

Validation data for HEK-Blue™ mNOD2 cells

<https://www.invivogen.com/hek-blue-mnod2>

For research use only

Version 23G24-AK

HEK-Blue™ mNOD2 cells were engineered from the human embryonic kidney HEK 293 cell line to assess the role of the Nucleotide-binding Oligomerization Domain-containing protein 2 (NOD2). These cells stably express the murine (m) *NOD2* gene and an NF- κ B-inducible SEAP reporter gene. A strong SEAP activation upon stimulation of NOD2 using the specific ligands L18-MDP (Muramyl dipeptide), MDP, and Murabutide can be readily monitored by performing the assay in HEK-Blue™ Detection medium (Figure 1). HEK293 cells express endogenous levels of various pattern recognition receptors (PRRs), including NOD1, TLR3, and TLR5. Therefore, HEK-Blue™ mNOD2 cells also respond to their cognate ligands Tri-DAP, Poly(I:C), and flagellin, respectively (Figure 2).

Dose-dependent NF- κ B responses of HEK-Blue™ mNOD2 to NOD2-specific ligands

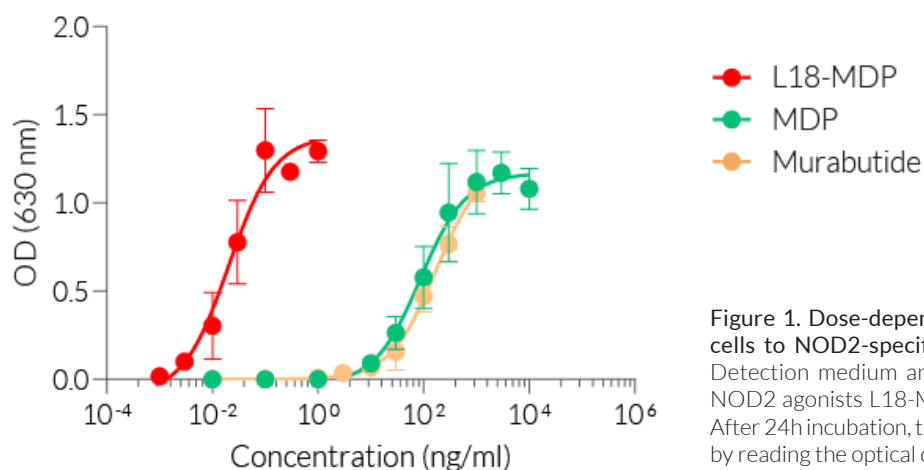


Figure 1. Dose-dependent NF- κ B responses of HEK-Blue™ mNOD2 cells to NOD2-specific ligands. Cells were incubated in HEK-Blue™ Detection medium and stimulated with increasing concentrations of NOD2 agonists L18-MDP (Muramyl dipeptide), MDP, and Murabutide. After 24h incubation, the levels of NF- κ B-induced SEAP were determined by reading the optical density (OD) at 630 nm (mean \pm SEM).

NF- κ B response of HEK-Blue™ mNOD2 cells to various PRR agonists and cytokines

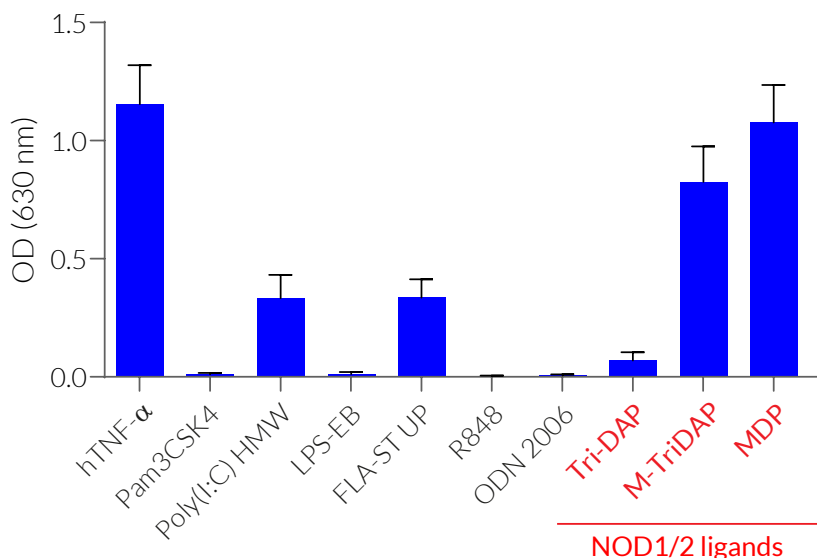


Figure 2. NF- κ B responses of HEK-Blue™ mNOD2 cells to various PRR agonists and cytokines. Cells were incubated in HEK-Blue™ Detection medium and stimulated for 24h with various cytokines and PRR agonists: Human (h)TNF- α (NF- κ B-positive control, 10 ng/ml), Pam3CSK4 (TLR2 ligand, 1 μ g/ml), Poly(I:C) HMW (TLR3 ligand, 1 μ g/ml), LPS-EK Ultrapure (UP) (TLR4 ligand, 10 μ g/ml), FLA-ST UP (TLR5 ligand, 1 μ g/ml), R848 (TLR7/8 ligand, 10 μ g/ml) ODN 2006 (TLR9 ligand, 10 μ g/ml), Tri-DAP (NOD1 ligand, 1 μ g/ml), M-Tri-DAP (NOD1/2 ligand, 1 μ g/ml), and MDP (NOD2 ligand, 1 μ g/ml). After 24h incubation, the levels of NF- κ B-induced SEAP were determined by reading the optical density (OD) at 630 nm (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