

# Validation data for ssRNA41/LyoVec™

<https://www.invivogen.com/ssrna41-lv>

For research use only

Version 24A10-AK

ssRNA41/LyoVec™ is a 20-mer single-stranded RNA oligonucleotide. It derives from ssRNA40 by replacement of all U nucleotides with adenosine and is used as a negative control for ssRNA40. 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. In comparison to ssRNA40, ssRNA41/LyoVec™ is unable to activate InvivoGen's HEK-Blue™ hTLR8 cells (**Figure 1**).

NF-κB response of HEK-Blue™ hTLR8 cells

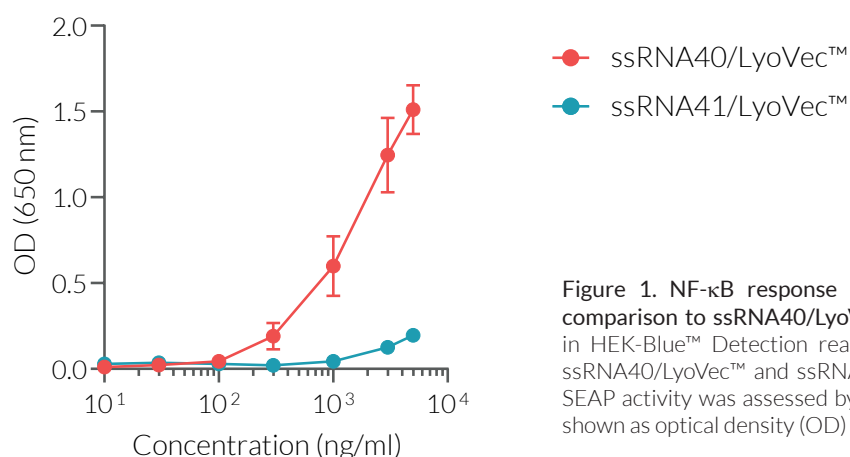


Figure 1. NF-κB response of HEK-Blue™ hTLR8 cells to ssRNA41/LyoVec™ in comparison to ssRNA40/LyoVec™. HEK-Blue™ cells expressing hTLR8 were cultured in HEK-Blue™ Detection reagent and stimulated with increasing concentrations of ssRNA40/LyoVec™ and ssRNA41/LyoVec™. After 24h incubation, the NF-κB-induced SEAP activity was assessed by measuring the SEAP level in the supernatant. Data are shown as optical density (OD) at 650 nm (mean ± 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](mailto:info@invivogen.com)