

Validation data for ADP-Heptose

<https://www.invivogen.com/adp-heptose>

For research use only

Version 19K26-ED

InvivoGen has synthesized and purified ADP-L-*glycero*-b-D-*manno*-heptose (ADP-Heptose), an intermediary sugar in the biosynthesis of lipopolysaccharide (LPS), an essential component of the outer membrane of Gram-negative bacteria. ADP-Heptose has been identified as a potent pathogen associated molecular pattern (PAMP) that binds to the cytosolic pattern recognition receptor (PRR) ALPK1 and activates the ALPK1-TIFA signaling pathway. Ultimately, ADP-Heptose induces a pro-inflammatory NF- κ B-dependent response in host cells (Figure 1).

Functional validation of ADP-Heptose

Stimulation of InvivoGen's HEK-Blue™ Null1-v cells (red), which express a NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP), with ADP-Heptose results in a clear dose-dependent response. Furthermore, this activation is dependent upon ALPK1 and TIFA, with no response observed in the derived knockout (KO) cell lines, HEK-Blue™ KO ALPK1 (yellow) and HEK-Blue™ KO-TIFA (blue) cells.

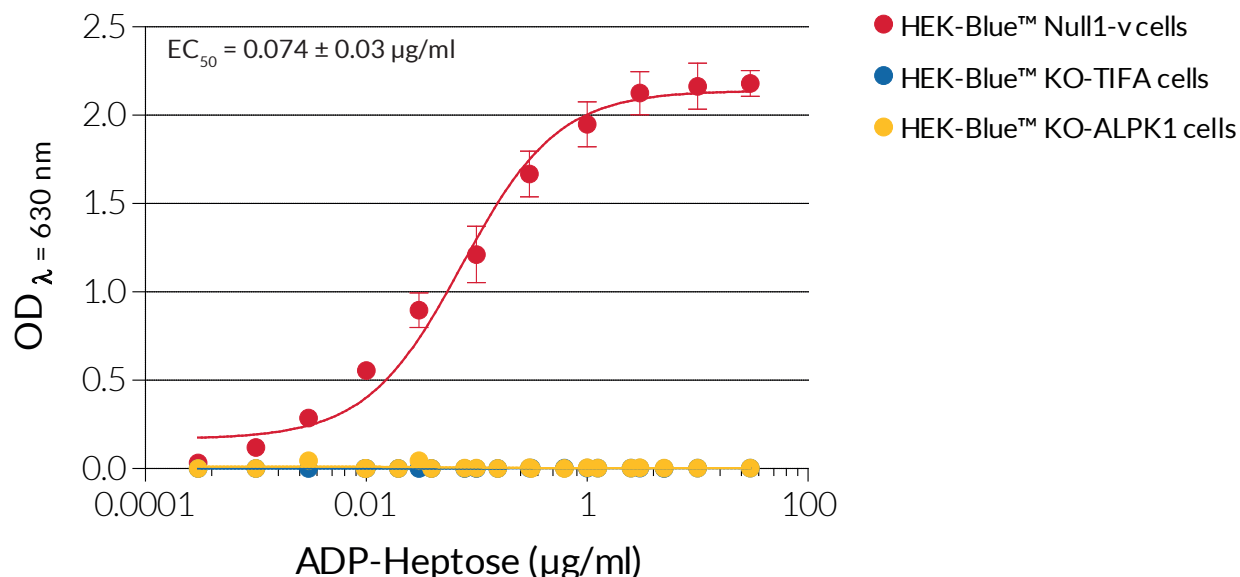


Figure 1: ADP-Heptose induced NF- κ B response. HEK-Blue™ Null1-v, HEK-Blue™ KO-ALPK1, and HEK-Blue™ KO-TIFA cells were incubated with increasing concentrations of ADP-Heptose (0 - 30 $\mu\text{g/ml}$). After overnight incubation, the NF- κ B response was assessed by measuring the activity of SEAP in the supernatant using QUANTI-Blue™ Solution, a SEAP detection reagent. Data are presented as optical density (OD) at 630 nm (mean \pm SEM). EC_{50} value is indicated (\pm std error).

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 3-622-34-80
E-mail: info@invivogen.com