropriate aliquots of cell suspension into new culture vessels. Culture at 37°C in a CO<sub>2</sub> incubator. At first passage switch to complete growth medium (**contains Blastidicin and Geneticin**). Cells should be split before they reach complete confluence.

**To passage the cells,** rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add complete growth medium and transfer to a tube. Spin down cells, resuspend cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ratio: 1:10 weekly.

## Induction of the target protein expression

Induce cells in MEM medium, 10% FBS, 1% non-essential amino acids, 1 mM Na-pyruvate, 1% Penicillin/Streptomycin, 0.2 µg/ml Doxycycline (MP Biomedicals #0219504401) for 24 hours before cell harvesting or assay.

## Vector and sequence

N'-terminal FLAG tagged human TRPM2 was cloned into a tetracycline regulated expression vector.

Polylinker: CMV-tetracycline operator (x 2)-BamHI-**TRPM2**-XhoI---SV40-neomycin<sup>R</sup>

hTRPM2 sequence (accession number NM\_003307)

MEPSALRKAGSEQEEGFEGLP RRVTDLGMVSNLRRS NSSLFKSWRLQCPFGNNDKQ  
ESLSSWIPENIKKKECVYFV ESKLSDAGKVVCQCGYTHEQHLEEATKPHTFQGTQWD  
PKKHVQEMPTDAFGDIVFTGLSQVKVKKYVRVSQDTPSSVIYHLMTQHWGLDVPNLLISV

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** info@bpsbioscience.com

TGGAKNFNMKPRLLKSIFRRGLVKVAQTTGAWIITGGSHTGVMKQVGEAVRDFSLSSSY  
KEGELITIGVATWGTVHRREGLIHPTGSFPAEYILDEDGQGNLTCLDSNHSHFILVDDGT  
HGQYGV EIPLRTRLEKFISEQTKERGGVAIKIPIVCVVLEGGPGTLHTIDNATTNGTPCVV  
VEGSGRVADVIAQVANLPVSDITISLIQKLSVFFQEMFETFTESRIVEWTKKIQDIVRRR  
QLLTVFREGKDGQDQVDVAILQALLKASRSQDHFHGHENWDHQLKLA VAWNRVDIARS  
EIFMDEWQWKPSDLHPTMTAALISNKPEFVKLFLENGVQLKEFVTWDTLLYLYENLDPS  
CLFHSLKQLKVLVEDPERPACAPAAPRLQMHHVAQVLRRELLGDFTQPLYPRPRHNDRLR  
LLLVPVHVKLVNQGVSLRSLYKRSSGHVFTTMDPIRDLLIWAIVQNRRELAGIWAQSQD  
CIAAALACSKILKELSKEEEDTDSSEMLALAEYEHRAIGVFTECYRKDEERAQKLLTR  
VSEAWGKTTCLQLALEAKDMKFVSHGGIQAFLTKVWWGQLSVDNGLWRVTLCMLAFP  
LLLTLGLISFREKRLQDVGTAAARARAFFTAPVVVFHLNLSYFAFLCLFAYVLMVDFQPV  
PSWCECAIYLWLFSLVCEEMRQLFYDPDECGLMKKAALYFSDFWNKLDVGAILLFVAG  
LTCRLIPATLYPGRVILSLDFILFCLRLMHIFTISKTLGPKIIIVKMMMVDVFFFLAVVWV  
SFGVAKQAILIHNERRVDWLFRGAVYHSYLTIFGQIPGYIDGVNFNPEHCSPNGTDPYK  
PKCPESDATQQRPAFPEWLTVLLLCLYLLFTNILLNLLIAMFNFTFQQVQEHTDQIWK  
QRHDLIEEYHGRPAAPPPFILLSHLQLFIKRVVLTAKRHKQLKNKLEKNEEAALLSWEI  
YLKENYLQNRQFQQKQRPEQKIEDISNKVDAMVDLLDPLKRSMSMEQRLASLEEQV  
AQTARALHWIVRTLRSAGFSSEADVPTLASQKAAEPPDAEPGGRKKTEEPGDSYHVNA  
RHLLYPNCPVTRFPVPNEKVPWETEFLIYDPPFYTAERKDAAMDPMGDTLEPLSTIQY  
NVVDGLRDRRSFHGPYTVQAGLPLNPMGRTGLRGRGSLSCFGPNHTLYPMVTRWRR  
NEDGAICRKSIIKMLEVLVVKLPLSEHWALPGGSREPGEMLPKRLKRILRQEHWPSFE  
NLLKCGMEVYKGYMDDPRNTDNAWIETVAVSVHFQDQNDVELNRLNSNLHACDSGAS  
IRWQVVDRIPLYANHKTLQKAAAEFGAHY

## References

1. Moran, M.M., *et al.* (2011) Transient receptor potential channels as therapeutic targets. *Nat. Rev. Drug Discov.* **10(8)**:601-620
2. Song, Y., *et al.* (2008) Development and validation of a cell-based high-throughput screening assay for TRPM2 channel modulators. *J. Biomol. Screen.* **13(1)**:54-61
3. Fonfria, E., *et al.* (2004) TRPM2 channel opening in response to oxidative stress is dependent on activation of poly(ADP-ribose) polymerase. *Br. J. Pharmacol.* **143(1)**:186-192

## License Disclosure

Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit use of this cell line in humans or for therapeutic or drug use. The license does not permit modification of the cell line in any way.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. BPS does not warrant the suitability of the cell line for any particular use, and does not accept any liability in connection with the handling or use of the cell line. Modifications of this cell line, transfer to another facility, or commercial use of the cells may require a separate license and additional fees; contact [sales@bpsbioscience.com](mailto:sales@bpsbioscience.com) for details. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.  
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)