

Validation data for CRX-527

<https://www.invivogen.com/crx527>

For research use only

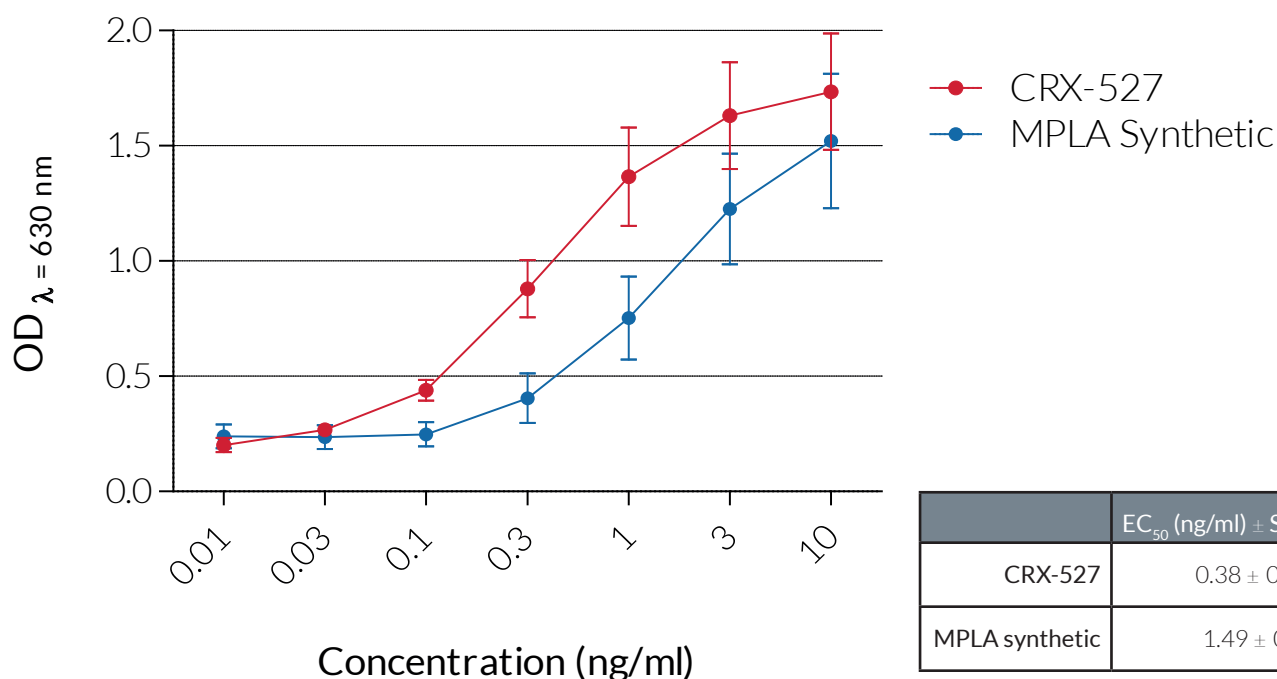
Version 19C18-ED

CRX-527 is a synthetic lipid A mimic belonging to the family of aminoacyl glucoaminide 4-phosphates (AGPs). Lipid A is the immunostimulatory structure of the bacterial-derived Toll-like receptor 4 (TLR4) agonist, lipopolysaccharide (LPS). CRX-527 is a highly-specific TLR4 agonist and is described as the most powerful synthetic lipid A mimic of the AGP family^{1,2}. Stimulation of InvivoGen's HEK-Blue™ human (h)TLR4 cells with CRX-527 results in a clear dose-dependent activation of TLR4, which is comparable to another LPS-like molecule, synthetic monophosphoryl lipid A (MPLA).

1. Stover, A.G. *et al.*, 2003. Structure-activity relationship of synthetic Toll-like receptor 4 agonists. *J. Biol. Chem.* 279:4440-4449.

2. Legat, A. *et al.*, 2010. CD14-independent responses induced by a synthetic lipid A mimetic. *Eur. J. Immunol.* 40:792-802.

Evaluation of TLR4 activation by CRX-527



CRX-527 induces a dose-dependent response in HEK-Blue™ hTLR4 cells. The cells were incubated with increasing concentrations of CRX-527 and synthetic monophosphoryl lipid A (MPLA). After overnight incubation in HEK-Blue™ detection medium, a SEAP detection growth medium, the activation of TLR4 was assessed by determining the presence of SEAP in the supernatant. Data are expressed as OD (λ=630nm) (±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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com