

Validation data for CL075

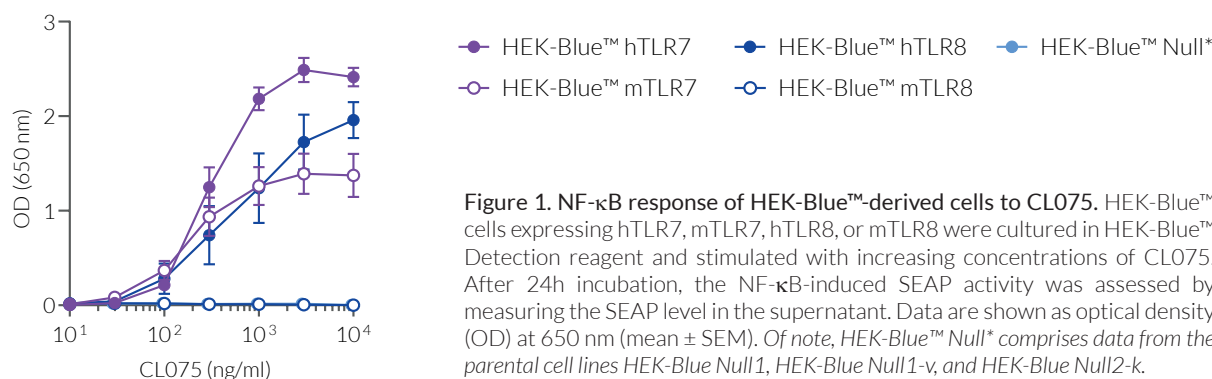
<https://www.invivogen.com/cl075>

For research use only

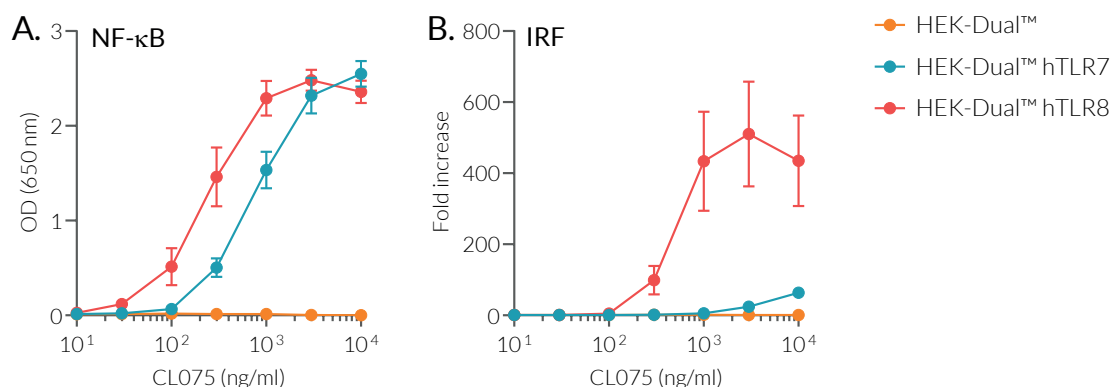
Version 23L18-AK

CL075 is a thiazoloquinolone derivative and a potent agonist of the Toll-like receptors 7 and 8 (TLR7 and TLR8). The ability of CL075 to activate TLR7 and TLR8 signaling was validated using a panel of InvivoGen's reporter cell lines. CL075 efficiently activates human (h)TLR7, murine (m)TLR7 and hTLR8, but not mTLR8, as assessed by the expression of an NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter in HEK-Blue™-derived cell lines (**Figure 1**). The induction of the NF- κ B and IRF pathways by CL075 has been tested using InvivoGen's HEK-Dual™ cells featuring two reporter genes, the NF- κ B-inducible SEAP and IRF-inducible Lucia luciferase, as well as the overexpression of TLR7 or TLR8 (**Figure 2**).

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



Dose-dependent NF- κ B and IRF responses in HEK-Dual™-derived cells



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