

# HSV-60/LyoVec™

Viral DNA motif complexed with LyoVec™; CDS Ligand

Catalog code: tlr1-hsv60c

<https://www.invivogen.com/hsv60>

For research use only

Version 21L23-MM

## PRODUCT INFORMATION

### Contents

- 100 µg HSV-60/LyoVec™

Note: Each vial contains 25 µg of HSV-60 complexed with 50 µg LyoVec™.

- 10 ml endotoxin-free water

### Sequence

5'-TAAGACACGATGCGATAAAATCTGTTTGTAAATTTATTA-  
3'-ATTCTGTGCTACGCTATTTTAGACAAACATTTTAAATAAT-

-AGGGTACAAATTGCCCTAGC-3'

-TCCCATGTTTAACGGGATCG-5'

### Storage and stability

- Product is shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension, store at 4°C. Resuspended product is stable for 1 week at 4°C when properly stored.

### Quality control

- The biological activity has been verified using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

Intracellular DNA from pathogens is recognized by multiple cytosolic DNA sensors (CDSs), which display contextual preferences for the recognition of DNA<sup>1</sup>. HSV-60 is a double-stranded 60 bp oligonucleotide containing viral DNA motifs<sup>2</sup>. HSV-60 derives from the herpes simplex virus 1 genome. Transfected HSV-60 was shown to potently induce interferon-beta (IFN-β) in a TLR-, DAI and RNA Pol III-independent, but STING-, TBK1- and IRF3-dependent manner. HSV-60 is recognized by the CDSs, DDX41<sup>3</sup> and IFI16<sup>2</sup>. HSV-60 is complexed with the cationic lipid LyoVec™ to facilitate its uptake. CDS ligands, including transfected HSV-60, trigger type I IFN production and the induction of IFN stimulated genes (ISG) through IFN regulatory factors (IRFs). In order to facilitate their study, InvivoGen has developed stable reporter cells in two well established immune cell models, the human monocytic THP-1 cell line and the murine RAW 264.7 macrophages. These cells express a reporter gene, secreted embryonic alkaline phosphatase or luciferase Lucia, under the control of an IRF-inducible promoter. For more information, visit <https://www.invivogen.com/cds-cell-lines>.

1. Sharma S. & Fitzgerald KA. 2011. Innate immune sensing of DNA. PLoS Pathog. 7(4):e1001310. 2. Unterholzner L. et al., 2010. IFI16 is an innate immune sensor for intracellular DNA. Nat Immunol. 11(11):997-1004. 3. Zhang Z. et al., 2011. The helicase DDX41 senses intracellular DNA mediated by the adaptor STING in dendritic cells. Nat Immunol. 12(10):959-65. 4. Arakawa R. et al., 2010. Characterization of LRRFIP1. Biochem Cell Biol. 88(6):899-906. 5. Lippmann J. et al., 2010. IFNbeta responses induced by intracellular bacteria or cytosolic DNA in different human cells do not require ZBP1 (DLM-1/DAI). Cell Microbiol. 10(12):2579-88.

### 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 Asia: +852 3622-3480

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

## METHODS

### Preparation of stock solution (50 µg/ml)

Stimulation of CDS can be achieved with 300 ng-10 µg/ml transfected HSV-60/LyoVec™.

- Add 500 µl of endotoxin-free water (provided) per vial of 25 µg HSV-60/LyoVec™. Mix gently. Allow at least 15 minutes for complete solubilization.
- Store at 4°C. Do not store for more than 1 week.

### Induction of type I IFNs in THP1-Lucia™ ISG cells

Induction of type I IFNs with HSV-60/LyoVec™ can be studied in a variety of cells. The human monocytic cell line THP-1 has been shown to express all the CDSs<sup>2-4</sup>, with the exception of DAI<sup>5</sup>. A protocol for the induction of type I IFNs using THP1-Lucia™ ISG cells, an IRF-Lucia luciferase reporter cell line, is given below:

1. Resuspend HSV-60/LyoVec™, as described above.
2. Stimulate cells with 300 ng/ml-10 µg/ml HSV-60/LyoVec™ complex for 16-48 hours.
3. Monitor induction of type I IFNs by measuring the levels of IRF-induced Lucia luciferase in the cell culture supernatant using QUANTI-Luc™, a Lucia luciferase detection reagent.

## RELATED PRODUCTS

Product	Description	Cat. Code
ISD/LyoVec™	CDS ligand	tlr1-isdc
QUANTI-Luc™	Lucia detection reagent	rep-qlc1
RAW-Lucia™ ISG Cells	Murine macrophages	rawl-isg
THP1-Lucia™ ISG Cells	Human monocytes	thp1-isg
VACV-70/LyoVec™	CDS ligand	tlr1-vav70c