

Validation data for Bay11-7082

<https://www.invivogen.com/bay11-7082>

For research use only

Version 22F07-MM

Bay11-7082 is a potent inhibitor of the transcription factor NF- κ B (nuclear factor kappa-light-chain-enhancer of activated B cells) and the NLRP3 (NOD-like receptor (NLR) pyrin domain-containing protein 3) inflammasome. NF- κ 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.

The ability of Bay11-7082 to inhibit the NLRP3 inflammasome was validated using InvivoGen's THP-1/HEK-Blue™ IL-1 β assay. This assay uses the secretion of IL-1 β by THP1-Null2 cells as an indicator of NLRP3 inflammasome induction. The IL-1 β production by these cells is measured using HEK-Blue™ IL-1 β cells. Treatment with Bay11-7082 inhibited IL-1 β secretion in a dose-dependent manner (Figure 1).

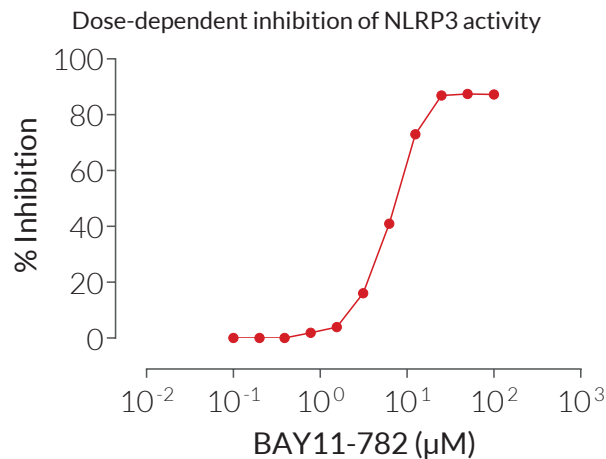


Figure 1: Bay11-7082 inhibits the NLRP3 inflammasome response in a dose-dependent manner.

THP1-Null2 cells were primed with LPS-EK (1 μ g/ml) for 3 h and then stimulated with MSU (150 μ g/ml) and increasing concentrations of Bay11-7082. After overnight incubation, IL-1 β secretion was analyzed by adding 50 μ l of supernatant from treated THP1-Null2 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 655 nm. Data are shown as percentage (%) inhibition.

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