

Validation data for Cholesterol Crystals

<https://www.invivogen.com/cholesterol-crystals-nlrp3-inflammasome-inducer>

For research use only

Version 25D04-NJ

Cholesterol crystals are the hallmark of atherosclerosis. They strongly stimulate the NLRP3 inflammasome-dependent induction of pro-inflammatory cytokines, including interleukin 1 β (IL-1 β). The ability of InvivoGen's cholesterol crystals to induce the NLRP3 inflammasome has been validated using THP1-Null2 and THP1 KO-NLRP3 cells. The production and release of IL-1 β was measured using HEK-Blue™ IL-1 β cells. The treatment with cholesterol crystals induced IL-1 β secretion in a dose-dependent manner (Figure 2).

Microscopic evaluation of cholesterol crystals

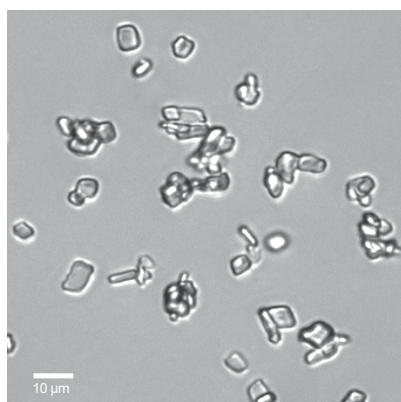


Figure 1. Microscopic analysis of Cholesterol Crystals. Cholesterol crystal particles resuspended in sterile water.

NLRP3 inflammasome activation by cholesterol crystals

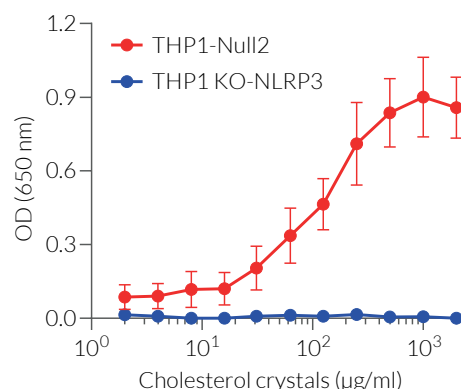


Figure 2. NLRP3-dependent IL-1 β secretion upon treatment with cholesterol crystals. THP1-Null2 and THP1 KO-NLRP3 cells were primed for 3 hours with LPS-EK (1 μ g/ml) prior to incubation with increasing concentrations of cholesterol crystals. After overnight incubation, the IL-1 β secretion was assessed in the culture supernatant using HEK-Blue™ IL-1 β sensor cells and the SEAP detection reagent QUANTI-Blue™ Solution. Data is shown as optical density (OD) at 650 nm (mean \pm SD).

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