TL7-887 Conjugatable TLR7 ligand (CL307) - VacciGrade™ Catalog Code: vac-tl7887, vac-tl7887-1 https://www.invivogen.com/tlr7-conjugatable-ligands

For research use only Version 24J24-NJ

PRODUCT INFORMATION

Contents

TL7-887 is available in two pack sizes:

- 250 μg: vac-tl7887
- 1 mg (4 x 250 μg): vac-tl7887-1

Note: TL7-887 appears as a lyophilized pellet or a light-yellow droplet.

Storage and stability

TL7-887 is shipped at room temperature. Upon receipt, store at -20 °C. Lyophilized product is stable for 6 months when properly stored.
Upon resuspension, prepare aliquots and store at -20 °C. Resuspended product is stable for at least 6 months when properly stored. Avoid repeated freeze-thaw cycles.

Quality control

• Purity: ≥ 95% (UHPLC)

• TL7-887 is VacciGrade[™] (preclinical grade). It is prepared under strict aseptic conditions and is guaranteed sterile. Absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue[™] hTLR2 and HEK-Blue[™] hTLR4 cells, and its endotoxin level is ≤5 EU/mg (measurement by kinetic chromogenic LAL assay).

• Biological activity has been confirmed using cellular assays.

• Conjugation to an Anti-HER2-hlgG1 and Anti-TROP2-hlgG1 mAbs and subsequent activation of TLR7 has been validated using cellular assays.

PRODUCT DESCRIPTION

TL7-887 is a conjugatable TLR7 ligand, synthesized from the base molecule CL307, a well-known specific TLR7 agonist. TL7-887 features a maleimide functional group that has been added to CL307 via a spacer. The bioconjugation to a protein of interest (POI) is driven by the maleimide reaction with free thiols on the POI, which generates a covalent thioether bond.

TL7-887 is a <u>ready-to-use</u>, "<u>pre-linked</u>" reagent, provided with one example of thioether conjugation protocol (see next page).

APPLICATIONS

TL7-887 conjugatable TLR7 ligand has been designed to allow the generation of a bioconjugate after attachment to a biomolecule with a chemical linker. Examples of therapeutic bioconjugates include immunostimulatory antibody-drug conjugates (ADCs) and antigenadjuvant conjugates (AACs).

• ADCs allow localized TLR activation and antibody-mediated effector functions ^{1,2}.

• AAC vaccines allow the codelivery of antigen and TLR agonist to antigen presenting cells (APCs) and thus, better antigen processing and presentation for the induction of adaptative immune responses³.



Formula: $C_{45}H_{69}N_{13}O_7$ Molecular weight: 904 g/mol Mol. ext. coefficient ε_{280nm} : 1140 M⁻¹ cm⁻¹ Solubility: 10 mg/ml (11 mM) in H₂O



1. Gadd A.J, et al. 2015. Targeted activation of toll-like receptors: conjugation of a toll-like receptor 7 agonist to a monoclonal antibody maintains antigen binding and specificity. Bioconjug. Chem. 26(8):1743. 2. Ackerman S.E, et al. 2021. Immune-stimulating antibody conjugates elicit robust myeloid activation and durable antitumor immunity. Nat Cancer. 2(1):18. 3. Xu Z. & Moyle P.M, 2018. Bioconjugation approaches to producing subunit vaccines composed of protein or peptide antigens and covalently attached Toll-Like receptor ligands. Bioconjug. Chem. 29(3):572.

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



Below is a protocol for cysteine-based thioether coupling of TL7-887 to a protein of interest (POI).

Samples are harvested throughout the procedure for calculation of the ligand to POI ratio (DAR) using a spectrophotometer.

<u>Note:</u> for more information on conjugation methods and DAR calculation, visitourwebsite at <u>https://www.invivogen.com/faq-conjugatable-ligands</u>. <u>Note:</u> lysine-based thioether coupling protocols are available in the literature².

CYSTEINE-BASED THIOETHER COUPLING

This strategy is widely used to generate monoclonal antibody (mAb) conjugates. It aims at targeting the cysteine residues forming the four inter-chain disulfides of the IgG.

Materials required (not supplied):

- Conjugation Buffer (pH 7.5): 0.1 M Sodium Phosphate, 0.1 M NaCl, 1 mM EDTA

- Reducing Agent: 0.2 M TCEP (tris(2-carboxyethyl) phosphine), freshly prepared

- **Stopping Buffer** (pH 8.0): 0.1 M Disodium Phosphate (Na₂HPO₄), 0.1 M NaCl, 50 mM Sodium Tetraborax (Na₂B₄O₇)
- Final Buffer (pH 6.5-7.5): 20 mM Sodium Phosphate, 0.15 M NaCl
- **OD Buffer**: Final Buffer + 0.2 M Cysteine
- Desalting spin columns 0.5 ml (refer to manufacturer's instructions)
- PVDF Syringe filters 0.22 μm, 4 mm or13 mm

Preparation of the mAb of interest

1. Prepare a solution of antibody in **Conjugation Buffer** at ~5 mg/ml. <u>IMPORTANT: the mAb solution must contain no tris</u>, cysteine, glycine, nor any other amine or thiol function.

2. <u>Sample 1:</u> Harvest and dilute 1 μ l of mAb solution into 9 μ l of **OD** Buffer. Store at 4°C until required.

3. Add Reducing Agent (1 mM final) to the mAb solution.

4. Vortex briefly and incubate for 90 to 120 min at 37°C using a plate or rotation shaker.

Preparation of TL7-887 (6.25 mg/ml; 6.9 mM)

Note: Ensure you see the lyophilized pellet or light-yellow droplet before resuspension.

- 1. Add 40 μL sterile water to 250 μg TL7-887 vial.
- 2. Mix gently by pipetting until the product is completely dissolved.
- 3. Do a quick spin down to collect all the material from the tube wall.
- 4. <u>Sample 2</u>: Harvest and dilute 1 μ l of TL7-887 solution into 49 μ l of **OD Buffer**. Store at 4°C until required.

Conjugation

Below is a protocol to obtain a TL7-887-ADC with a DAR ~8 using a TL7-887 to Anti-HER2-hlgG1 molar ratio of 12:1.

<u>Note:</u> this example has been used to validate the bioconjugate's potency at stimulating TLR7 and targeting HER2 using in-house cellular assays.

<u>IMPORTANT</u>: the user should assess different molar ratios to optimize the DAR depending on the immunoglobulin of interest and the following applications (in vitro or in vivo).

- 1. Add 12 μI of TL7-887 to 200 μI of TCEP-treated mAb solution.
- 2. Vortex briefly and incubate for 4 hours at 4°C using a plate or rotation shaker.
- 3. Desalt the conjugate against $\ensuremath{\textbf{Stopping Buffer}}$ using a spin column.
- 4. Incubate at room temperature for ~30 min.
- 5. Desalt the conjugate against $\ensuremath{\mathsf{Final}}\xspace$ Buffer using a spin column.
- 6. <u>Sample 3</u>: Harvest and dilute 2 μl of the ADC and dilute into 8 μl of the **OD Buffer**. Store at 4°C until required.
- 7. Filter sterilize the ADC using a 0.22 μ m PVDF syringe filter.
- 8. Store ADC at 4°C (short-term) or -20°C (long-term).

Below is a summary table of calculated volumes of TL7-887 depending on desired DAR with a standard 150 kDa mAb.

| Theoretical DAR | TL7-887/mAb Molar Ratio | Volume of TL7-887 per mg of mAb |
|-----------------|----------------------------|------------------------------------|
| ~ 8 | 12 | 12 μL |
| ~ 5 | 9 | 9 μL |
| ~ 4 | 6 | 6 µL |
| ~ 2 | 4 | 4 μl |

RELATED PRODUCTS

| Product | Description | Cat. Code |
|---|---|--|
| Anti-HER2-hlgG1 Anti-TROP2-hlgG1 Anti-β-Gal-hlgG1* CL307 | Monoclonal antibody Monoclonal antibody Monoclonal antibody TLR7 agonist tlrl-c307 | her2-mab1-1 trop2-mab1-1 bgal-mab1-1 |
| HEK-Blue™ hTLR7 cells HEK-Blue™ Detection | Human TLR7 reporter cells Detection reagent | hkb-htlr7 hb-dt2 |

