

Validation data for ODN 2395 Control

<https://www.invivogen.com/odn2395-control>

For research use only

Version 24E06-AK

ODN 2395 Control is a synthetic oligonucleotide (ODN) designed as a negative control for the Toll-like receptor 9 (TLR9) agonist ODN 2395. ODN 2395 is a Class C CpG ODN specific for the human (h) and mouse (m) TLR9. In HEK-Blue™-derived reporter cells, ODN 2395 efficiently activates hTLR9 and mTLR9. As expected, ODN 2395 Control does not activate mTLR9 when compared to the immunostimulatory ODN 2395 (Figure 1). In addition, ODN 2395 Control does not activate hTLR9 (data not shown).

Dose-dependent NF- κ B response in HEK-Blue™ mTLR9 cells

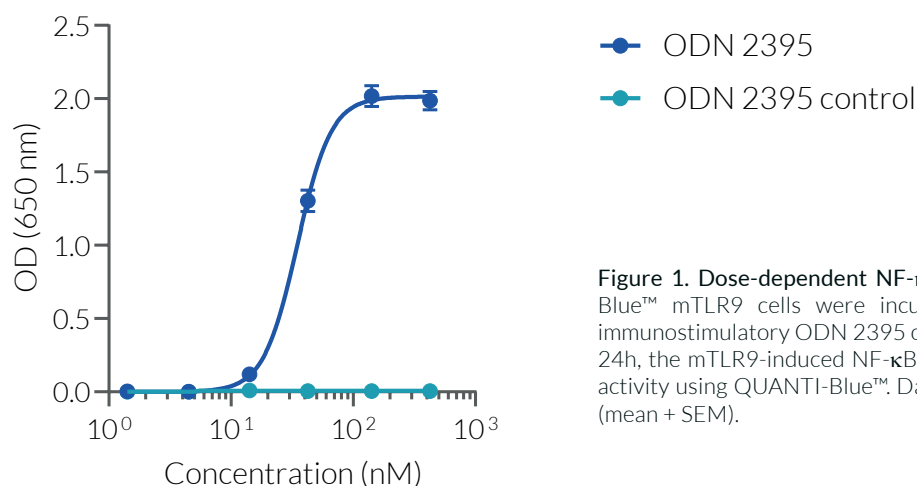


Figure 1. Dose-dependent NF- κ B response in HEK-Blue™ mTLR9 cells. HEK-Blue™ mTLR9 cells were incubated with increasing concentrations of the immunostimulatory ODN 2395 or the negative control ODN 2395 Control. After 24h, the mTLR9-induced NF- κ B response was assessed by measuring the SEAP activity using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean + SEM).

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