

# Validation data for THP1-Dual™ hTLR3 cells

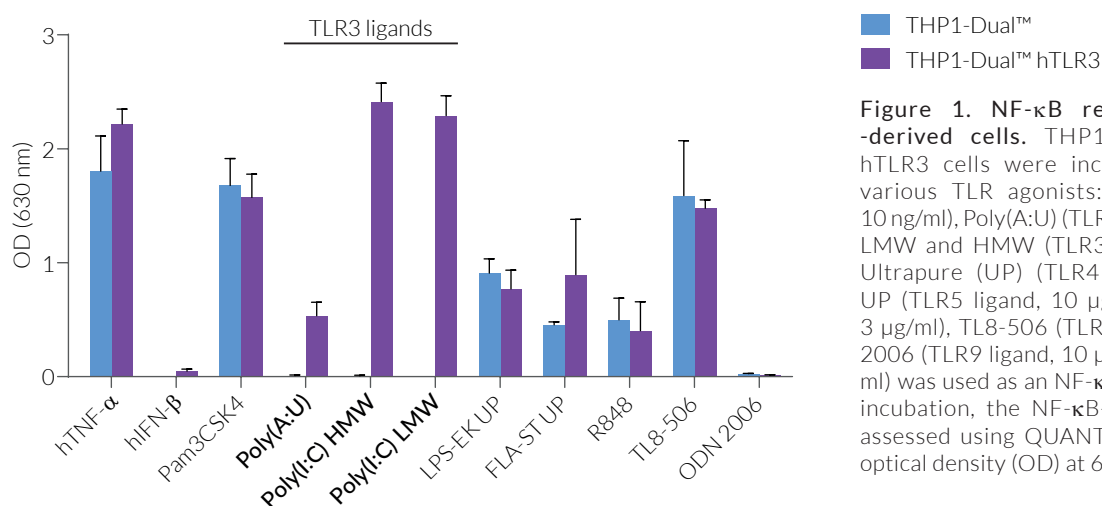
<https://www.invivogen.com/thp1-dual-hltr3>

For research use only

Version 23C24-AK

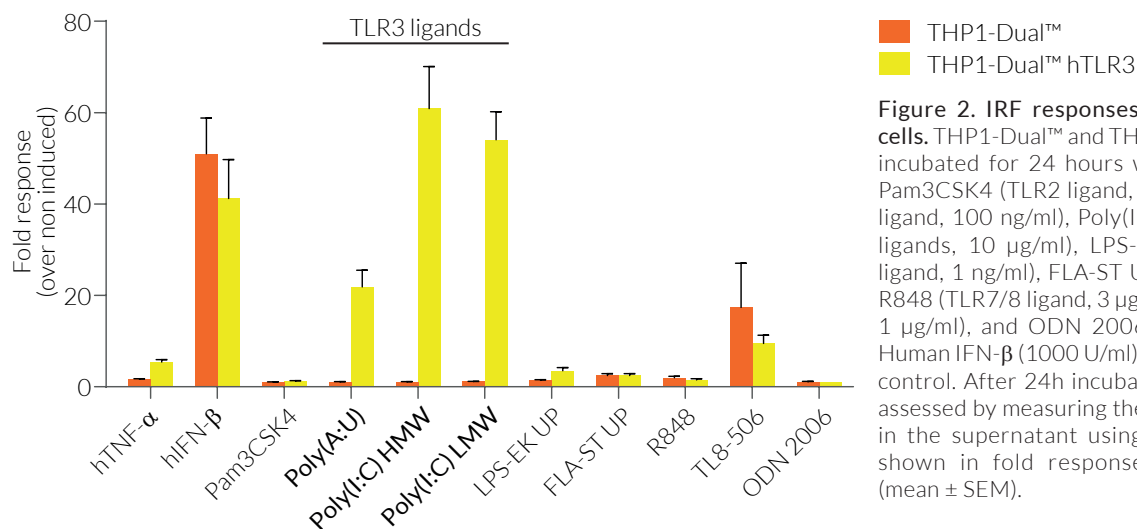
THP1-Dual™ hTLR3 cells were generated from the THP1-Dual™ cell line through the stable expression of the human Toll-like receptor 3 (hTLR3) as well as elements important for its signaling. These cells also feature two reporter genes allowing the simultaneous study of NF-κB- and IRF-induced responses, by monitoring the SEAP (secreted embryonic alkaline phosphatase) and Lucia luciferase activities, respectively. Due to the stable expression of hTLR3, these cells show strong NF-κB and IRF responses upon incubation with synthetic double-stranded (ds)RNA analogs, such as Poly(A:U) (polyadenylic-polyuridylic acid) or Poly(I:C) (polyinosinic-polycytidylic acids), when compared to their parental cells (Figures 1 & 2). Of note, as THP-1 cells express endogenous levels of various TLRs, THP1-Dual™ - derived cells respond to the cognate ligands including Pam3CSK4, LPS, or flagellin.

## Functional validation of THP1-Dual™-derived cells (NF-κB responses)



**Figure 1. NF-κB responses in THP1-Dual™-derived cells.** THP1-Dual™ and THP1-Dual™ hTLR3 cells were incubated for 24 hours with various TLR agonists: Pam3CSK4 (TLR2 ligand, 10 ng/ml), Poly(A:U) (TLR3 ligand, 100 ng/ml), Poly(I:C) LMW and HMW (TLR3 ligands, 10 µg/ml), LPS-EK Ultrapure (UP) (TLR4 ligand, 10 ng/ml), FLA-ST UP (TLR5 ligand, 10 µg/ml), R848 (TLR7/8 ligand, 3 µg/ml), TL8-506 (TLR8 ligand, 1 µg/ml), and ODN 2006 (TLR9 ligand, 10 µg/ml). Human TNF-α (10 ng/ml) was used as an NF-κB-positive control. After 24h incubation, the NF-κB-induced SEAP activity was assessed using QUANTI-Blue™. Data are shown as optical density (OD) at 630 nm (mean ± SEM).

## Functional validation of THP1-Dual™-derived cells (IRF response)



**Figure 2. IRF responses in THP1-Dual™-derived cells.** THP1-Dual™ and THP1-Dual™ hTLR3 cells were incubated for 24 hours with various TLR agonists: Pam3CSK4 (TLR2 ligand, 10 ng/ml), Poly(A:U) (TLR3 ligand, 100 ng/ml), Poly(I:C) LMW and HMW (TLR3 ligands, 10 µg/ml), LPS-EK Ultrapure (UP) (TLR4 ligand, 1 ng/ml), FLA-ST UP (TLR5 ligand, 10 µg/ml), R848 (TLR7/8 ligand, 3 µg/ml), TL8-506 (TLR8 ligand, 1 µg/ml), and ODN 2006 (TLR9 ligand, 10 µg/ml). Human IFN-β (1000 U/ml) was used as an IRF-positive control. After 24h incubation, 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 ± 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)