

Validation data for Poly(I:C) HMW

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

For research use only

Version 23J20-MM

Poly(I:C) HMW is a synthetic analog of double-stranded RNA (dsRNA). It is a potent activator of TLR3. Upon agonist recognition, TLR3 interacts with the adaptor molecule TRIF and triggers downstream signaling pathways that lead to the activation of interferon regulatory factor (IRF3), NF- κ B and AP-1. The biological activity of Poly(I:C) HMW has been tested using InvivoGen's HEK-Blue™ hTLR3 cells which stably express human TLR3 and an NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter (Figure 1).

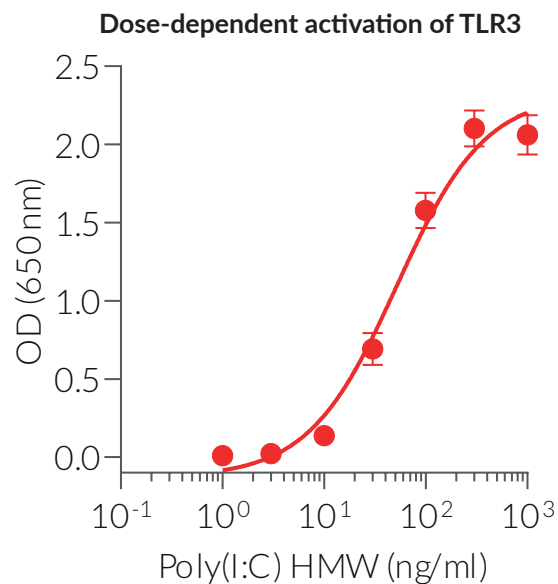


Figure 1. Poly(I:C) HMW is a potent activator of human (h)TLR3. HEK-Blue™ hTLR3 cells were incubated in HEK-Blue™ Detection medium and stimulated with increasing concentrations of Poly(I:C) HMW. After 24h incubation, the levels of NF- κ B-induced SEAP were determined by reading the optical density (OD) at 630 nm (mean \pm SEM).

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