

Poly(I:C) HMW Rhodamine

High Molecular Weight

Rhodamine labeled synthetic analog of dsRNA; TLR3 ligand

Catalog code: tlrl-picr

<https://www.invivogen.com/polyic-rhodamine>

For research use only

Version 21A27-MM

PRODUCT INFORMATION

Contents

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

Storage and stability

- Poly(I:C) HMW Rhodamine is shipped at room temperature. Lyophilized product can be stored at -20°C for up to 6 months. Protect from light.
- Upon resuspension, Poly(I:C) HMW Rhodamine should be aliquoted and stored at -20°C. Protect from light. Resuspended product is stable for 3 months at -20°C. Avoid repeated freeze-thaw cycles.

Quality control

- Human TLR3 (hTLR3) activity tested using HEK-Blue™ hTLR3 cells
- Rhodamine fluorescence evaluated on RAW-Blue™ cells using FACS analysis.

Spectral properties of Rhodamine

Excitation λ max: 546 nm

Emission λ max: 576 nm

DESCRIPTION

Poly(I:C) HMW Rhodamine was chemically labeled by covalent coupling of a rhodamine probe containing a reactive alkylating group. This confers fluorescent properties to poly(I:C) HMW with a slight reduction of TLR3 recognition (see figure 2 on the next page).

Polyinosinic-polycytidylic acid (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 interferons (IFN) and other cytokines production. Poly(I:C) is recognized by Toll-like receptor 3 (TLR3)^{1,2}. Upon poly(I:C) recognition, TLR3 activates the transcription factor interferon regulatory factor 3 (IRF3), through the adapter protein Toll-IL-1 receptor (TIR) domain-containing adapter inducing IFN-β (TRIF, also known as TICAM-1)³. Activation of IRF3 leads to the production of type I IFNs, especially IFN-β. A second pathway involves the recruitment of TNF receptor-associated factor 6 (TRAF6) or receptor interacting protein 1 (RIP1), with the subsequent activation of the transcription factors NF-κB and AP-1⁴. Activation of this pathway triggers the production of inflammatory cytokines and chemokines such as TNF-α, IL-6 and CXCL10. Poly(I:C) is also recognized by the cytosolic RNA helicases retinoic acid-inducible protein 1 (RIG-I) and melanoma differentiation-associate gene 5 (MDA-5)⁵.

1. Alexopoulou L. *et al.*, 2001. Recognition of double-stranded RNA and activation of NF-κB by Toll-like receptor 3. *Nature*, 413:732-8. 2. Matsumoto M. *et al.*, 2002. Establishment of a monoclonal antibody against human Toll-like receptor 3 that blocks double-stranded RNA-mediated signaling. *BBRC* 293:1364-9. 3. Yamamoto M. *et al.*, 2003. Cutting edge: A novel Toll/IL-1 receptor domain-containing adapter that preferentially activates the IFN-β promoter in the Toll-like receptor signaling. *Science* 301:640. 4. Kawai T. & Akira S., 2008. Toll-like receptor and RIG-I-like receptor signaling. *Ann N Y Acad Sci.* 1143:1-20. 5. Kato H. *et al.*, 2006. Small interfering RNAs mediate sequence-independent gene suppression and induce immune activation by signaling through toll-like receptor 3. *Nature* 441:101-5.

APPLICATIONS

Poly(I:C) Rhodamine can be used for various applications:

- flow cytometry
- fluorescent and confocal microscopy.

METHODS

Preparation of sterile stock solution (100 µg/ml)

Stimulation of TLR3 can be achieved with 10 ng-10 µg/ml Poly(I:C) Rhodamine.

1. Add 100 µl of the endotoxin-free water provided to the 10 µg Poly(I:C) HMW Rhodamine vial to obtain a solution at 100 µg/ml.
2. Homogenize the solution by pipetting up and down until complete solubilization. Protect from light.

Fluorescent *in vitro* labeling with Poly(I:C) HMW Rhodamine

The following protocol describes a method to detect Poly(I:C) HMW Rhodamine in murine macrophages RAW-Blue™ cells.

1. Prepare a Raw-Blue™ cell suspension (500,000 cells/ml) in DMEM with 10% (v/v) heat-inactivated fetal bovine serum.
2. In a 96-well plate, add 180 µl of the Raw-Blue™ cell suspension per well.
3. Stimulate cells with 10 ng-10 µg/ml Poly(I:C) HMW Rhodamine for 16 h.
4. Rinse cells with 1X PBS in order to remove free Poly(I:C) Rhodamine.
5. Analyze fluorescent labeling using one of the applications listed. Typical results are shown in figure 1.

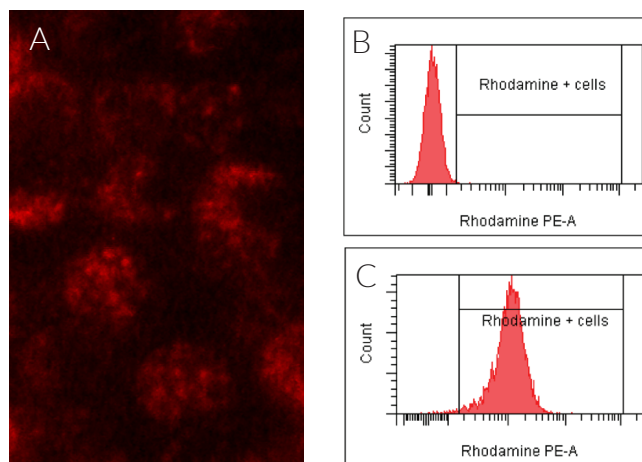


Figure 1. Intracellular fluorescent labeling of RAW-Blue™ (A) cells following a 16 h incubation with 10 µg/ml Poly(I:C) HMW Rhodamine (100,000 cells/well of a 96-well plate). FACS analysis of RAW-Blue™ cells following a 16 h incubation with (B) 10 µg/ml Poly(I:C) HMW and (C) 10 µg/ml Poly(I:C) HMW Rhodamine (100,000 cells/well of a 96-well plate).

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

TLR3 activation with Poly(I:C) HMW Rhodamine

Poly(I:C) HMW Rhodamine-induced TLR3 activation can be monitored using TLR3 reporter cell lines, such as [HEK-Blue™ hTLR3 cells](#). These cells are designed for studying the stimulation of hTLR3 by monitoring the activation of NF-κB. Stimulation with a TLR3 ligand activates NF-κB and AP-1 which induces the production of SEAP. Levels of SEAP can be easily determined with [QUANTI-Blue™ Solution](#) (a detection medium that turns purple/blue in the presence of alkaline phosphatase).

1. Prepare a HEK-Blue™ hTLR3 cell suspension (250,000 cells/ml) in DMEM with 10% (v/v) heat-inactivated fetal bovine serum.
2. In a 96-well plate, add 180 µl of the HEK-Blue™ hTLR3 cell suspension per well.
3. Stimulate cells with 10 ng-10 µg/ml Poly(I:C) HMW Rhodamine for 6 to 24 h.
4. Determine Poly(I:C) HMW Rhodamine stimulation on TLR3 by assessing SEAP reporter expression using a SEAP detection medium, such as [QUANTI-Blue™ Solution](#) or [HEK-Blue™ Detection](#).

A typical stimulation curve is given in figure 2.

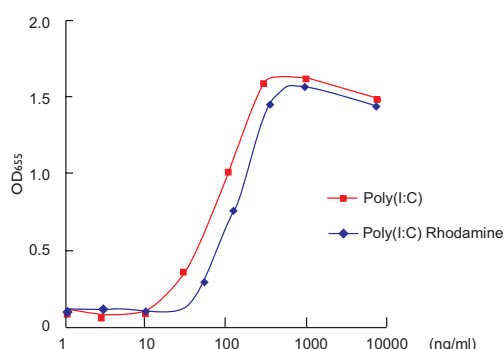


Figure 2. HEK-Blue™ hTLR3 cells were stimulated with increasing concentrations of Poly(I:C) HMW and Poly(I:C) HMW Rhodamine (50,000 cells/well of a 96-well plate). After 18h incubation, NF-κB-induced SEAP activity was assessed using QUANTI-Blue™ Solution.

RELATED PRODUCTS

Product	Description	Cat. Code
RAW-Blue™ cells	Macrophage reporter cells	raw-sp
HEK-Blue™ hTLR3 cells	Human TLR3 reporter cells	hkb-htrl3
HEK-Blue™ Detection	SEAP detection medium	hb-det2
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs
Poly(I:C) HMW	TLR3 ligand	tlrl-pic
Poly(I:C) HMW Fluorescein	Fluorescein labeled Poly(I:C)	tlrl-picf
Poly(I:C) HMW Biotin	Biotin labeled Poly(I:C)	tlrl-picb
Poly(A:U)	TLR3 ligand	tlrl-pau

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