

Validation data for TL7-975

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

For research use only

Version 23J12-NJ

TL7-975 is a “click-chemistry compatible” conjugatable TLR7 ligand featuring an azido group added to the N-terminus of the base molecule CL307, a well-known TLR7 agonist. TL7-975 efficiently triggers a cellular response upon recognition by either human or murine TLR7 (**Figure 1**). TL7-975 can be used to generate immunostimulatory antibody-drug conjugates (ADCs) as conjugation to a Anti-HER2-hIgG1 and subsequent activation of TLR7 has been validated using cellular assays (**Figure 2**). The Anti-HER2/TL7-975 ADC is more potent at inducing a TLR7-mediated response than TL7-975 only in cells expressing HER2 (**Figure 2A**). Of note, at high concentrations, TL7-975, Anti-HER2/TL7-975 ADC and Anti- β -Gal/TL7-975 ADC, 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 TL7-975

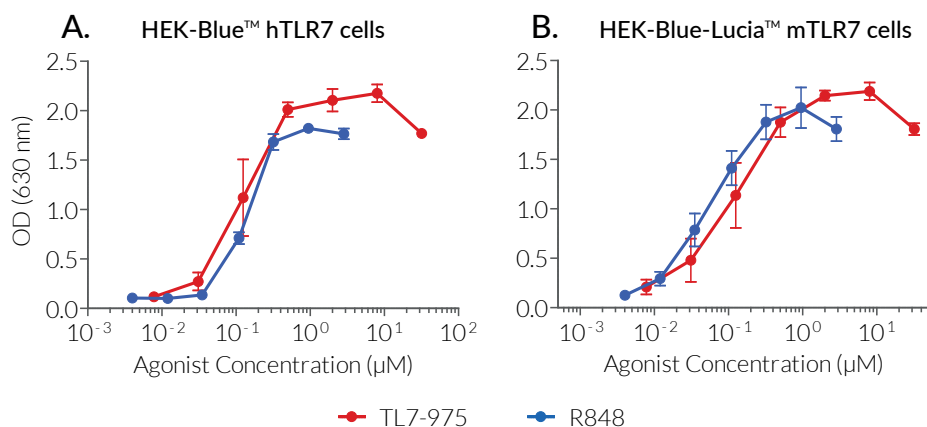


Figure 1: Dose-response of human and murine TLR7 reporter cells to TL7-975.

$\sim 4 \times 10^5$ HEK-Blue™ hTLR7 (**A**) or HEK-Blue-Lucia™ mTLR7 (**B**) cells were stimulated with increasing concentrations of TL7-975 (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 \pm SEM.

Biological activity of TL7-975 conjugated to Anti-HER2-hIgG1

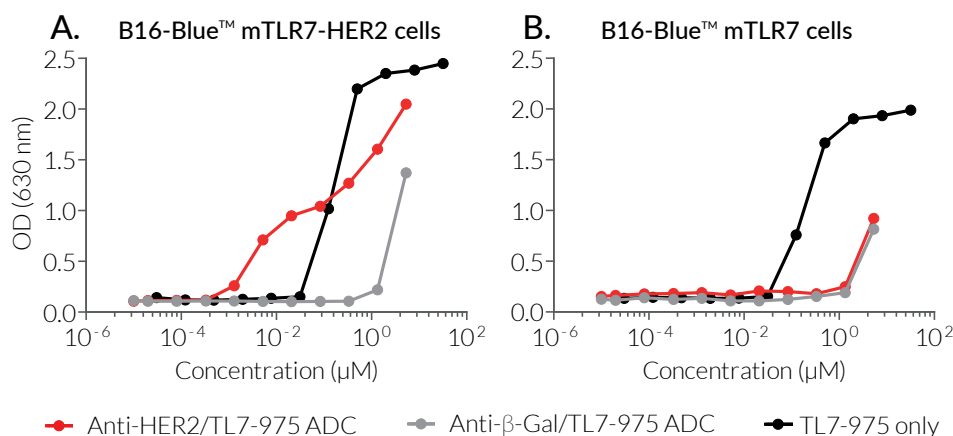


Figure 2: Dose-response of HER2-expressing and murine TLR7 reporter cells to Anti-HER2/TL7-887 ADC.

$\sim 5 \times 10^5$ B16-Blue™ mTLR7-HER2 cells (**A**) or B16-Blue™ mTLR7 control cells (**B**) were stimulated with increasing concentrations of Anti-HER2/TL7-975 ADC (Ratio 1:12), Anti- β -Gal/TL7-975 ADC (Ratio 1:12), or TL7-975 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