

Validation data for ssRNA40/LyoVec™

<https://www.invivogen.com/ssrna40-lv>

For research use only

Version 24A10-AK

ssRNA40/LyoVec™ is a 20-mer single-stranded RNA oligonucleotide containing a GU-rich sequence. It is complexed with the cationic lipid LyoVec™ to protect it from degradation and facilitate its uptake. Moreover, phosphorothioate linkages were incorporated in order to extend the effective molecular lifetime by minimizing extra and intracellular nuclease degradation. The biological activity of ssRNA40 has been tested using InvivoGen's HEK-Blue™ cells expressing human or murine 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

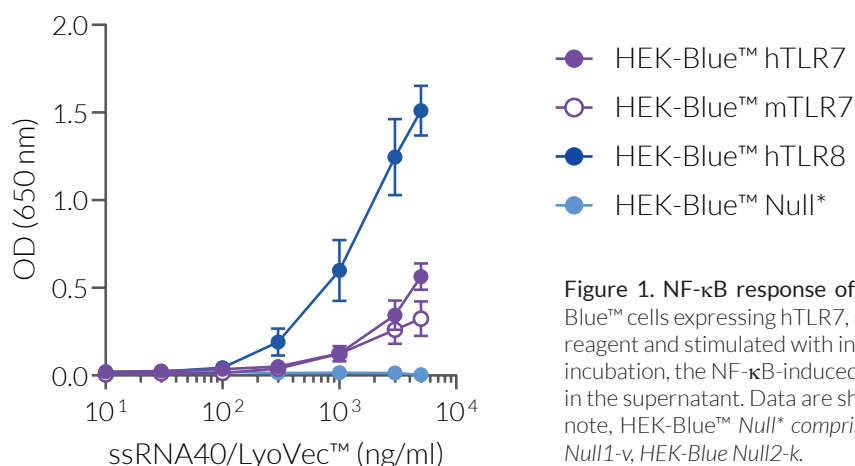


Figure 1. NF- κ B response of HEK-Blue™-derived cells to ssRNA40/LyoVec™. HEK-Blue™ cells expressing hTLR7, mTLR7, or hTLR8 were cultured in HEK-Blue™ Detection reagent and stimulated with increasing concentrations of ssRNA40/LyoVec™. After 24h incubation, the NF- κ B-induced SEAP activity was assessed by measuring the SEAP level in the supernatant. 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, HEK-Blue Null1-v, HEK-Blue Null2-k.

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