

Validation data for Alum Hydroxide

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

For research use only

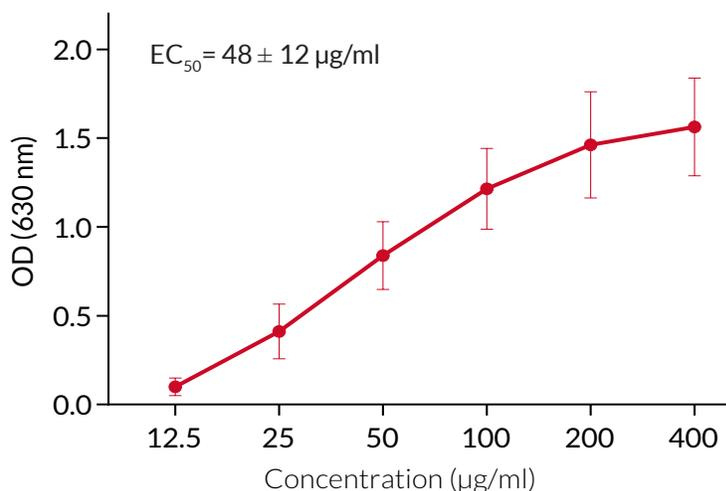
Version 19L06-MM

Aluminium Hydroxide (alum) is an inducer of the NLRP3 inflammasome, a large intracellular multiprotein complex that plays a central role in innate immunity^{1,2}. NLRP3 inflammasome activation requires an initial signal ('priming'), provided by microbial molecules, such as lipopolysaccharide (LPS), and a secondary signal, provided by a wide array of stimuli including bacterial toxins, endogenous molecules or crystalline substances such as alum. This triggers the multimerization of the NLRP3 inflammasome and caspase-1 activation with the subsequent maturation and secretion of IL-1 β and IL-18.

InvivoGen's Alum Hydroxide is designed for *in vitro* assays. Its ability to induce the NLRP3 inflammasome has been validated using THP1-Null cells. The production of IL-1 β by THP1-Null cells was measured using HEK-Blue™ IL-1 β cells. Treatment with Alum Hydroxide induced IL-1 β secretion, an indicator of NLRP3 inflammasome activation, in a dose-dependent manner.

1. Schroder K. & Tschopp J., 2010. The inflammasomes. Cell 140(6):821-32. 2. Franchi L. et al., 2012. Sensing and reacting to microbes through the inflammasomes. Nat Immunol 13(4):325-32.

Evaluation of NLRP3 inflammasome activation



IL-1 β production in THP1-Null cells. THP1-Null cells, primed with LPS (1 µg/ml for 3h), were stimulated with increasing concentrations of Alum Hydroxide. After overnight incubation, IL-1 β secretion was analyzed by adding 50 µl of supernatant from treated THP1-Null cells to HEK-Blue™ IL-1 β cells. IL-1 β -induced activation of NF- κ B was assessed by measuring the levels of SEAP in the supernatant of HEK-Blue™ IL-1 β cells using QUANTI-Blue™ Solution, a SEAP detection reagent, and by reading the optical density (OD) 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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com