

# Validation data for Isoliquiritigenin

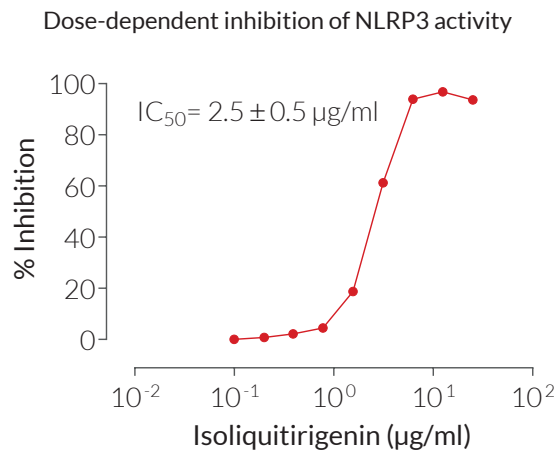
<https://www.invivogen.com/isoliquiritigenin>

For research use only

Version 21A25-MM

Isoliquiritigenin is an inhibitor of NF- $\kappa$ B and the NLRP3 (NOD-like receptor (NLR) pyrin domain-containing protein 3) inflammasome. NF- $\kappa$ B regulates multiple aspects of innate and adaptive immune functions and serves as a pivotal mediator of inflammatory responses. While the NLRP3 inflammasome is an innate immune sensor that is activated by a two-step process; a first signal ('priming') is provided by microbial molecules such as lipopolysaccharide (LPS), and the second signal is provided by a wide array of stimuli including endogenous molecules or crystalline substances such as monosodium urate (MSU) crystals. Inflammasome activation triggers caspase-1-mediated interleukin-1 $\beta$  (IL-1 $\beta$ ) production and secretion.

The ability of isoliquiritigenin to inhibit the NLRP3 inflammasome was validated using InvivoGen's THP-1/HEK-Blue™ IL-1 $\beta$  assay. This assay uses the secretion of IL-1 $\beta$  by THP1-Null2 cells as an indicator of NLRP3 inflammasome induction. The IL-1 $\beta$  production by these cells is measured using HEK-Blue™ IL-1 $\beta$  cells. Treatment with isoliquiritigenin inhibited IL-1 $\beta$  secretion in a dose-dependent manner (**Figure 1**).



**Figure 1:** Isoliquiritigenin inhibits the NLRP3 inflammasome response in a dose-dependent manner.

THP1-Null2 cells, primed with LPS-EK (1 µg/ml for 3 h), were stimulated with MSU (150 µg/ml) and increasing concentrations of isoliquiritigenin. After overnight incubation, IL-1 $\beta$  secretion was analyzed by adding 50 µl of supernatant from treated THP1-Null2 cells to HEK-Blue™ IL-1 $\beta$  cells. IL-1 $\beta$ -induced activation of NF- $\kappa$ B was assessed by measuring the levels of SEAP in the supernatant of HEK-Blue™ IL-1 $\beta$  cells using QUANTI-Blue™ Solution, a SEAP detection reagent, and by reading the optical density (OD) at 655 nm. Data are shown as a percentage (%) inhibition of the maximal response for the ligand with no inhibitor.

## 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](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)