

# Poly(I:C) (LMW)/LyoVec™

RIG-I/MDA-5 Ligand

Catalog code: tlr1-picwlv

<https://www.invivogen.com/polyic-lmw-lyovec>

For research use only

Version 25C25-MM

## PRODUCT INFORMATION

### Contents

- 4 x 25 µg lyophilized poly(I:C) (LMW)/LyoVec™ 1:6 ratio (w/w)  
*Note: Each vial contains 25 µg of poly(I:C) LMW complexed with 150 µg LyoVec™. Poly(I:C) LMW (low molecular weight) is a preparation of poly(I:C) with an average size is 0.2-1 kb.*
- 10 ml endotoxin-free water

### Storage and stability

- Poly(I:C) (LMW)/LyoVec™ is shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension, store product at 4°C. Resuspended product is stable for 1 week at 4°C.

### Quality control

- The biological activity has been tested using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using cellular assays.

## DESCRIPTION

Polyinosinic-polycytidylic acid (poly(I:C)) is a synthetic analog of double-stranded RNA (dsRNA), a molecular pattern associated with viral infection. Poly(I:C) induces a strong innate immune response initiated by two types of pattern recognition receptors (PRRs): the Toll-like receptors (TLRs) and the RIG-I-like receptors (RLRs)<sup>1</sup>. The TLR family consists of more than 10 members expressed on the cell surface membrane or endosomes. The RLRs form a family of cytoplasmic RNA helicases that includes RIG-I and MDA-5. Naked poly(I:C) is recognized by TLR3 whereas transfected poly(I:C) is sensed by RIG-I/MDA-5 in a cell-type-specific manner<sup>2,3</sup>.

Poly(I:C)(LMW)/LyoVec are preformed complexes between poly(I:C) LMW and the transfection reagent LyoVec™. These complexes induce the activation of the RIG-I/MDA-5 signaling pathway at concentrations ranging from 100 ng to 1 µg/ml in InvivoGen's RLR reporter cells.

**1. Kawai T. & Akira S., 2007.** Antiviral signaling through pattern recognition receptors. *J Biochem.* 141(2):137-45. **2. Gitlin L. et al., 2006.** Essential role of mda-5 in type I IFN responses to polyriboinosinic:polyribocytidylic acid and encephalomyocarditis picornavirus. *PNAS* 103(22):8459-8464. **3. Kato H. et al., 2005.** Cell type-specific involvement of RIG-I in antiviral response. *Immunity.* 23(1):19-28.

## METHODS

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

- Add 500 µl endotoxin-free water (provided) and mix gently. Allow at least 15 minutes to resuspend the product.

*Note: The suspension may contain floating fine particles.*

### RIG-I/MDA-5 stimulation

**A549-Dual™ cells** express RIG-I, MDA-5, and an IFN regulatory factor-inducible Lucia luciferase reporter gene which provides a simple method to monitor activation of IFN signaling.

1. Resuspend poly(I:C) (LMW)/LyoVec™ as described above.
2. Add 20 µl of poly(I:C)(LMW)/LyoVec™ at different concentrations (100 ng to 1 µg/ml) per well of a flat-bottom 96-well plate.

#### *Notes:*

- At final concentrations higher than 1 µg/ml, some cytotoxicity may be observed.
  - Naked poly(I:C) LMW may be used as negative control.
3. To each well containing poly(I:C) (LMW)/LyoVec™, add 180 µl of an **A549-Dual™ cell** suspension (50,000 cells per well).
  4. Incubate for 18-24 hours at 37°C.
  5. Determine poly(I:C) (LMW)/LyoVec™ stimulation of RIG-I/MDA-5 by assessing Lucia luciferase reporter gene expression using **QUANTI-Luc™ 4 Lucia/Gaussia**.

## RELATED PRODUCTS

Products	Catalog Code
3p-hpRNA	tlr1-hprna
5'ppp-dsRNA	tlr1-3prna
A549-Dual™ Cells	a549d-nfis
LyoVec™	lyec-1
Poly(I:C) LMW	tlr1-picw
QUANTI-Luc™ 4 Lucia/Gaussia	rep-qlc4lg1

## 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)