

PRODUCT INFORMATION



STING A162 variant (human, recombinant)

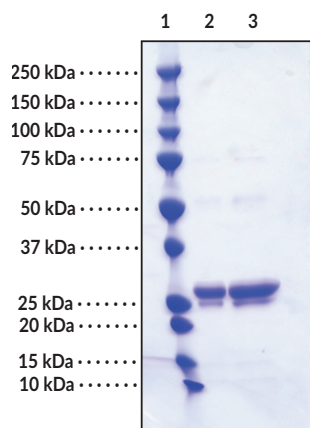
Item No. 25306

Overview and Properties

Source:	N-terminal histidine-tagged human recombinant STING R232, S162A mutant expressed in <i>E. coli</i>
Amino Acids:	138-379
Uniprot No.:	Q86WV6
Molecular Weight:	28.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥2 years
Purity:	batch specific (≥75% estimated by SDS-PAGE)
Supplied in:	50 mM HEPES, pH 8.0, 150 mM sodium chloride, 1 mM DTT, and 10% glycerol
Protein Concentration:	batch specific mg/ml
Activity:	Serial dilutions of canonical 3'3'-cGAMP were incubated with 5 µg recombinant human STING A162 variant in 50 mM HEPES, pH 7.5, 150 mM NaCl, 10% glycerol, and SYPRO® Orange dye at 4°C. ¹ The reaction was read on a BioRad CFX96 Touch™ Real-Time PCR Detection System at 4-100°C. ² The binding of the ligand stabilizes the protein structure, increasing the melting temperature (T_m), which is detected via a thermal shift assay (TSA), also known as a differential scanning fluorimetry (DSF) assay. ²

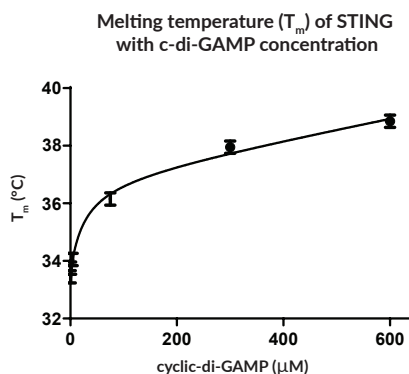
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: STING A162 variant (2 µg)
Lane 3: STING A162 variant (4 µg)

Representative gel image shown; actual purity may vary between each batch.



Binding Activity of STING A162 variant (Item No. 25306). STING A162 variant (5 µg) was incubated with serial dilutions of 3'3'-cGAMP (Item No. 17966) and SYPRO® Orange dye. The detected increase in T_m indicates binding.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/27/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

STING A162 variant (human recombinant) contains amino acids 138-379 of the wild-type variant (R232) with an alanine substituted for serine at position 162. Stimulator of interferon genes (STING) is a component of the innate immune response that binds to cyclic dinucleotides, which are bacterial second messengers, leading to activation of NF- κ B and transcription of immunomodulatory genes, including type I interferon (IFN).³⁻⁶ The R232 variant of STING is the most common variant in the human population, found at a frequency of 57.9% in the 1000 Genome Project.⁷ The SNP variant H232 is found at a 13.7% frequency. The S162A point mutation is located in the cyclic dinucleotide binding site of the human STING variants R232 and H232 and allows human STING to bind to DMXAA, a compound previously known to bind mouse, but not human, STING.^{8,9} When STING S162A is bound to DMXAA, it adopts the closed conformation, similar to the conformation it has when bound to the second messenger 2'3'-cGAMP (Item No. 19887), and activates the IFN pathway similarly to mouse STING.⁶

References

1. Diner, E.J., Burdette, D.L., Wilson, S.C., *et al.* The innate immune DNA sensor cGAS produces a noncanonical cyclic dinucleotide that activates human STING. *Cell Rep.* **3(5)**, 1355-1361 (2013).
2. Niesen, F.H., Berglund, H., and Vedadi, M. The use of differential scanning fluorimetry to detect ligand interactions that promote protein stability. *Nat. Protoc.* **2(9)**, 2212-2221 (2007).
3. Sun, L., Wu, J., Du, F., *et al.* Cyclic GMP-AMP synthase is a cytosolic DNA sensor that activates the type I interferon pathway. *Science* **339(6121)**, 786-791 (2013).
4. Wu, J., Sun, L., Chen, X., *et al.* Cyclic GMP-AMP is an endogenous second messenger in innate immune signaling by cytosolic DNA. *Science* **339(6121)**, 826-830 (2013).
5. Konno, H., Konno, K., and Barber, G.N. Cyclic dinucleotides trigger ULK1 (ATG1) phosphorylation of STING to prevent sustained innate immune signaling. *Cell* **155(3)**, 688-698 (2013).
6. Burdette, D.L., Monroe, K.M., Sotelo-Troha, K., *et al.* STING is a direct innate immune sensor of cyclic-di-GMP. *Nature* **478(7370)**, 515-518 (2011).
7. Yi, G., Brendel, V.P., Shu, C., *et al.* Single nucleotide polymorphisms of human STING can affect innate immune response to cyclic dinucleotides. *PLoS One* **8(10)**, e77846 (2013).
8. Gao, P., Ascano, M., Zillinger, T., *et al.* Structure-function analysis of STING activation by c[G(2',5')pA(3',5')p] and targeting by antiviral DMXAA. *Cell* **154(4)**, 748-762 (2013).
9. Conlon, J., Burdette, D.L., Sharma, S., *et al.* Mouse, but not human STING, binds and signals in response to the vascular disrupting agent 5,6-dimethylxanthenone-4-acetic acid. *J. Immunol.* **190(10)**, 5216-5225 (2013).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM