

# Polyinosinic- polycytidylic acid . potassium salt (TLRGRADE®) (synthetic)

TLR3 is a key effector of the immune response against viruses by recognizing double-stranded RNA. Polyinosinic-polycytidylic acid (Poly(I:C)), a synthetic double-stranded homopolymer, binds TLR3 and activates the transcription factor interferon regulator factor 3 (IRF3) following the initiation of TIR domain-containing adaptor protein (TRIF)-dependent TLR signaling. Activation of IRF3 by poly(I:C) leads to the production of type I interferons. Poly(I:C) also activates the RNA helicases MDA-5 and RIG-1. Poly(I:C) can be used to generate stable mature dendritic cells *in vitro* and is considered a potent adjuvant in vaccine formulations, especially those targeting dendritic cells, thanks to the synergy between MDA5 and TLR3 activation.

Citations: 23

[View Online »](#)

## Ordering Information

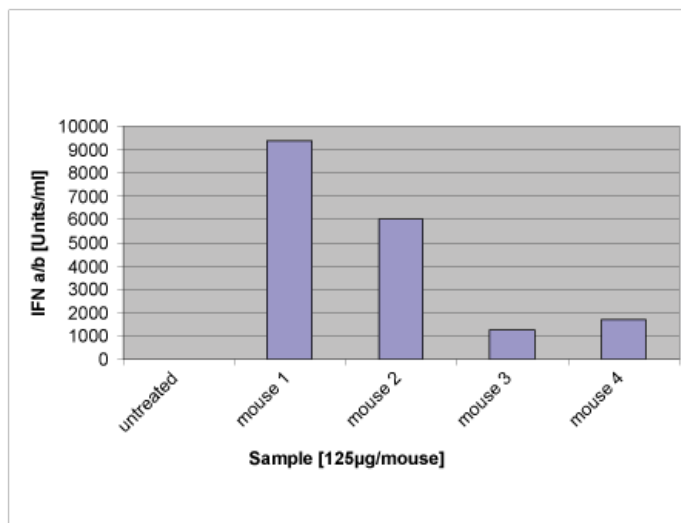
[Order Online »](#)

ALX-746-021-M002	2mg
ALX-746-021-M005	5mg

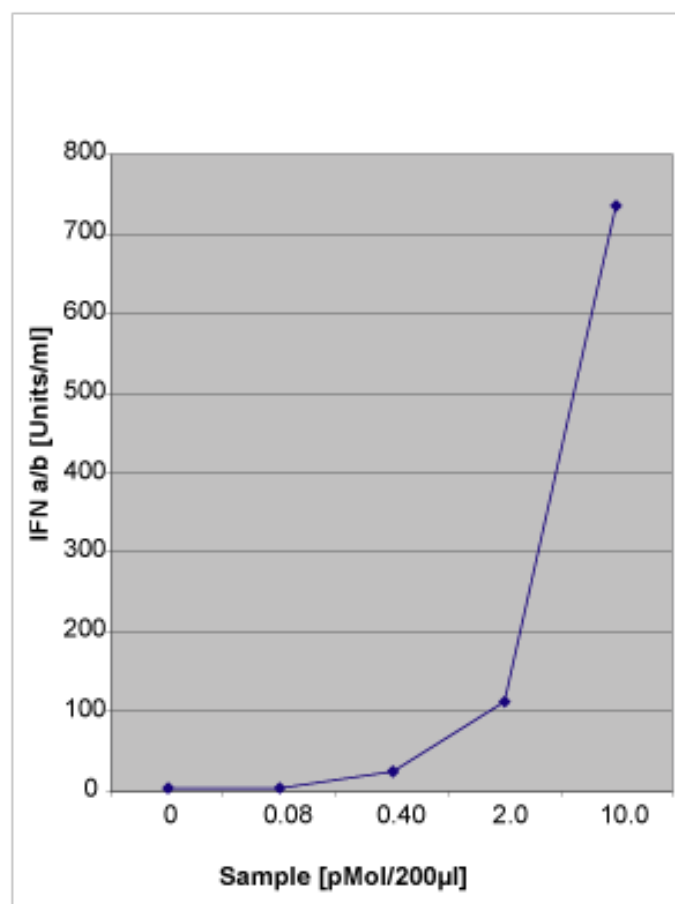
## Manuals, SDS & CofA

[View Online »](#)

- Specific activator of TLR3 and MDA5
- Used to generate stable mature dendritic cells
- Potential adjuvant in cancer vaccines



**Figure 2:** Poly(I:C) . K (TLRGRADE®) (synthetic) (Prod. No. [ALX-746-021](#)) induces cytokine release in mouse *in vivo*. **Method:** Poly(I:C) was injected i.v. and plasma levels of IFN α/β were determined 2 hours later by cytokine ELISA.



**Figure 1:** Poly(I:C) . K (TLRGRADE®) (synthetic) (Prod. No. [ALX-746-021](#)) induces cytokine release in murine BM-derived macrophages. **Method:** Poly(I:C) was added to bone marrow-derived murine macrophages in a 96-well plate at the indicated concentrations, cell supernatants were harvested after 24 hours and IFN α/β levels analyzed by cytokine ELISA.

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Aqueous stock solution is stable for 1 day when stored at +4°C.
Handling	Protect from light. For maximum product recovery after thawing, centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	poly(I:C) . K, poly(I:C)
CAS	31852-29-6
Couple Target	MDA, TLR, TLR3
Couple Type	Activator, Ligand
Endotoxin Content	<0.002EU/μg (LAL test; BioWhittaker)
Formula	$(C_{10}H_{10}N_4NaO_7P)_x \cdot (C_9H_{11}N_3NaO_7P)_x$
Formulation	Lyophilized. Sterile.
Purity Detail	Activity and endotoxin tested – TLRGRADE®.
Quantity	Aliquoted by weight of dry material. Contains polymer, residual salt and water. Sufficient for at least 200 cellular activation assays at 10-25μg/ml.
Reconstitution	For a 1mg/ml stock solution, dissolve the total vial content in 2ml (ALX-746-021-M002) or 5ml (ALX-746-021-M005) endotoxin-free buffer or water. To obtain optimal dissolving we recommend the following procedure: – Add 50% of the solvent and let dissolve for 10 min. – Add remaining 50% of the solvent and mix thoroughly. – Moderate warming may aid dissolving.

Source

Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)