

**Human Recombinant ACE2 Stable Cell Line**  
**Cat. No. M00770****Version 04242020**

---

I Introduction .....	1
II Background.....	1
III Representative Data.....	2
IV Thawing and Subculturing.....	2
V References .....	3
Limited Use License Agreement.....	4

---

**I. INTRODUCTION**

Catalog Number: M00770

Cell Line Name: HEK293/ACE2

Gene Synonyms: Angiotensin-converting enzyme 2, ACEH

Expressed Gene: Codon Optimized from NM\_021804.2; no expressed tags

Host Cell: HEK293

Quantity: Two vials of frozen cells (>1×10<sup>6</sup> cells/vial)

Stability: 15 passages

Application: Binding assay or use as immunogen

Freeze Medium: 95% Culture Medium, 5% (V/V) DMSO

Complete Growth Medium: DMEM, 10% FBS

Culture Medium: DMEM, 10% FBS, 250 µg/ml Hygromycin

Mycoplasma Status: Not detected\*

Storage: Liquid nitrogen immediately upon receipt

*\* The mycoplasma test was performed with MycoAlert™ PLUS Mycoplasma Detection Kit of Lonza.*

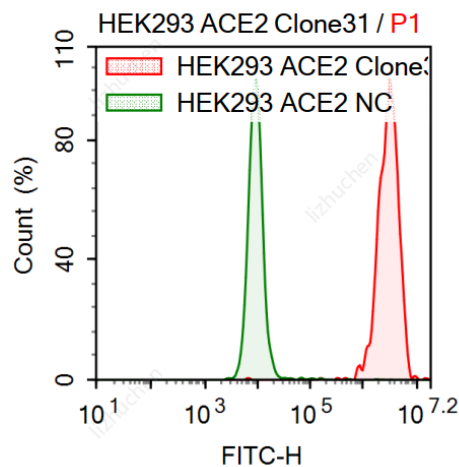
**II. BACKGROUND**

Angiotensin-converting enzyme 2 (ACE2), discovered as a homologue of ACE, acts as its physiological counterbalance providing homeostatic regulation of circulating angiotensin II (Ang II) levels. ACE2 is a zinc metalloenzyme and carboxypeptidase located as an ectoenzyme on the surface of endothelial and other cells. While its primary substrate appears to be Ang II, it can

hydrolyze a number of other physiological substrates. Additionally, ACE2 functions in other noncatalytic cellular roles including the regulation of intestinal neutral amino acid transport. It also serendipitously acts as the receptor for the severe acute respiratory syndrome virus. Upregulation of ACE2 expression and function is increasingly recognized as a potential therapeutic strategy in hypertension and cardiovascular disease, diabetes, lung injury, and fibrotic disorders. ACE2 is regulated at multiple levels including transcriptional, posttranscriptional (miRNA and epigenetic), and posttranslational through its shedding from the cell surface.

### III. REPRESENTATIVE DATA

#### Protein Expression Validation



**Figure 1.** FACS analysis of ACE2 expression in HEK293/ACE2 clone 31.

### IV. THAWING AND SUBCULTURING

#### Thawing Protocol

1. Remove the vial from liquid nitrogen tank and thaw cells quickly in a 37°C water-bath.
2. Just before the cells are completely thawed, decontaminate the outside of the vial with 70% ethanol and transfer the cells to a 15 ml centrifuge tube containing 9 ml of complete growth medium.
3. Pellet cells by centrifugation at 800 rpm for 4 min, and remove the medium.
4. Resuspend the cells in complete growth medium.
5. Transfer the cell suspension to a 10 cm dish with 10 ml of complete growth medium.
6. Grow the cells in incubator with 37°C, 5 % CO<sub>2</sub>.
7. Add antibiotic the following day.

#### Sub-culturing Protocol

1. Remove the culture medium from cells.
2. Wash cells with PBS (pH=7.4) to remove all traces of serum that contains trypsin inhibitor.
3. Add 2.0 ml of 0.25% (w/v) Trypsin- EDTA (GIBCO, Cat No. 25200-072) solution into 10 cm dish and observe the cells under an inverted microscope until cell layer is dispersed (usually within 3 to 5 minutes).  
Note: To avoid cells clumping, do not agitate the cells by hitting or shaking the dish while waiting for the cells detach. If cells are difficult to detach, please place the dish in 37°C incubator for ~2 min.
4. Add 6.0 to 8.0 ml of complete growth medium into dish and aspirate cells by gently pipetting.
5. Centrifuge the cells at 800 rpm for 4 min, and remove the medium.
6. Resuspend the cells in culture medium and add the cells suspension to new culture dish.
7. Grow the cells in incubator with 37°C, 5% CO<sub>2</sub>.

Subcultivation Ratio: 1:3 to 1:6.

Medium Renewal: Every 2 to 3 days

## V. REFERENCES

1. Anthony J. Turner, *The Protective Arm of the Renin Angiotensin System (RAS)*, 2015
2. Stertz S, *The intracellular sites of early replication and budding of SARS-coronavirus*, 2007

## Limited Use License Agreement

This is a legal agreement between you (Licensee) and GenScript USA Inc. governing use of GenScript's stable cell line products and protocols provided to licensee. By purchasing and using the stable cell line, the buyer agrees to comply with the following terms and conditions of this label license and recognizes and agrees to such restrictions:

- 1) The products are not transferable and will be used at the site where they were purchased. Transfer to another site owned by buyer will be permitted only upon written request by buyer followed by subsequent written approval by GenScript.
- 2) The purchaser 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.
- 3) The products sold by GenScript are for laboratory and animal research purposes only. The products are not to be used on humans, for consumption, or for any unlawful uses.

GenScript USA Inc. will not assert against the buyer a claim of infringement of patents owned or controlled by GenScript USA Inc. and claiming this product based upon the manufacture, use or sale of a clinical diagnostic, therapeutic and vaccine, or prophylactic product developed in research by the buyer in which this product or its components has been employed, provided that neither this product nor any of its components was used in the manufacture of such product. For information on the use of this product for other purposes, contact Marketing Department, GenScript USA Inc., 120 Centennial Avenue, Piscataway, New Jersey 08840, U.S.A. Phone: 1-732-885-9188. Fax: 1-732-210-0262. Email: [marketing@genscript.com](mailto:marketing@genscript.com).

**GenScript USA Inc,**  
860 Centennial Ave.  
Piscataway, NJ 08854  
Toll-Free: 1-877-436-7274  
Tel: 1-732-885-9188, Fax: 1-732-210-0262  
Email: [product@genscript.com](mailto:product@genscript.com)  
Web: <http://www.genscript.com>  
**For Research Use Only.**