

Validation data for Brefeldin A

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

For research use only

Version 19G30-ED

Brefeldin A (BFA) is a small, potent, and reversible inhibitor of intracellular protein trafficking and vesicle formation between the endoplasmic reticulum (ER) and the Golgi apparatus. As ER-Golgi vesicle trafficking plays a pivotal role in the secretion of many proteins, BFA is commonly utilized as an inhibitor of protein secretion in cellular assays (Figure 1). Additionally, it has been established that BFA inhibits type I interferon (IFN) production by blocking the dissociation of activated STING from the ER (Figure 2). Of note, while morphological changes are seen in THP1 cells when BFA is used at concentrations above 0.6 μM , there is no change to the cellular viability.

Inhibition of STING-induced IRF activity by Brefeldin A

Treatment of THP1-Dual™ cells with BFA for 1 hour results in reduced levels of an IRF-inducible Lucia luciferase in a dose-dependent manner upon incubation with the STING ligands 2'3'-cGAMP and cAIM(PS)2 Difluor (Rp/Sp). Similarly, reduced NF- κ B-inducible SEAP (secreted alkaline phosphatase) is also observed in these cells (data not shown).

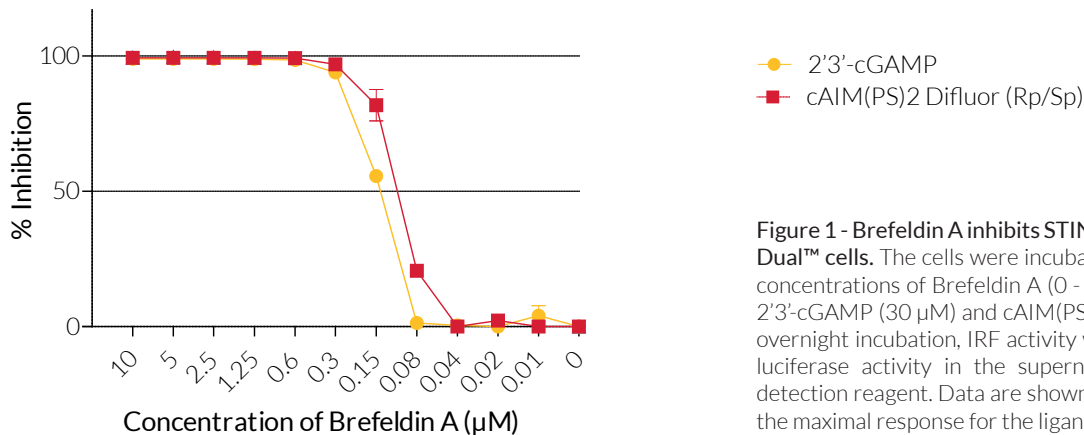


Figure 1 - Brefeldin A inhibits STING-induced IRF activity in THP1-Dual™ cells. The cells were incubated in the presence of increasing concentrations of Brefeldin A (0 - 10 μM) for 1 hour before adding 2'3'-cGAMP (30 μM) and cAIM(PS)2 Difluor (Rp/Sp) (10 μM). After overnight incubation, IRF activity was assessed by measuring Lucia luciferase activity in the supernatant, using the QUANTI-Luc™ detection reagent. Data are shown as a percentage (%) inhibition of the maximal response for the ligand with no BFA.

Inhibition of constitutive STING-induced IRF activity by Brefeldin A

BFA prevents the movement of STING to the ER-Golgi intermediate compartment, where it activates the TBK1-IRF3 signaling axis, and ultimately triggers expression of IFNs. Upon incubation with BFA for 18 hours, reduced levels of an IRF-inducible Lucia luciferase in a dose-dependent manner is observed in THP1-Dual™ KI-hSTING-S154 (constitutively active STING variant) cells. This cell line displays constitutive activation of STING and expression of an IRF-dependent Lucia luciferase in the absence of an inducer, due to the knock-in of a 'gain-of-function' mutation.

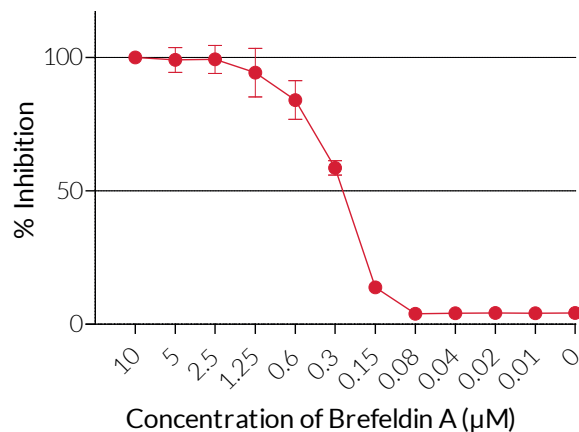


Figure 2 - Brefeldin A inhibits constitutively activated STING in THP1-Dual™ KI-hSTING-S154 cells. The cells were incubated in the presence of increasing concentrations of Brefeldin A (0 - 10 μM) for 18 hours. Inhibition of STING signaling was assessed by measuring the activity of the constitutively expressed Lucia luciferase in the supernatant, using the QUANTI-Luc™ detection reagent. Data are shown as a percentage (%) inhibition of the maximal response of the cell line with no BFA.

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