

Validation data for ssPolyU Naked

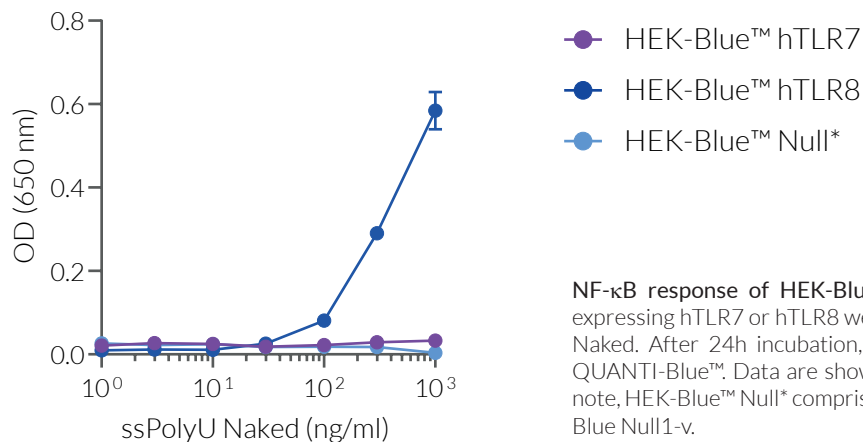
<https://www.invivogen.com/sspolyu>

For research use only

Version 24A11-AK

ssPolyU Naked is a lyophilized and non-complexed preparation of single-stranded poly-uridine (ssPolyU). It is not complexed with the cationic lipid LyoVec™; thus, it is not protected from degradation. The biological activity of ssPolyU Naked has been tested using InvivoGen's HEK-Blue™ cells expressing human TLR7 or TLR8 together with an NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter (Figure 1).

Dose-dependent NF- κ B response in HEK-Blue™-derived cells



NF- κ B response of HEK-Blue™-derived cells to ssPolyU Naked. HEK-Blue™ cells expressing hTLR7 or hTLR8 were stimulated with increasing concentrations of ssPolyU Naked. After 24h incubation, the NF- κ B-induced SEAP activity was assessed using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean \pm SEM). Of note, HEK-Blue™ Null* comprises data from parental cell lines HEK-Blue Null1 and HEK-Blue Null1-v.

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