

# Validation data for CL264

<https://www.invivogen.com/cl264>

For research use only

Version 23L18-AK

CL264 is a 9-benzyl-8 hydroxyadenine derivative containing a glycine on the benzyl group (*in para*) and a potent agonist of the Toll-like receptor 7 (TLR7). The ability of CL264 to activate TLR7, but not TLR8 signaling was validated using a panel of InvivoGen's reporter cell lines. CL264 efficiently activates human (h) and mouse (m)TLR7, but not h/mTLR8, as assessed by the expression of an NF- $\kappa$ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter in HEK-Blue™-derived cell lines (Figure 1). The induction of the NF- $\kappa$ B and IRF pathways by CL264 has been tested using InvivoGen's HEK-Dual™ cells featuring two reporter genes, the NF- $\kappa$ B-inducible SEAP and IRF-inducible Lucia luciferase, as well as the overexpression of TLR7 or TLR8 (Figure 2).

## Dose-dependent NF- $\kappa$ B response in HEK-Blue™-derived cells

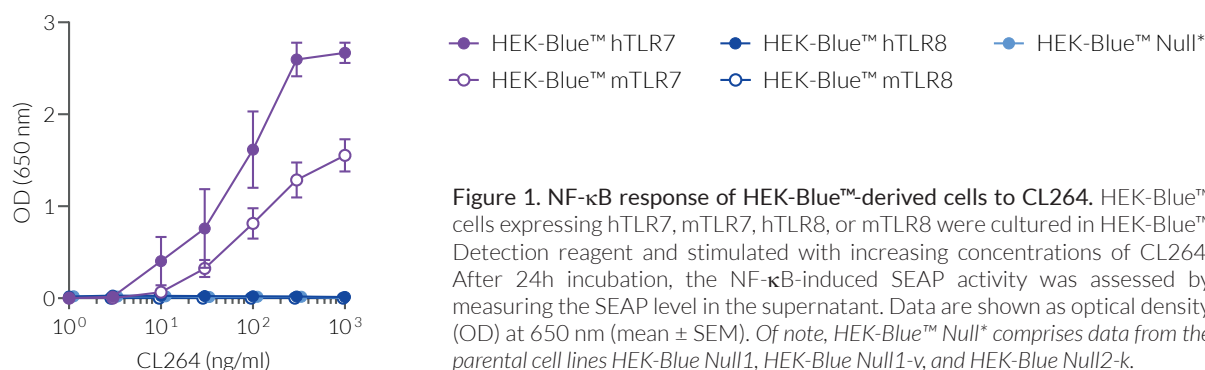


Figure 1. NF- $\kappa$ B response of HEK-Blue™-derived cells to CL264. HEK-Blue™ cells expressing hTLR7, mTLR7, hTLR8, or mTLR8 were cultured in HEK-Blue™. Detection reagent and stimulated with increasing concentrations of CL264. After 24h incubation, the NF- $\kappa$ B-induced SEAP activity was assessed by measuring the SEAP level in the supernatant. Data are shown as optical density (OD) at 650 nm (mean  $\pm$  SEM). Of note, HEK-Blue™ Null\* comprises data from the parental cell lines HEK-Blue Null1, HEK-Blue Null1-v, and HEK-Blue Null2-k.

## Dose-dependent NF- $\kappa$ B and IRF responses in HEK-Dual™-derived cells

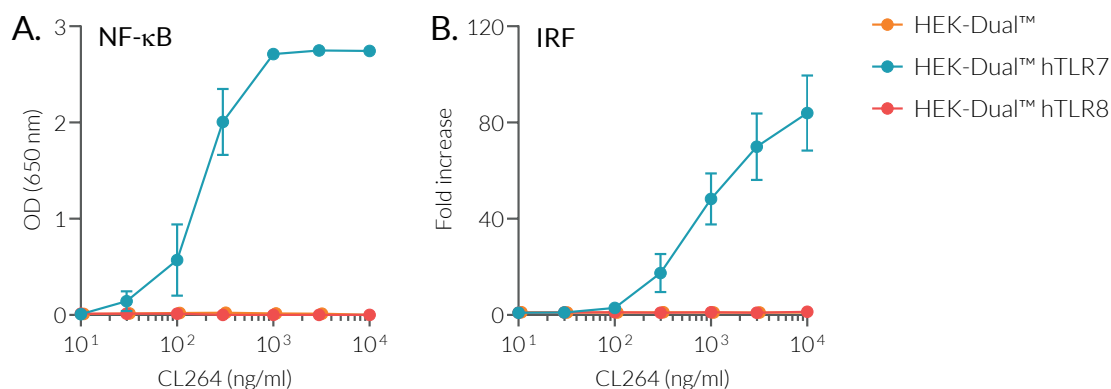


Figure 2. NF- $\kappa$ B and IRF responses of HEK-Dual™-derived cells to CL264. HEK-Dual™, HEK-Dual™ hTLR7, and HEK-Dual™ hTLR8 cells were incubated for 24 hours with increasing concentrations of CL264. After 24h incubation, the (A) NF- $\kappa$ B-induced SEAP activity was assessed using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean  $\pm$  SEM). (B) The IRF response was assessed by measuring the activity of Lucia luciferase in the supernatant using QUANTI-Luc™. Data are shown in fold response over non-induced cells (mean  $\pm$  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](mailto:info@invivogen.com)