

Validation data for LPS-RS (Standard)

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

For research use only

Version 23F14-AK

Lipopolysaccharide (LPS)-RS from the photosynthetic bacterium *Rhodobacter sphaeroides* is a potent antagonist of LPS from pathogenic bacteria. It does not induce TLR4 signaling but is detected by the LAL assay, the standard endotoxin detection assay. The standard LPS-RS preparation is extracted by a phenol-water mixture. It contains other bacterial components, such as lipoproteins, and therefore stimulates TLR2 (Figure 1) while inhibiting LPS-induced TLR4 activation (Figure 2), as verified using InvivoGen's HEK-Blue™ hTLR2 and HEK-Blue™ hTLR4 cells.

Dose-dependent activation of TLR2 and TLR4

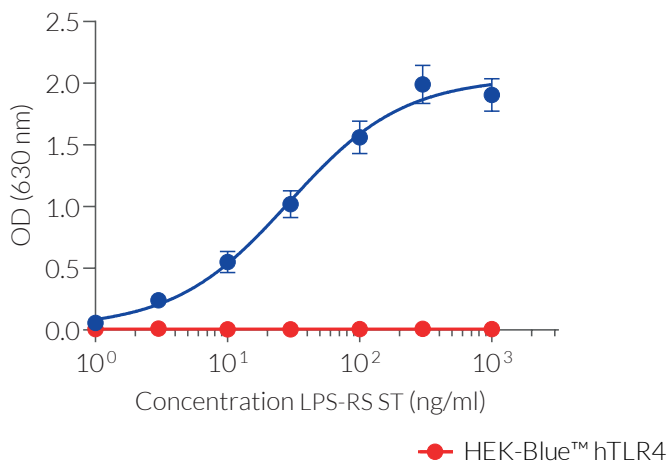


Figure 1. LPS-RS Standard (ST) is a potent activator of human (h)TLR2. The cells were incubated with increasing concentrations of LPS-RS ST. After overnight incubation in HEK-Blue™ detection medium, a SEAP detection growth medium, the response of hTLR2 and hTLR4 was assessed by determining the presence of SEAP in the supernatant. Data are expressed as optical density at 630 nm (\pm SEM).

Dose-dependent inhibition of LPS-induced TLR4 activation

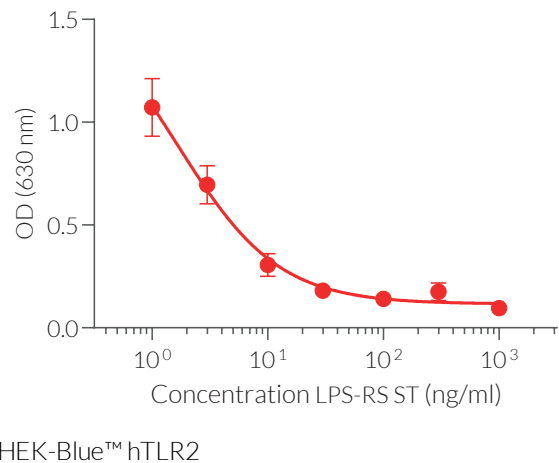


Figure 2. LPS-RS ST is a potent antagonist of LPS-EK Ultrapure (UP). The cells were treated with 0.3 ng/ml LPS-EK UP and increasing concentrations of LPS-RS ST. After overnight incubation, the inhibition of LPS-induced hTLR4 activation was assessed by determining the presence of SEAP in the supernatant using QUANTI-Blue™ Solution. 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