

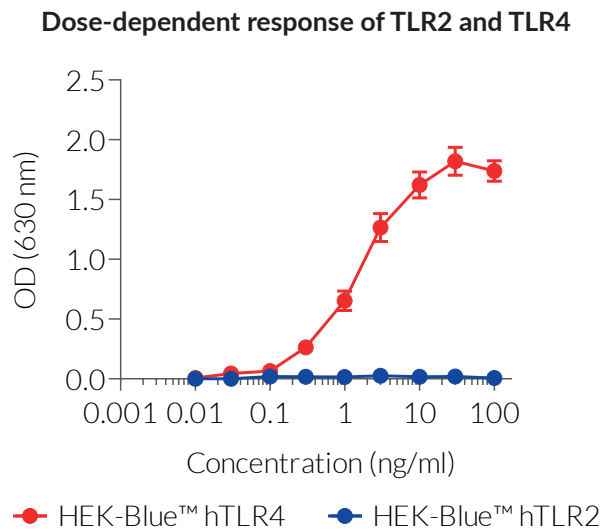
# Validation data for LPS-EK Ultrapure

<https://www.invivogen.com/lps-ek>

For research use only

Version 23F14-AK

LPS-EK is a preparation of a rough (r)-form of lipopolysaccharide (LPS) purified from the Gram-negative *E. coli* K12, a prototypical laboratory strain. It is the preferred model in biochemical genetics, molecular biology, and biotechnology. LPS-EK Ultrapure (UP) is extracted by successive enzymatic hydrolysis steps and purified by the previously described phenol-TEA-DOC extraction protocol. This process removes contaminating lipoproteins. Therefore LPS-EK UP only activates TLR4, as verified using InvivoGen's HEK-Blue™ hTLR2 and HEK-Blue™ hTLR4 cells (**Figure 1**).



**Figure 1. LPS-EK UP is a potent activator of human (h)TLR4.** The cells were incubated with increasing concentrations of LPS-EK UP. After overnight incubation in HEK-Blue™ detection medium, a SEAP detection growth medium, the response of hTLR2 and hTLR4 was assessed by determining the presence of SEAP in the supernatant. Data are expressed as optical density at 630 nm ( $\pm$ 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](mailto:info@invivogen.com)