

# hACE2-Fc

Soluble human Angiotensin-Converting Enzyme 2 (ACE2) protein fused to a human IgG1 Fc tag  
Catalog code: fc-hace2

<https://www.invivogen.com/human-ace2-proteins>

For research use only, not for diagnostic or therapeutic use

Version 20F23-NJ

## PRODUCT INFORMATION

### Contents:

- 50 µg of lyophilized hACE2-Fc protein
- 1.5 ml endotoxin-free water

### Protein construction:

Human ACE2 extracellular domain [M1-S740] with a human IgG1 Fc tag in C-terminus

**Accession sequence:** AAQ89076

**Species:** Human

**Tag:** C-terminal human IgG1 Fc

**Total protein size:** 967 a.a. (secreted form)

**Molecular weight:** ~130 kDa

**Purification:** protein G affinity chromatography

**Purity:** >95% (SDS-PAGE)

### Formulation:

0.2 µm filtered solution in a sodium phosphate buffer with glycine, saccharose, and stabilizing agents

### Storage:

- Product is shipped at room temperature. Store lyophilized product at -20°C. Lyophilized product is stable for at least 1 year.
- Reconstituted protein is stable for 1 month when stored at 4°C and for 1 year when aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles.

### Quality control:

- The size and purity of the protein has been confirmed by SDS-PAGE.
- hACE2-Fc has been functionally validated by ELISA using a coated Spike-RBD-His fusion protein
- Absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and TLR4 cellular assays.

## BACKGROUND

Human ACE2 (angiotensin I-converting enzyme-2) is a type I surface transmembrane protein expressed in arteries, heart, kidneys, and epithelia of the lung and small intestine<sup>1,2</sup>. It plays a critical role in the pathogenesis of the Coronavirus Disease-19 (COVID-19) caused by the severe acute respiratory syndrome coronavirus-2 (SARS-CoV-2). Indeed, hACE2 is now established as being the receptor for the Spike (S) protein of SARS-CoV and SARS-CoV-2, facilitating viral entry into target cells<sup>3-5</sup>. The blockade of ACE2 and the delivery of an excessive soluble form of ACE2 are among the strategies being investigated to treat COVID-19<sup>5,6</sup>. Of note, treatment with a soluble form of ACE2 not only slows viral entry into target cells, but also rescues the endogenous cellular ACE2 activity, protecting the lung from injury<sup>6</sup>.

## PRODUCT DESCRIPTION

hACE2-Fc is a soluble human ACE2 fusion protein generated by fusing the ACE2 extracellular domain (a.a. 1-740) to a C-terminal human IgG1 Fc tag with a TEV (Tobacco Etch Virus) sequence linker. This fusion protein has a molecular weight of ~130 kDa on a SDS-PAGE gel. hACE2-Fc has been generated by recombinant DNA technology, produced in CHO cells, and purified by protein G affinity chromatography.

## APPLICATIONS

- SARS-CoV and SARS-CoV-2 neutralization assays.
- Screening of small molecule inhibitors or of neutralizing antibodies able to block Spike-RBD and ACE2 interaction.

## METHODS

### hACE2-Fc resuspension (100 µg/ml)

*Note: Ensure you see the lyophilized pellet before resuspension.*

- Add 500 µl of endotoxin-free water to the vial and gently pipette until completely resuspended.
- Prepare aliquots and store at -20°C or 4°C.

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

## PROTEIN SEQUENCE

MSSSSWLLLLSLVAVTAAQSTIEEQAKTFLDKFNHE  
AEDLFYQSSLASWNYNTNITEENVQNMNNAQDK  
WSAFLKEQSTLAQMYPLQEIQNLTVKLQLQALQQ  
NGSSVLSSEDKSKRLNTILNTMSTIYSTGKVCNPDN  
PQECLELLEPGLNEIMANSLDYNERLWAWESWRSE  
VGKQLRPLYYEYVVLKNEMARANHYEDYGDYWR  
GDYEVNGVDGYDYSRGLIEDVEHTFEEIKPLYE  
HLHAYVRAKLMNAYPSYISPIGCLPAHLLGDMWG  
RFWNTNLYSLTVPFGQKPNIDVTDAMVDQAWDAQ  
RIFKEAEKFFVSVGLPNMTQGFWENSMLTDPGNV  
QKAVCHPTAWDLGKGDFRILMCTKVTMDDFLTA  
HHEMGHIQYDMAYAAQPFLLRNGANEGFHEAVG  
EIMSLSAATPKHLKLSIGLLSPDFQEDNETEINFLL  
KQALTI VGTLPFTYMLEKWRWMVFKGEIPKQW  
MCKWVEMKREIVGVVEPVPHDETYCDPASLFHV  
SNDYSFIRYYTRTLTYQFQFQEAALCQAAKHEGPLH  
KCDISNSTEAGQKLFNMLRRLGKSEPWTALLENVV  
GAKNMNVRPLLNYFEPLFTWLKDKQNKNSFVGWS  
TDWSPYADQSIKVRISLKSALGDKAYEWNNDNEM  
LFRSSVAYAMRQYFLKVKKNQMILFGEEDVRVANL  
KPRISFNFFVTAPKNVSDIIPRTEVEKAIRMSRSRI  
NDAFRLNDNSLEFLGIQPTLGPNNQPPVSRTE  
NLYFQGSSEPKSSDKTHTCPCPAPEAEGGSPVFL  
FPPKPKDQLMISRTPEVTCVVVDVSHEDPEVKFN  
WYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLH  
QDWLNGKEYKCKVSNKALPASIEKTISKAKGQPR  
EPQVYTLPPSREEMTKNQVSLTCLVKGFYPSDIAV  
EWESNGQPENNYKTTPPVLDSDGSFFLYSKLTVD  
KSRWQQGNVFSQSVLHEALHNHYTQKLSLSLSPGK

Green: signal sequence

Blue: ACE2 sequence

Black: TEV cleavage sequence

Red: Human IgG1 Fc sequence

## REFERENCES

1. Li F., 2016. Structure, function, and evolution of coronavirus spike proteins. *Annu. Rev. Virol.* 3:237-261. 2. Li F. *et al.*, 2005. Structure of SARS coronavirus spike receptor-binding domain complexed with receptor. *Science.* 309:1864-1868. 3. Wang N. *et al.*, 2020. Subunit vaccines against emerging pathogenic human coronaviruses. *Front. Microbiol.* 11:298. DOI: 10.3389/fmicb.2020.00298. 4. Padron-Regalado E., 2020. Vaccines for SARS-CoV-2: Lessons from other coronavirus strains. *Infect. Dis. Ther.* DOI: 10.1007/s40121-020-00300-x. 5. Monteil V. *et al.*, 2020. Inhibition of SARS-CoV-2 infections in engineered human tissues using clinical-grade soluble human ACE2. *Cell.* 181:1-9. 6. Zhang *et al.*, 2020. Angiotensin-converting enzyme 2 (ACE2) as a SARS-CoV-2 receptor: molecular mechanisms and potential target. *Intensive Care Medicine.* 46(4):586-590.

## RELATED PRODUCTS

Product	Catalog Code
Anti-Spike-RBD-hIgG1	srbd-mab1
Anti-Spike-RBD-hIgM	srbd-mab5
Anti-Spike-RBD-hIgA2	srbd-mab6
Spike-RBD-Fc	fc-sars2-srbd
Spike-RBD-His	his-sars2-srbd
Spike-S1-Fc	fc-sars2-s1
Spike-S1-His	his-sars2-s1
pDUO2-hACE2-TMPRSS2a	pduo2-hace2tpsa

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)