

Validation data for THP1-Dual™ KO-MyD cells

<https://www.invivogen.com/thp1-dual-komyd>

For research use only

Version 23A10-AK

THP1-Dual™ KO-MyD cells are engineered from the THP1-Dual™ cell line through the verified stable knockout of the human (h)MyD88 gene (Figure 1). These cells also feature two reporter genes, allowing the simultaneous study of NF-κB- and IRF-induced responses by monitoring the SEAP (secreted embryonic alkaline phosphatase) and Lucia luciferase activities. Lucia luciferase and SEAP activities are readily assessable in the cell culture supernatant using QUANTI-Luc™ and QUANTI-Blue™ Solution detection reagents, respectively. THP1-Dual™ KO-MyD cells display significantly reduced responses to ligands that activate receptors whose signaling is MyD88-dependent, such as TLR1/2, TLR2/6, TLR4, TLR5, TLR8 and the interleukin-1 receptor (IL-1R). These cells remain responsive to ligands that activate receptors whose signaling is MyD88-independent, such as the tumor necrosis factor (TNF) receptor or NOD1 (Figure 2).

Validation of MyD88 Knockout

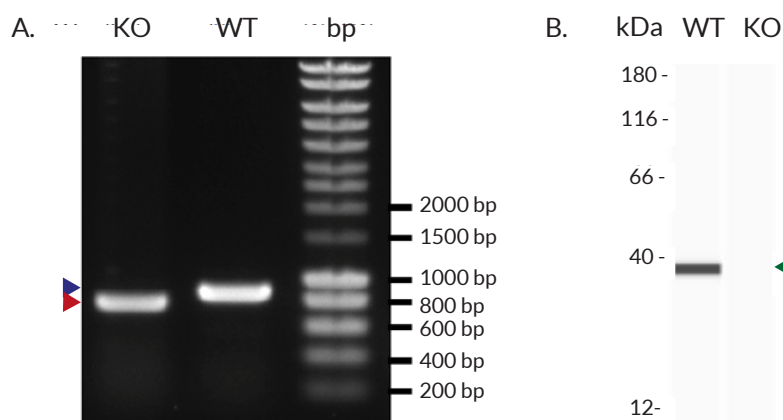


Figure 1: Validation of MyD88 knockout in THP1-Dual™ KO-MyD cells. (A) The targeted MyD88 region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-MyD (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-MyD cells feature a biallelic deletion. (B) Lysates from THP1-Dual™ (WT) and THP1-Dual™ KO-MyD (KO) cells were analyzed by Western blot (Wes™) using an anti-human MyD88 antibody, followed by an HRP conjugated anti-rabbit secondary antibody. The green arrow indicates the expected band for the MyD88 protein (33 kDa).

Functional validation of MyD88 knockout (NF-κB response)

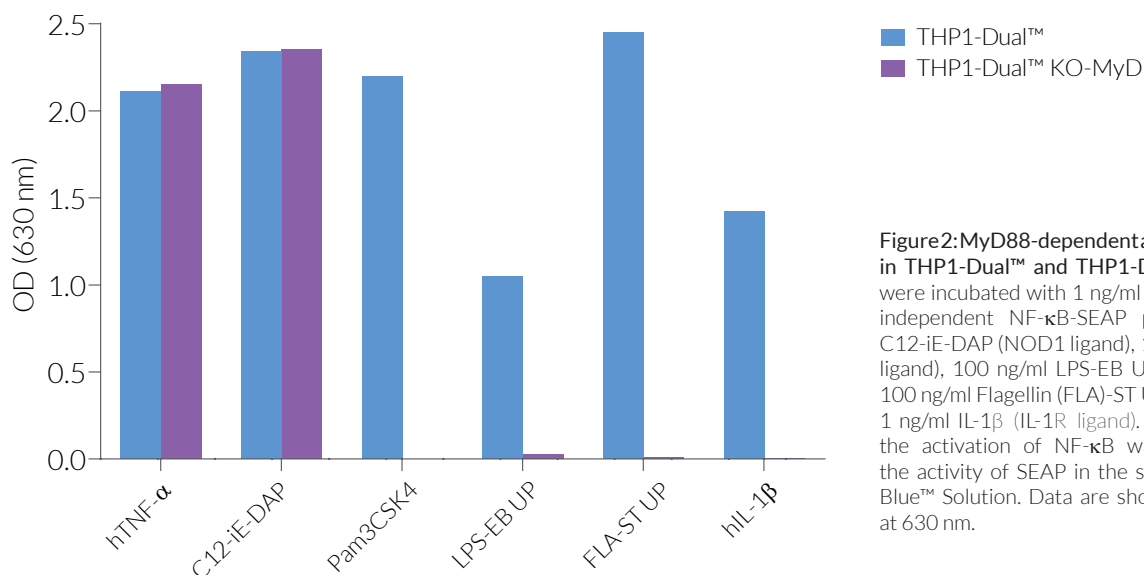


Figure 2: MyD88-dependent and independent responses in THP1-Dual™ and THP1-Dual™ KO-MyD cells. Cells were incubated with 1 ng/ml human (h) TNF-α (MyD88-independent NF-κB-SEAP positive control), 1 μg/ml C12-iE-DAP (NOD1 ligand), 1 ng/ml Pam3CSK4 (TLR1/2 ligand), 100 ng/ml LPS-EB Ultrapure (UP; TLR4 ligand), 100 ng/ml Flagellin (FLA)-ST Ultrapure (TLR5 ligand), and 1 ng/ml IL-1β (IL-1R ligand). After overnight incubation, the activation of NF-κB was assessed by measuring the activity of SEAP in the supernatant using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 nm.

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