

Validation data for MPLA-SM*

<https://www.invivogen.com/mpla>

For research use only

Version 22K24-NJ

Monophosphoryl Lipid A (MPLA-SM) is a TLR4 agonist derived from Lipid A, the immunostimulatory structure of lipopolysaccharide (LPS). This natural compound is extracted from the LPS of *Salmonella minnesota* Re595 (Re mutant). The preparation is a mix of MPLA congeneric forms differing in the number of acyl chains, and possibly responsible for the partial TLR4 agonist function of some preparations. MPLA-SM* results from an improved process of MPLA-SM extraction. While MPLA-SM* and MPLA-SM have the same ability to activate murine TLR4 (Figure 1), MPLA-SM* is more potent than MPLA-SM at inducing human TLR4 responses (Figure 2).

Mouse TLR4 activation by *S. minnesota* monophosphoryl lipid A

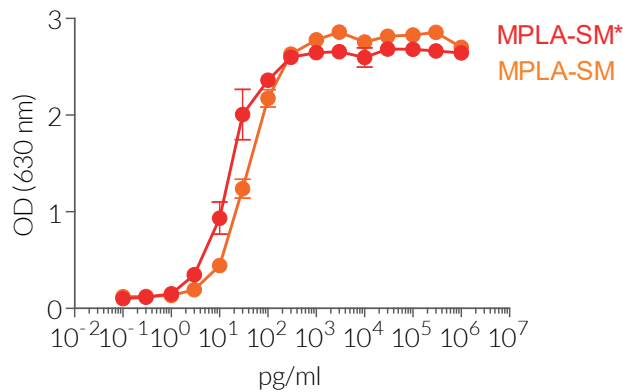


Figure 1. MPLA-SM and MPLA-SM* induce a similar dose-dependent response in HEK-Blue™ mTLR4 cells.

The cells were incubated with increasing concentrations of two preparations of *S. minnesota* monophosphoryl lipid A, MPLA-SM and MPLA-SM*. After overnight incubation in HEK-Blue™ detection medium, a SEAP detection growth medium, the activation of mouse (m)TLR4 was assessed by determining the presence of SEAP in the supernatant. Data are expressed as optical density at 630 nm (±SEM).

Human TLR4 activation by *S. minnesota* monophosphoryl lipid A

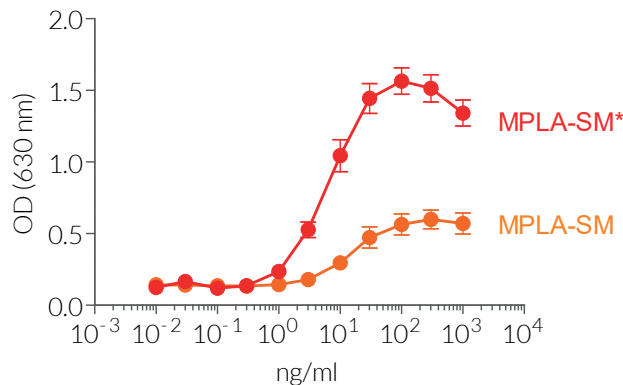


Figure 2. MPLA-SM* is more potent than MPLA-SM at inducing a dose-dependent response in HEK-Blue™ hTLR4 cells.

The cells were incubated with increasing concentrations of two preparations of *S. minnesota* monophosphoryl lipid A, MPLA-SM and MPLA-SM*. After overnight incubation in HEK-Blue™ detection medium, a SEAP detection growth medium, the activation of human (h)TLR4 was assessed by determining the presence of SEAP in the supernatant. Data are expressed as optical density at 630 nm (±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