

Validation data for LPS-EB (Standard)

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

For research use only

Version 23F14-AK

LPS-EB is a preparation of a smooth (s)-form lipopolysaccharide (LPS) purified from the Gram-negative *E. coli* O111:B4, a pathogenic serotype of *E. coli* known to cause significant gastric disease. The standard LPS-EB preparation is extracted by a phenol-water mixture. It contains other bacterial components, such as lipoproteins, and therefore stimulates both TLR4 and TLR2, as verified using InvivoGen's HEK-Blue™ hTLR2 and HEK-Blue™ hTLR4 cells (Figure 1).

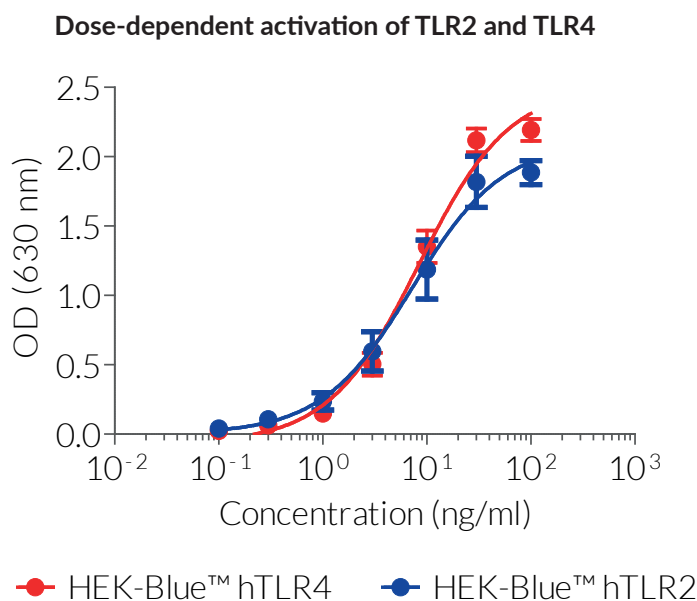


Figure 1. LPS-EB (Standard) is a potent activator of human (h)TLR2 and hTLR4. The cells were incubated with increasing concentrations of LPS-EB (Standard). After overnight incubation in HEK-Blue™ detection medium, a SEAP detection growth medium, the activation of hTLR2 and hTLR4 was assessed by determining 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