

# Validation data for HEK-Blue™ STAT6-hSTING-R232 Cells

<https://www.invivogen.com/sting-r232-stat6-cells>

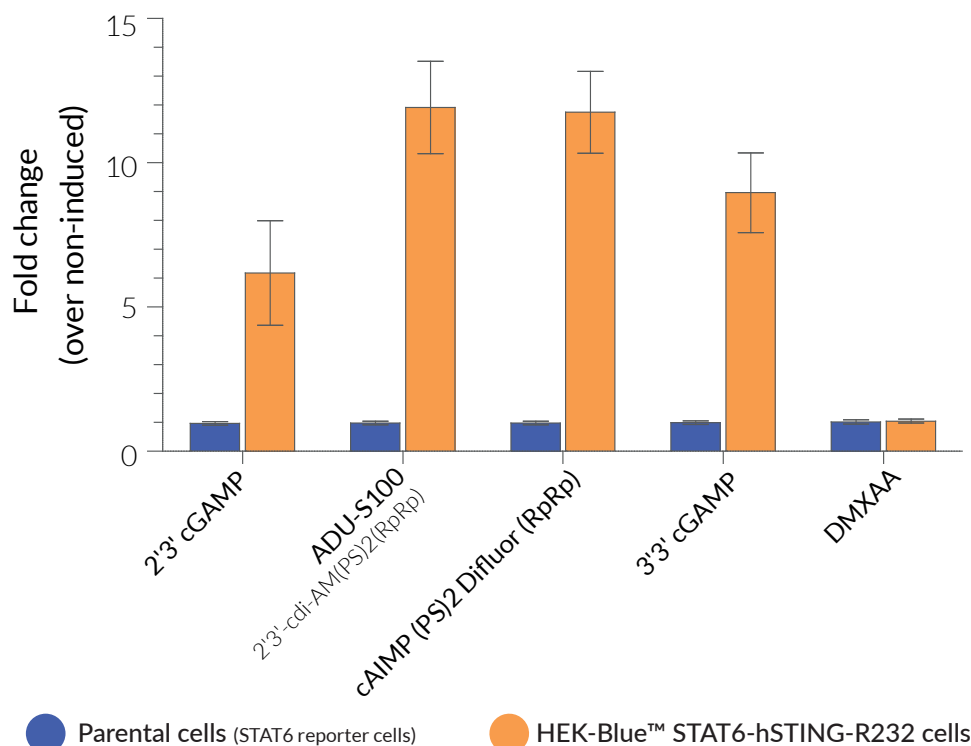
For research use only

Version 20J09-ED

HEK-Blue™ STAT6-hSTING-R232 cells were generated by stable transfection and overexpression of the wild-type human (h)STING R232 variant in a human embryonic kidney (HEK)293-derived cell line that expresses human STAT6. Additionally, these cells feature a STAT6-inducible SEAP (secreted embryonic alkaline phosphatase) reporter to monitor the activation of the STAT6 signaling pathway (Figure 1).

## Functional validation of STAT6 activation by hSTING (R232 variant)

Activation of the STAT6 signaling pathway in response to known STING ligands, such as cyclic dinucleotides (CDNs), has been assessed in HEK-Blue™ STAT6-hSTING-R232 cells. These cells exhibit a robust STAT6-dependent response to a range of STING agonists compared to the parental cell line. As expected, these cells exhibit no detectable response to the specific murine STING ligand DMXAA.



**Figure 1: STAT6 activation upon STING induction.** HEK-Blue™ STAT6-hSTING-R232 and their parental cells (STAT6 reporter cells) were incubated with 2'3'-cGAMP (30 µg/ml), 2'3'-cGAM(PS)2 (Rp/SP) (ADU-S100; 30 µg/ml), cAIMP(PS)2 Difluor (Rp/Sp) (30 µg/ml), 3'3'-cGAMP (30 µg/ml), or DMXAA (30 µg/ml) in cell culture medium. After overnight incubation, the STAT6 response was assessed by measuring the activity of SEAP in the supernatant using QUANTI-Blue™ Solution, a SEAP detection reagent. Data are presented fold change over non-induced cells (mean ± 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 Hong Kong: +852 3-622-34-80

E-mail: [info@invivogen.com](mailto:info@invivogen.com)