

# Validation data for THP1-Dual™ cells

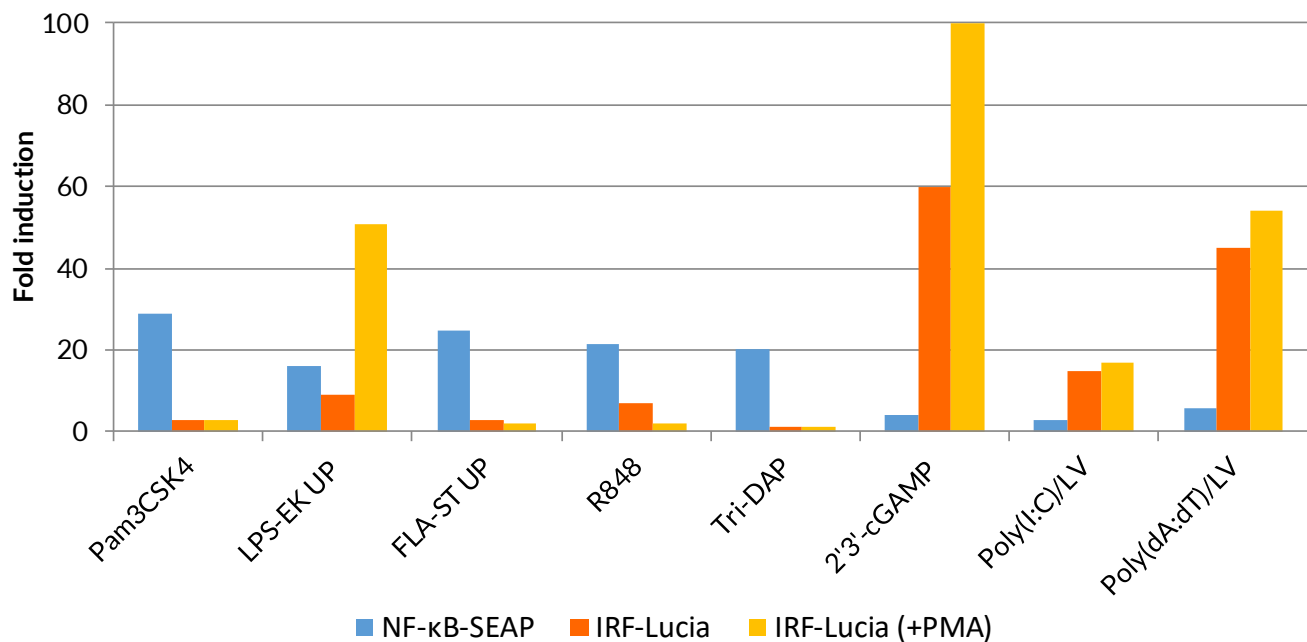
<https://www.invivogen.com/thp1-dual>

For research use only

Version 20B26-MM

THP1-Dual™ cells are derived from the human THP-1 monocyte cell line by stable integration of two inducible secreted reporter genes: Lucia luciferase and SEAP (secreted embryonic alkaline phosphatase). As a result, they allow the simultaneous study of the NF-κB pathway, by monitoring the activity of SEAP, and the IRF (interferon regulatory factor) pathway, by assessing the activity of Lucia luciferase. THP1-Dual™ cells induce activation of NF-κB in response to certain TLR agonists, such as Pam3CSK4 and flagellin. They trigger the IRF pathway upon stimulation with type I interferons (IFNs) and RLR (RIG-I-like receptor), CDS (cytosolic dsDNA sensor) or STING agonists, such as transfected poly(I:C) or poly(dA:dT) and 2'3'cGAMP respectively.

## NF-κB and IRF responses of THP1-Dual™ Cells to PRR ligands



**Figure 1: NF-κB/IRF dual response of THP1-Dual™ cells.** Cells were pretreated or not pretreated with PMA (Phorbol 12-myristate 13-acetate; 20 ng/ml for 3 hours) and stimulated with 1 ng/ml Pam3CSK4 (TLR2), 100 ng/ml LPS-EK UP (TLR4), 100 ng/ml FLA-ST UP (TLR5), 10 µg/ml R848 (TLR7/8), 10 µg/ml Tri-DAP (NOD1), 3 µg/ml 2'3'-cGAMP (STING), 1 µg/ml poly(I:C)/LyoVec™ (RLR) or 100 ng/ml poly(dA:dT)/LyoVec™ (CDS). After 24h incubation, NF-κB and IRF activation were assessed by measuring the levels of SEAP and Lucia luciferase using QUANTI-Blue™ and QUANTI-Luc™, respectively. With QUANTI-Blue™ the levels of SEAP were determined by reading the optical density (OD) at 655 nm. With QUANTI-Luc™ the levels of Lucia luciferase were determined by measuring the relative light units (RLUs) in a luminometer.

### 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 Hong Kong: +852 3622-34-80  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)