

Validation data for TL7-887 VacciGrade™

<https://www.invivogen.com/tlr7-conjugatable-ligands>

For research use only

Version 22D21-NJ

TL7-887 is a ready-to-use “pre-linked” conjugatable TLR7 ligand, synthesized from the base molecule CL307, a well-known TLR7 agonist. TL7-887 efficiently triggers a cellular response upon recognition by either human or murine TLR7 (Figure 1). TL7-887 can be used to generate immune-stimulating antibody conjugates (ISACs) as conjugation to a Anti-HER2-hlgG1 and subsequent activation of TLR7 has been validated using cellular assays. The Anti-HER2/TL7-887 ISAC is more potent at inducing a TLR7-mediated response than TL7-887 only in cells expressing HER2 (Figure 2A). Of note, at high concentrations, TL7-887, Anti-HER2/TL7-887 ISAC and Anti-βGal/TL7-887 ISAC, induce a TLR7-mediated cellular response, independently of HER2 expression (Figure 2A and B). This observation could be explained by cellular uptake through endocytosis/pinocytosis.

Biological activity of TL7-887

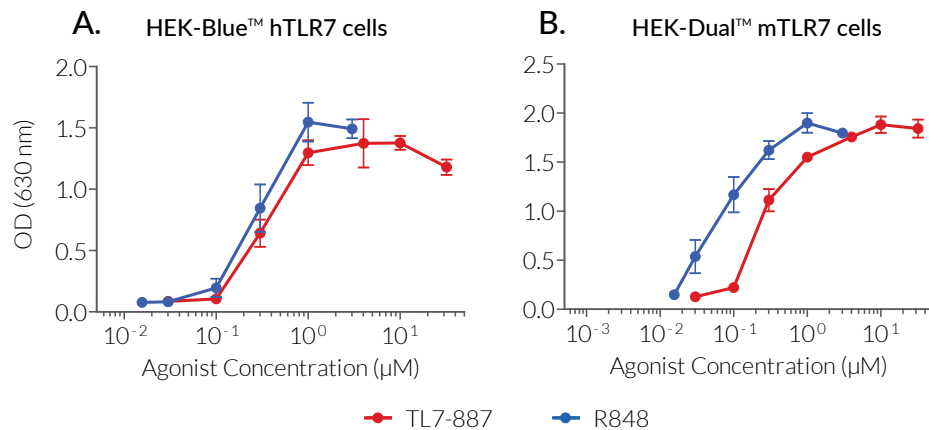


Figure 1: Dose-response of human and murine TLR7 reporter cells to TLR7 conjugatable ligand TL7-887.

~ 4x10⁵ HEK-Blue™ hTLR7 (A) or HEK-Dual™ mTLR7 (B) cells were stimulated with increasing concentrations of TL7-887 (TLR7 agonist), or R848 (TLR7/8 agonist) as a control. Cells were incubated overnight in HEK-Blue™ Detection, a cell culture medium that allows real-time detection of SEAP activity in the supernatant. The optical density (OD) at 630 nm is shown as mean ± SEM.

Biological activity of TL7-887 conjugated to Anti-HER2-hlgG1

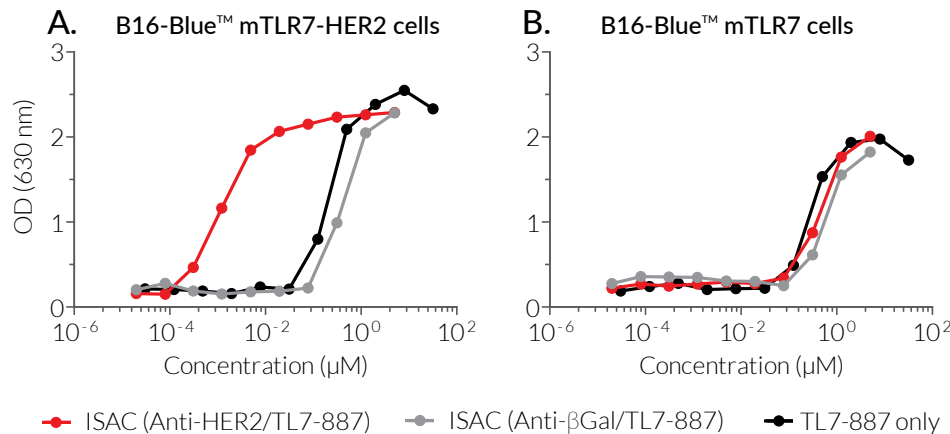


Figure 2: Dose-response of HER2-expressing and murine TLR7 reporter cells to Anti-HER2/TL7-887 ISAC.

~ 5x10⁵ B16-Blue™ mTLR7-HER2 cells (A) or B16-Blue™ mTLR7 control cells (B) were stimulated with increasing concentrations of Anti-HER2/TL7-887 ISAC (DAR ~8), Anti-βGal/TL7-887 ISAC (DAR ~8), or TL7-887 only. After overnight incubation, the TLR7 response was determined using QUANTI-Blue™ Solution, a SEAP detection reagent. The optical density (OD) at 630 nm is shown.

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