

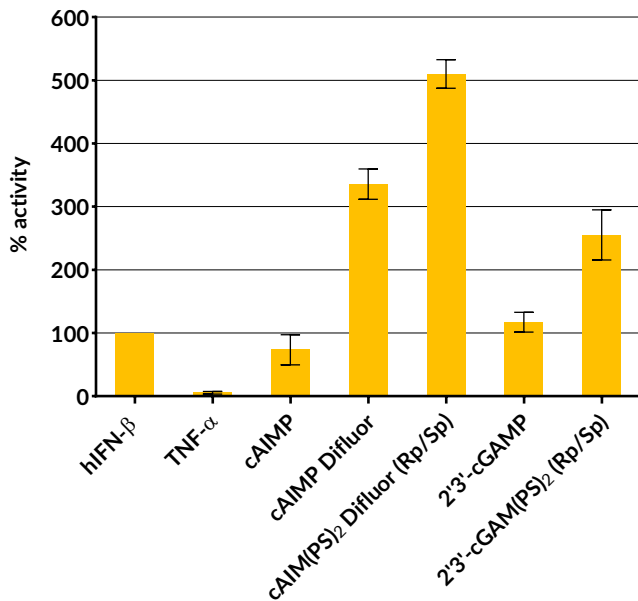
Validation data for cAIMP and its analogs

For research use only

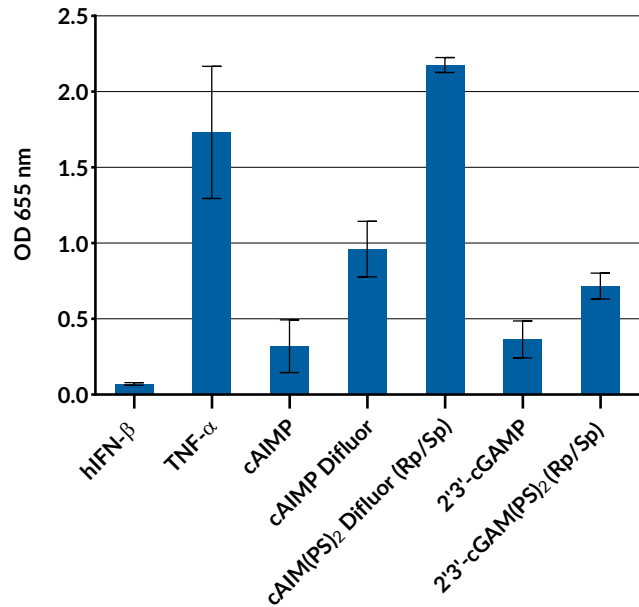
Version # 17107-MM

cAIMP and its analogs are novel STING-activating cyclic dinucleotides (CDNs) that induce NF- κ B and interferon (IFN) signaling. The activity of these CDNs has been compared to the reference agonist for human STING, 2'3'-cGAMP (see figures 1 & 2) using the THP1-Dual™ cells, a cell line derived from the human THP-1 monocyte cell line by stable integration of two inducible reporter constructs. As a result, these cells allow the simultaneous study of the IFN regulatory factor (IRF) pathway, by assessing the activity of a secreted Lucia luciferase, and the NF- κ B pathway, by monitoring the activity of SEAP. Interestingly, cAIMP exhibits similar potency to 2'3'-cGAMP, while the difluorinated (cAIMP Difluor) and bisphosphorothioate Difluor (cAIM(PS)₂ Difluor (Rp/Sp)) analogs are more potent than cAIMP and 2'3'-cGAMP.

IRF INDUCTION (Lucia luciferase reporter)



NF- κ B INDUCTION (SEAP reporter)



THP1-Dual™ cells were stimulated for 24 hours with human IFN- β (1×10^4 U/ml), TNF- α (300 pg/ml), cAIMP, cAIMP Difluor, cAIM(PS)₂ Difluor (Rp/Sp), 2'3'-cGAMP and 2'3'-cGAM(PS)₂ (Rp/Sp). All CDNs were used at 10 μ g/ml.

Figure 1: IRF induction was determined by measuring the relative light units (RLUs) in a luminometer using QUANTI-Luc™, a Lucia luciferase detection reagent. The IRF induction of each ligand is expressed relative to that of hIFN- β at 1×10^4 U/ml (taken as 100%).

Figure 2: NF- κ B induction was determined using QUANTI-Blue™, a SEAP detection reagent, and by reading the optical density (OD) at 655 nm. TNF- α has been included as a positive control to activate the NF- κ B signaling pathway.

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 3-622-34-80
E-mail: info@invivogen.com