

Validation data for CLI-095

<https://www.invivogen.com/cli-095>

For research use only

Version 22F14-AK

CLI-095 is a small-molecule inhibitor of the TLR4 signaling. In detail, CLI-095 attaches to cysteine 747 in the intracellular sphere of TLR4, thus blocking the intracellular domain. Subsequently, both MyD88-dependent and TRIF dependent pathways stimulated by lipopolysaccharide (LPS) are impaired. The ability of CLI-095 to suppress human (h)TLR4 signaling was validated using InvivoGen's HEK-Blue™ hTRL4 reporter cells (**Figure 1**). These cells stably express the TLR4, MD2 and CD14 genes as well as an NF- κ B-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene.

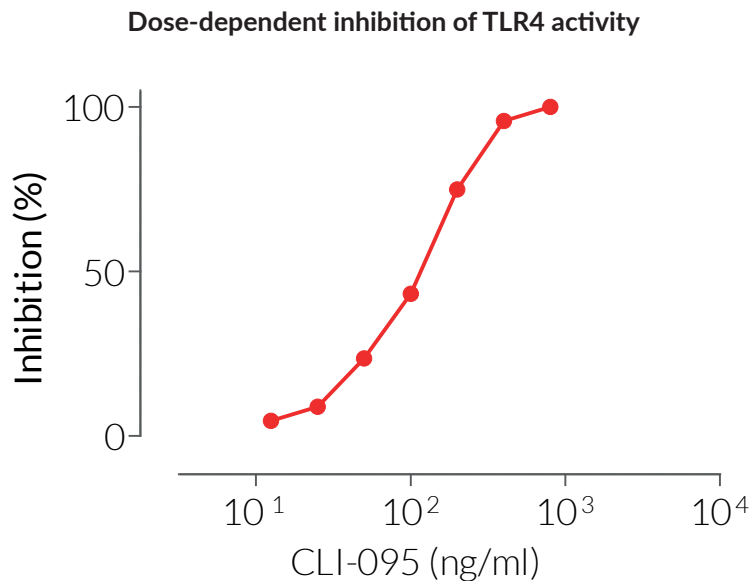


Figure 1: CLI-095 is a potent inhibitor of hTLR4 signaling pathway.

HEK-Blue™ hTLR4 cells were incubated in the presence of increasing concentrations of CLI-095 for 3 hours before adding 10 ng/ml LPS-EK Ultrapure (TLR4 agonist). After overnight incubation at 37°C, the neutralizing activity of CLI-095 was determined by measuring the reduction of SEAP production in the supernatant using the QUANTI-Blue™ Solution detection reagent. Data are shown in percentage (%) of inhibition.

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