

Validation data for Bafilomycin A1

<https://www.invivogen.com/BafilomycinA1>

For research use only

Version 21F28-MM

Bafilomycin A1 (BafA1) is a specific vacuolar H⁺ ATPase (V-ATPase) inhibitor. Due to its ability to specifically target V-ATPase and hence disrupt autophagic flux, BafA1 is frequently used to study autophagy and endosomal acidification. Notably, by inhibiting V-ATPase, BafA1 can be used to block the activation of nucleic acid sensing endosomal Toll-Like receptors (TLRs), such as TLR9, by neutralizing endosomal pH.

The ability of BafA1 to inhibit V-ATPase was validated using InvivoGen's HEK-Blue™ hTLR9 cells. These cells were specifically designed for the study of human TLR9-induced NF-κB signaling by monitoring the activity of secreted embryonic alkaline phosphatase (SEAP) reporter activity. The SEAP production by these cells was measured using QUANTI-Blue™ Solution, a SEAP detection reagent. Treatment with BafA1 inhibited SEAP activity in a dose-dependent manner (Figure 1).

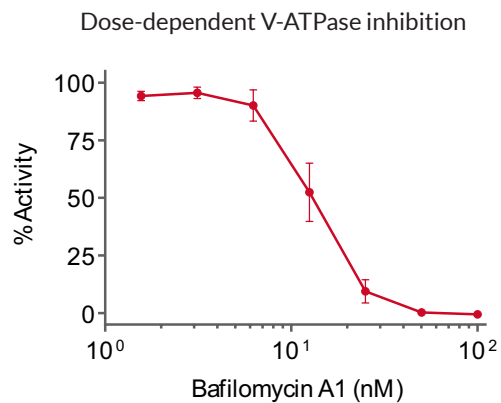


Figure 1: Bafilomycin A1 inhibits the TLR9 response in a dose-dependent manner.

HEK-Blue™ hTLR9 cells were stimulated with ODN 2006 (0.3 µg/ml) and increasing concentrations of BafA1. After overnight incubation, BafA1-induced inhibition of TLR9-NF-κB signaling was assessed by measuring the levels of SEAP using QUANTI-Blue™ Solution and by reading the optical density (OD) at 630 nm. Data are shown as percentage (%) activity ± SEM.

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