Poly(I:C) HMW Biotin High Molecular Weight (HMW)

# Biotin-labeled synthetic analog of dsRNA; TLR3 ligand

Catalog code: tlrl-picb https://www.invivogen.com/polyic-biotin

### For research use only

Version 20114-MM

### PRODUCT INFORMATION Contents

- 10 µg Poly(I:C) HMW Biotin
- 1.5 ml endotoxin-free water

#### Storage and storage

- Poly(I:C) HMW Biotin is shipped at room temperature and can be stored at -20°C for up to 6 months.

- Upon resuspension, prepare aliquots of Poly(I:C) HMW Biotin and store at -20°C. Resuspended product is stable for 3 months at -20°C. Avoid repeated freeze-thaw cycles.

### Quality control:

- TLR3 activity has been verified using HEK-Blue<sup>™</sup> hTLR3 cells.
- The absence of bacterial contamination (e.g. lipoproteins and
- endotoxins) has been confirmed using cellular assays.
- Biotin coupling has been validated by flow cytometry

### DESCRIPTION

Poly(I:C) HMW Biotin is a biotin-labeled Toll-like receptor 3 (TLR3) agonist. It was chemically produced by covalent coupling of biotin with high molecular weight polyinosinic-polycytidylic acid (poly(I:C) HMW). Poly(I:C) HMW, with an average size of 1.5-8 kb, comprises long strands of inosine poly(I) homopolymer annealed to strands of cytidine poly(C) homopolymer.

Poly(I:C) is a synthetic analog of double stranded RNA (dsRNA), a molecular pattern associated with viral infection. Both natural and synthetic dsRNAs are known to induce type I interferon (IFN) production<sup>1</sup>. Depending on its location in the cell, poly(I:C) can activate distinct pattern recognition receptor (PRR) pathways: endosomal poly(I:C) activates TLR3<sup>2</sup>, whereas cytosolic poly(I:C) activates retinoic acid-inducible protein I (RIG-I)<sup>3</sup>, melanoma differentiation-associate gene 5 (MDA-5)<sup>4</sup>, and protein kinase RNAactivated (PKR)<sup>5</sup>. Activation of these PRRs results in the induction of multiple signaling pathways, including NF-κB and IFN regulatory factors (IRFs). Importantly, the labeled ligand retains the biological activity of poly(I:C). Each lot is thoroughly tested to ensure the absence of lot-to-lot variation.

1. Kawasaki T. & Kawai T., 2014. Toll-like receptor signaling pathways. Front Immunol. 5:461. 2. Alexopoulou L. *et al.*, 2001. Recognition of double-stranded RNA and activation of NF-kB by Toll-like receptor 3. Nature, 413:732-8. 3. Kawai T. & Akira S., 2008. Toll-like receptor and RIG-I-like receptor signaling. Ann N Y Acad Sci. 1143:1-20. 4. McCartney S. *et al.*, 2009. Distinct and complementary functions of MDA5 and TLR3 in poly(I:C)-mediated activation of mouse NK cells. J Exp Med. 206(13):2967-76. 5. Lemaire P.A. *et al.*, 2008. Mechanism of PKR Activation by dsRNA. J Mol Biol. 381(2):351-60. 6. Hasan M. *et al.*, 2011. Antimicrobial peptides inhibit polyionsinic-polycytidylic acid-induced immune responses. J Immunol. 187(11):5653-9. 7. Sugimoto N. *et al.*, 2014. Helicase proteins DHX29 and RIG-10 cosense cytosolic nucleic acids in the human airway system. PNAS. 111(21):7747-52.

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 3622-34-80 E-mail: info@invivogen.com

## **APPLICATIONS**

Poly(I:C) HMW Biotin can be used for various applications; including histochemical staining, cytometry, fluorescence microscopy, and ligand binding assays<sup>6.7</sup>.

### **METHODS**

#### Preparation of sterile stock solution (100 $\mu$ g/ml)

 Add 100 μl of the endotoxin-free water provided to the 10 μg Poly(I:C) HMW Biotin vial to obtain a solution at 100 μg/ml.
Homogenize the solution by pipetting up and down until the product is completely dissolved.

#### Working concentration: 30 ng -1 µg/ml

### TLR3 activation with Poly(I:C) HMW Biotin

Poly(I:C) HMW Biotin-induced TLR3 activation can be monitored using TLR3 reporter cell lines, such as HEK-Blue<sup>™</sup> hTLR3 cells. These cells were transfected with the human TLR3 gene and an NF-κB-inducible SEAP (secreted alkaline phosphatase) reporter gene. Levels of SEAP can be easily determined with QUANTI-Blue<sup>™</sup> Soultion, a SEAP detection reagent.

For more information regarding the HEK-Blue hTLR3 cells please visit: <u>https://www.invivogen.com/hek-blue-htlr3</u>.

1. Prepare a HEK-Blue<sup>™</sup> TLR3 cell suspension (280,000 cells/ml) in growth medium without Normocin<sup>™</sup>.

2. In a 96-well plate, add 180 µl of the HEK-Blue<sup>™</sup> TLR3 cell suspension per well.

3. Stimulate cells with 30 ng -1  $\mu\text{g/ml}$  Poly(I:C) HMW Biotin for 6 to 24 h.

4. Determine Poly(I:C) HMW Biotin stimulation of TLR3 by assessing SEAP expression using QUANTI-Blue<sup>™</sup> Solution.

<u>Note:</u> Alternatively, HEK-Blue<sup> $\sim$ </sup> Detection can be used to assess SEAP expression.

# **RELATED PRODUCTS**

Product	Description	Cat. Code
HEK-Blue <sup>™</sup> hTLR3 cells	Human TLR3 reporter cells	hkb-htlr3
HEK-Blue <sup>™</sup> mTLR3 cells	Murine TLR3 reporter cells	hkb-mtlr3
HEK-Blue <sup>™</sup> Detection	SEAP detection reagent	hb-det2
QUANTI-Blue <sup>™</sup> Solution	SEAP detection reagent	rep-qbs
Poly(A:U)	TLR3 ligand	tlrl-pau
Poly(I:C) HMW	TLR3 ligand	tlrl-pic
Poly(I:C) HMW Fluorescein	Fluorescein-labeled poly(I:C)	tlrl-picf
Poly(I:C) HMW Rhodamine	Rhodamine-labeled poly(I:C)	tlrl-picr

