

Validation data for CU-CPT9a

<https://www.invivogen.com/cucpt9a>

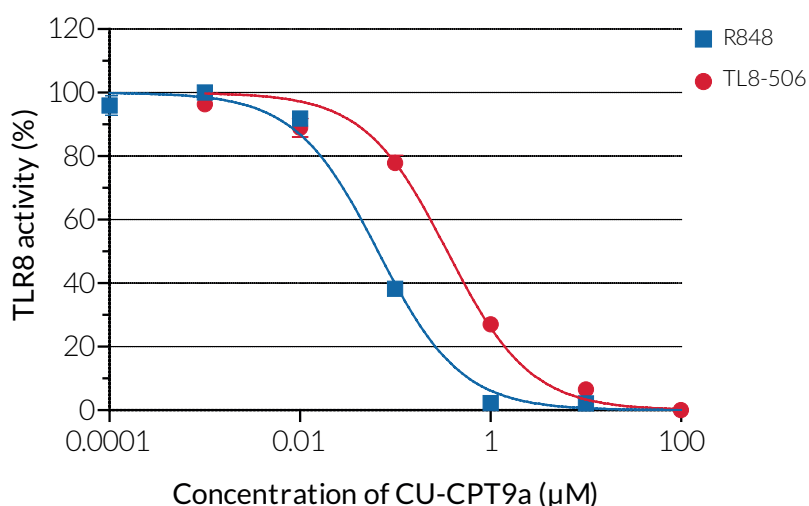
For research use only

Version 19J25-ED

CU-CPT9a is a potent and selective inhibitor of human Toll-like receptor 8 (hTLR8) (Figure 1 and 2). CU-CPT9a stabilizes the TLR8 dimer in its resting state and not only prevents TLR8 activation but also antagonizes any binding of TLR8 ligands.

Evaluation of hTLR8 inhibition by CU-CPT9a

Treatment of HEK-Blue™ hTLR8 cells with CU-CPT9a results in the inhibition of the inducible NF-κB response in a dose-dependent manner upon incubation with either R848, a TLR7/TLR8 agonist or TL8-506, a specific TLR8 agonist. Please note that at high concentrations (i.e. 100 mM) crystal-like structures may be observed in the resuspended CU-CPT9a product.



CU-CPT9a IC ₅₀ (μM) ± Std Error	
R848	0.066 ± 0.004
TL8-506	0.352 ± 0.09

Figure 1: CU-CPT9a inhibits hTLR8 in a dose-dependent response in HEK-Blue™ hTLR8 cells. The cells were incubated with increasing concentrations of CU-CPT9a for 3 hours. Following this, either TL8-506 (10 μg/ml), a specific TLR8 agonist, or R848 (10 μg/ml), a TLR7/TLR8 agonist, were added to the cells. After overnight incubation, activation of TLR8 (NF-κB activity) was assessed by measuring SEAP activity in the supernatant, using QUANTI-Blue™ Solution, a SEAP detection reagent. Data are shown as a percentage (%) of TLR8 activity (no inhibitor).

Specificity of CU-CPT9a

Treatment of HEK-Blue™ hTLR8, hTLR7, and murine (m)TLR7 cells with CU-CPT9a results in the inhibition of the inducible NF-κB response in the hTLR8 cells only upon incubation with R848, a TLR7/TLR8 agonist. There was minimal to no inhibition noted for both human and murine TLR7, highlighting the highly specific nature of CU-CPT9a for TLR8.

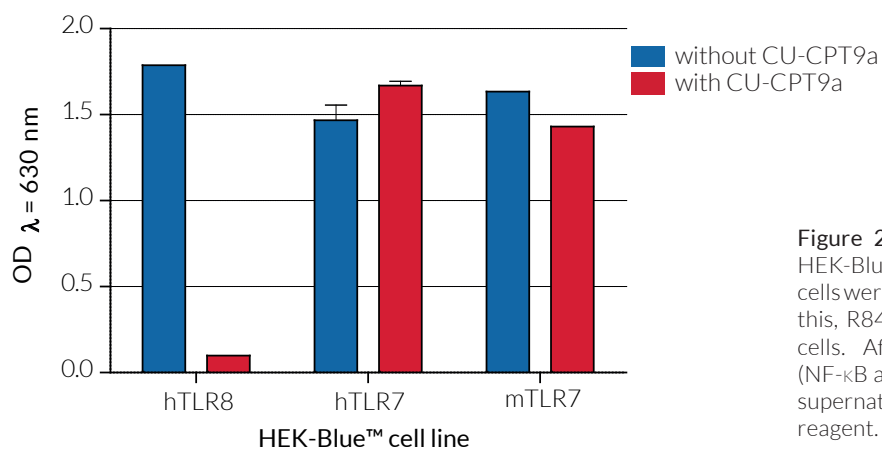


Figure 2: Specific inhibition of human TLR8 by Cu-CPT9a. HEK-Blue™ hTLR8, HEK-Blue™ hTLR7, and HEK-Blue™ mTLR7 cells were incubated with Cu-CPT9a (1 μM) for 3 hours. Following this, R848 (10 μg/ml), a TLR7/TLR8 agonist, was added to the cells. After overnight incubation, activation of TLR8 or TLR7 (NF-κB activity) was assessed by measuring SEAP activity in the supernatant, using QUANTI-Blue™ Solution, a SEAP detection reagent. Data are shown as optical density (OD) at 630nm.

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