

# 5'ppp-dsRNA Control/LyoVec™

Negative control for 5'ppp-dsRNA complexed with LyoVec™

Catalog code: t1rl-3prnac1v, t1rl-3prnac1v-100

<https://www.invivogen.com/5-ppp-dsrna-lyovec-control>

For research use only

Version 18J26-MM

## PRODUCT INFORMATION

### Contents

- 5'ppp-dsRNA Control/LyoVec™ is available in two quantities:
  - 25 µg 5'ppp-dsRNA Control/LyoVec™: cat. code t1rl-3prnac1v
  - 100 µg (4 x 25µg) 5'ppp-dsRNA Control/LyoVec™: cat. code t1rl-3prnac1v-100
- sterile endotoxin-free water, 1.5 ml with #t1rl-3prnac1v and 10 ml with #t1rl-3prnac1v-100

### Sequence of 5' ppp-dsRNA Control

5'- GCAUGCACCUCUGUUUGA -3' (19 mer)

3'- CGUACGCGGAGACAAACU -5' (19 mer)

The double stranded RNA Control is obtained by hybridization of non-triphosphatase single-stranded 19-mer phosphodiester RNA with its complementary strand (also non-triphosphatase). Each strand is chemically synthesized by solid-phase synthesis and purified by reverse phase HPLC.

### Storage and stability

- 5'ppp-dsRNA Control/LyoVec™ is provided lyophilized and shipped at room temperature. Store lyophilized product at -20°C. Lyophilized product is stable for 12 months when properly stored.
- Upon resuspension, store 5'ppp-dsRNA Control/LyoVec™ at 4°C. Resuspended product is stable for 1 week when properly stored.

### Quality control

- The absence of bacterial contamination, lipoproteins and endotoxins, has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

5' triphosphate double stranded RNA (5' ppp-dsRNA) is a synthetic ligand for retinoic acid-inducible protein I (RIG-I). RIG-I is a cytosolic pattern recognition receptor that senses pathogen-associated molecular patterns (PAMPs) on viral RNA and triggers an antiviral immune response by the activation of type I interferons (IFNs)<sup>1</sup>. RIG-I shares the identification of the dsRNA structure with another sensor MDA-5 (recognition of poly(I:C)/LyoVec™) and is specifically activated by the uncapped 5' triphosphate moiety on viral RNA<sup>2</sup>. A synthetic approach to the exact structure requirement to RIG-I recognition demonstrated that a short blunt double-stranded conformation containing a triphosphate at the 5' end is required<sup>3,4</sup>. 5'ppp-dsRNA Control does not contain the 5' triphosphate moiety, however, caution should be taken as dsRNA without 5' ppp can induce an IFN response in certain cell lines via non-RIG-I pathways, such as TLR7 in plasmacytoid dendritic cells<sup>5</sup>. 5' ppp-dsRNA Control is complexed with the cationic lipid LyoVec™ to facilitate its uptake.

1. Yoneyama M. & Fujita T., 2007. Function of RIG-I-like Receptors in Antiviral Innate Immunity. *J. Biol. Chem.* 282:15315-8. 2. Hornung V. *et al.*, 2006. 5'-Triphosphate RNA is the ligand for RIG-I. *Science.* 314:994-7. 3. Schlee M. *et al.*, 2009. Recognition of 5' triphosphate by RIG-I helicase requires short blunt double-stranded RNA as contained in panhandle of negative-strand virus. *Immunity.* 31(1):25-34. 4. Schmidt A. *et al.*, 2009. 5'-triphosphate RNA requires base-paired structures to activate antiviral signaling via RIG-I. *PNAS* 106(29):12067-72. 5. Schlee M. *et al.*, 2010. The chase for the RIG-I ligand - Recent advances. *Mol Ther.* 18(7):1254-62.

## METHODS

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

- Add 500 µl sterile endotoxin-free water (provided) per vial of 25 µg 5'ppp-dsRNA Control/LyoVec™. Mix gently. Allow at least 15 minutes for complete solubilization.
- Store at 4°C. Do not store for more than 1 week.

Below is a protocol for determining RIG-I stimulation with 5'ppp-dsRNA. Please note that 5'ppp-dsRNA Control has no activity on RIG-I. Use 5'ppp-dsRNA Control/LyoVec™ at the same concentration as 5'ppp-dsRNA/LyoVec™.

### Induction of RIG-I in B16-Blue™ ISG cells

Induction of RIG-I with 5'ppp-dsRNA/LyoVec™ can be studied in a variety of cells expressing RIG-I. If your cell line does not naturally express the RIG-I gene, transfect with a plasmid expressing a RIG-I gene, such as pUNO-hRIG-I or pUNO-mRIG-I.

RIG-I activation can be easily monitored using InvivoGen's B16-Blue™ ISG cells, an IFN regulatory factor (IRF)-inducible secreted embryonic alkaline phosphatase (SEAP) reporter cell line. In these cells, activation of RIG-I triggers the production of IFNs that results in the secretion of SEAP. A protocol for the induction of RIG-I using B16-Blue™ ISG cells is given below:

- Resuspend 5'ppp-dsRNA/LyoVec™, as described above.
- Stimulate cells with 1-10 µg/ml 5'ppp-dsRNA/LyoVec™ for 16-24 hours.

*Note: The use Poly(I:C) (as a positive control) is highly recommended.*

- Monitor induction of RIG-I by measuring the levels of SEAP in the cell culture supernatant using QUANTI-Blue™, a SEAP detection reagent.

## RELATED PRODUCTS

Product	Catalog Code
5'ppp-dsRNA/LyoVec™	t1rl-3prnac1v
B16-Blue™ ISG cells	bb-ifnabg
Poly(I:C) HMW/LyoVec™	t1rl-pic1v
Poly(I:C) LMW/LyoVec™	t1rl-picwlv
pUNO1-hRIG-I (human gene)	puno1-hrighi
pUNO-mRIG-I (mouse gene)	puno-mrighi
QUANTI-Blue™ Solution	req-qbs

## 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)

 **InvivoGen**  
www.invivogen.com