

pUNO1Fc-SARS2-N

Plasmid designed for the production of the SARS-CoV-2 Nucleocapsid::Fc fusion protein

Catalog code: p1fc-cov2-n

<https://invivogen.com/sars2-nucleocapsid-tag-production-vectors>

For research use only

Version 20H28-ED

PRODUCT INFORMATION

Contents

- 20 µg of lyophilized plasmid DNA
- 2 x 1 ml Blasticidin at 10 mg/ml

Storage and Stability

- Product is shipped at room temperature.
- Lyophilized DNA should be stored at -20°C.
- Resuspended DNA should be stored at -20°C and is stable for at least 1 year.
- Store Blasticidin at 4°C or -20°C. The expiry date is specified on the product label.

Quality control

- After purification by ion exchange chromatography, predominant supercoiled conformation is verified by electrophoresis.
- Plasmid construct is confirmed by restriction analysis and full-length open reading frame (ORF) sequencing.

METHODS

• Plasmid resuspension

- Quickly spin the tube containing the lyophilized plasmid to pellet the DNA.
- To obtain a plasmid solution at 1 µg/µl, resuspend the DNA in 20 µl of sterile water.
- Store resuspended plasmid at -20°C.

• Plasmid amplification and cloning

Plasmid amplification and cloning can be performed in *E. coli* GT116 or other commonly used laboratory *E. coli* strains, such as DH5α.

• Blasticidin usage

Blasticidin should be used at 25-100 µg/ml in bacteria and 1-30 µg/ml in mammalian cells.

PLASMID FEATURES

SARS-CoV-2 Nucleocapsid::hlgG1-Fc cassette

- **SV40 enhancer** is comprised of a 72-base-pair repeat and allows the enhancement of gene expression in a wide range of hosts. The enhancement varies from 2-fold in non-permissive cells to 20-fold in permissive cells. Furthermore, the SV40 enhancer is able to direct nuclear localization of plasmids¹.
- **EF-1α/HTLV composite promoter** is a composite promoter comprised of the Elongation Factor-1α (EF-1α) core promoter² and the 5' untranslated region of the Human T-Cell Leukemia Virus (HTLV)³. EF-1α utilizes a type 2 promoter that encodes for a «house keeping» gene. It is expressed at high levels in all cell cycles and lower levels during G0 phase. The promoter is also non-tissue specific; it is highly expressed in all cell types. The R segment and part of the U5 sequence (R-U5') of the HTLV Type 1 Long Terminal Repeat has been coupled to the EF-1α promoter to enhance stability of DNA and RNA. This modification not only increases steady state transcription, but also significantly increases translation efficiency possibly through mRNA stabilization.

• SARS-CoV-2 Nucleocapsid ORF

Nucleocapsid (N) is a phosphoprotein that associates with the viral RNA genome and forms the ribonucleoprotein core. N has a number of distinct domains, including an N-terminal RNA-binding domain, which captures the RNA, a C-terminal dimerization domain, which acts as an anchor to the replication-transcription complexes (RTCs) and, an intrinsically disordered central Ser/Arg (SR)-rich linker for primary phosphorylation^{4,5}. The N protein is highly immunogenic and abundantly expressed during infection, and is thus, capable of inducing protective immune responses against SARS-CoV-2. The pUNO1Fc-SARS2-N plasmid contains the nucleocapsid coding sequence from the Wuhan-Hu-1 isolate, with optimized signal sequence and codon usage.

• **TEV sequence** (Glu-Asn-Leu-Tyr-Phe-Gln-(Gly/Ser)) is cleavable by the TEV (Tobacco Etch Virus) protease between Gln and Gly/Ser residues.

• **Fc** is the Fc region of human IgG1 cloned at the C-terminus of the nucleocapsid and is followed by a stop codon. This tag comprises the CH2 and CH3 domains of the IgG heavy chain and the hinge region. The hinge serves as a flexible spacer between the two parts of the Fc fusion protein allowing each part of the molecule to function independently.

• **SV40 pAn** is the Simian Virus 40 late polyadenylation (pAn) signal enables efficient cleavage and polyadenylation reactions, resulting in high levels of steady-state mRNA⁶.

Antibiotic Selection cassette

• **hCMV (human cytomegalovirus) enhancer & promoter** drives the expression of the blasticidin resistance in mammalian cells.

• **EM7** is a bacterial promoter that enables the constitutive expression of the antibiotic resistance gene in *E. coli*.

• **bsr (blasticidin resistance gene)** from *Bacillus cereus* encodes a deaminase that confers resistance to the antibiotic blasticidin. The *bsr* gene is driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. Therefore, blasticidin can be used to select stable mammalian cells transfectants and *E. coli* transformants.

• **Human β-Globin pAn** is a strong polyadenylation (pAn) signal placed downstream of *bsr*. The use of beta-globin pAn minimizes interference⁷ and possible recombination events with the SV40 pAn signal.

General features of pUNO1Fc-SARS2-N

- **pMB1 ori** is a minimal *E. coli* origin of replication.

REFERENCES

1. Dean DA. *et al.* 1999. Sequence requirements for plasmid nuclear import. *Exp. Cell. Res.* 253:713-22.
2. Kim D. *et al.*, 1990. Use of the human elongation factor 1α promoter as a versatile and efficient expression system. *Gene* 91(2):217-23.
3. Takebe Y. *et al.*, 1988. SR alpha promoter: an efficient and versatile mammalian cDNA expression system composed of the simian virus 40 early promoter and the R-U5 segment of human T-cell leukemia virus type 1 long terminal repeat. *Mol Cell Biol.* 8(1):466-72.
4. Kang, S. *et al.* 2020. Crystal structure of SARS-CoV-2 nucleocapsid protein RNA binding domain reveals potential unique drug targeting sites. *Acta Pharm Sin B.*
5. Khan, M.T. *et al.* 2020. SARS-CoV-2 nucleocapsid and Nsp3 binding: an in silico study. *Arch Microbiol.*
6. Carswell S. & Alwine J., 1989. Efficiency of utilization of the simian virus 40 late polyadenylation site: effects of upstream sequences. *Mol Cell Biol.* 9(10):4248-58.
7. Yu J. & Russell J., 2001. Structural and functional analysis of an mRNP complex that mediates the high stability of human β-globin mRNA. *Mol Cell Biol.* 21(17):5879-88.

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

GENERAL PRODUCT USE

Please note: Due to the lack of a unique restriction site at the 3' end of the fusion gene, this plasmid does not facilitate the subcloning of the Fc-tagged nucleocapsid gene.

- **Stable gene expression in mammalian cells.** pUNO1 plasmids can be used directly in transfection experiments both *in vitro* and *in vivo*. pUNO1 plasmids contain the blasticidin-resistance gene (*bsr*) driven by the CMV promoter/enhancer in tandem with the bacterial EM7 promoter. This allows the amplification of the plasmid in *E. coli*, as well as the selection of stable clones in mammalian cells using the same selective antibiotic. pUNO1 plasmids facilitate high levels of expression and secretion of the gene product.

- **Detection and purification of the nucleocapsid protein.** The pUNO1Fc-SARS2-N plasmid has been designed to generate the nucleocapsid with a C-terminal human IgG1-Fc (Fc) tag to facilitate the detection of the secreted protein with an anti-Fc antibody, and its purification using [protein G](#) affinity resins.

RELATED PRODUCTS

Product	Description	Cat. Code
Blasticidin	Selection antibiotic	ant-bl-1
ChemiComp GT116	Competent <i>E. coli</i>	gt116-11
Protein G / Agarose	Ig binding protein	gel-agg-2
pUNO1-hACE2	Expression vector	puno1-hace2
pUNO1-hTMPRSS2a	Expression vector	puno1-htp2a
pUNO1-hTMPRSS2b	Expression vector	puno1-htp2b
pUNO1His-SARS2-S	Production vector	p1his-cov2-s
pUNO1Fc-SARS2-S	Production vector	p1fc-cov2-s
pUNO1His-SARS2-S1	Production vector	p1his-cov2-s1
pUNO1His-SARS2-RBD	Production vector	p1his-cov2-rbd
pUNO1Fc-SARS2-RBD	Production vector	p1fc-cov2-rbd
pUNO1His-SARS2-N	Production vector	p1his-cov2-n

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