

# Validation data for ATP

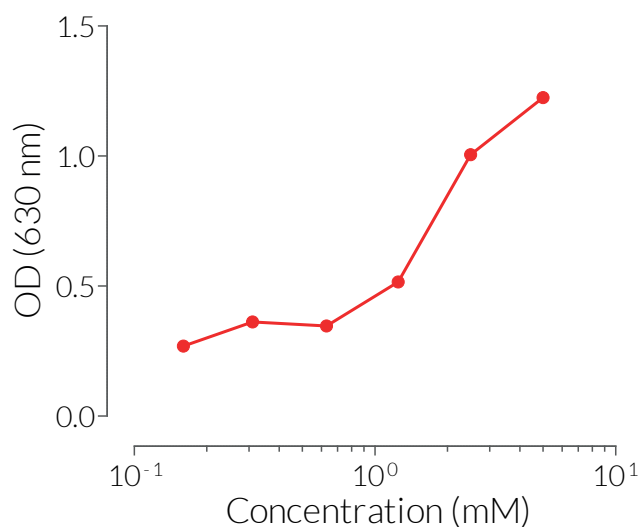
<https://www.invivogen.com/atp>

For research use only

Version 22H26-AK

The organic compound adenosine triphosphate (ATP) is a key potassium (K<sup>+</sup>) efflux agent, that induces NLRP3 inflammasome formation. The biological activity of ATP has been validated using THP1-Null cells. The production of IL-1 $\beta$  by THP1-Null cells was measured using HEK-Blue™ IL-1 $\beta$  cells. Treatment with ATP induced IL-1 $\beta$  secretion, an indicator of NLRP3 inflammasome activation, in a dose-dependent manner.

## Dose-dependent IL-1 $\beta$ induction



IL-1 $\beta$  production in THP1-Null cells. Human THP-1 monocytes were primed for 3 hours with LPS-EK(10  $\mu$ g/ml) prior to the incubation with increasing concentrations of ATP. The next day, the supernatant was incubated with HEK-Blue™ IL-1 $\beta$  cells for 24 hours and IL-1 $\beta$  was assessed using QUANTI-Blue™ Solution. The optical density (OD) was read at 630 nm.

### 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)