

# Validation data for ADP-L-Heptose

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

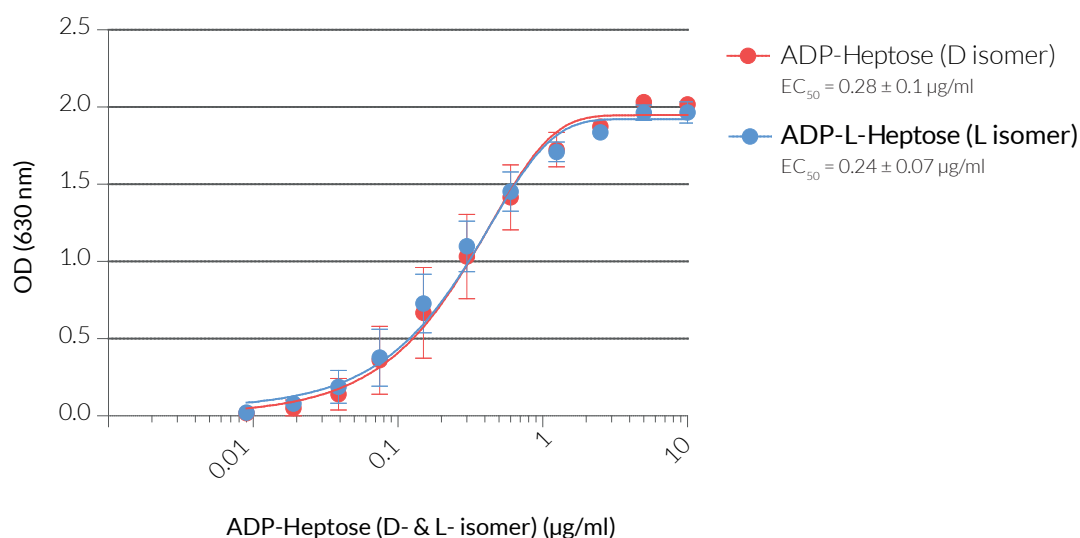
For research use only

Version 21F08-ED

InvivoGen has synthesized and purified ADP-L-Heptose, an intermediary sugar in the biosynthesis of lipopolysaccharide (LPS), an essential component of the outer membrane of Gram-negative bacteria. It is generated by a multi-step biosynthesis pathway, in which the final step is the interconversion between two isomers, ADP-D-glycero- $\beta$ -D-manno-heptose (ADP-Heptose) and ADP-L-glycero- $\beta$ -D-manno-heptose (ADP-L-Heptose). Both the D- and L- isomers activate the ALPK1-TIFA signaling cascade and induce a pro-inflammatory NF- $\kappa$ B-dependent response (Figure 1).

## Functional validation of ADP-L-Heptose

Induction of InvivoGen's HEK-Blue™ Null1-v cells, which express a NF- $\kappa$ B-inducible secreted embryonic alkaline phosphatase (SEAP), with either ADP-Heptose (D- isomer) or ADP-L-Heptose (L- isomer) results in comparable dose-dependent responses.



**Figure 1: ADP-L-Heptose induced NF- $\kappa$ B response.** HEK-Blue™ Null1-v were incubated with increasing concentrations of ADP-Heptose or ADP-L-Heptose (0 - 10 µg/ml). After overnight incubation, the NF- $\kappa$ 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<sub>50</sub> 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 Asia: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)