

PRODUCT INFORMATION



STING R232 variant (N242A mutant; human, recombinant)

Item No. 40241

Overview and Properties

Synonyms: Endoplasmic Reticulum Interferon Stimulator, ERIS, Mediator of IRF3 Activation, MITA, Mitochondrial Mediator of IRF3 Activation, MPYS, Stimulator of Interferon Genes, Stimulator of Interferon Response cGAMP Interactor 1, STING1, N-Terminal Methionine-Proline-Tyrosine-Serine Plasma Membrane Tetraspanner, TMEM173, Transmembrane Protein 173

Source: Active recombinant human N-terminal His-tagged STING R232 variant (N242A mutant) expressed in *E. coli*

Amino Acids: 138-379

Uniprot No.: Q86WV6

Molecular Weight: 28.7 kDa

Storage: -80°C (as supplied)

Stability: ≥1 year

Purity: *batch specific* (≥80% estimated by SDS-PAGE)

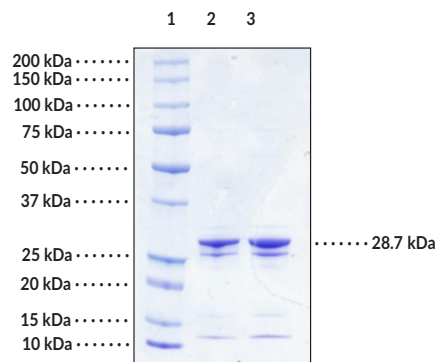
Supplied in: 50 mM HEPES, pH 8.0, with 100 mM sodium chloride, and 10% glycerol

Protein

Concentration: *batch specific* mg/ml

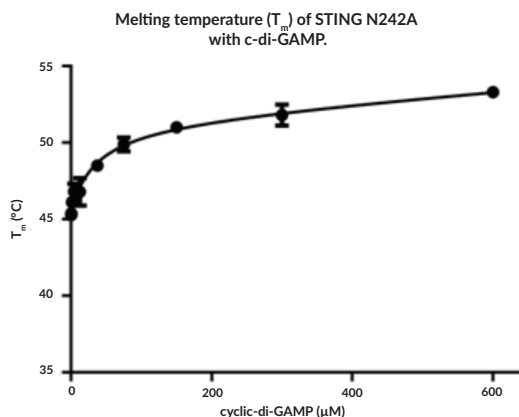
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: STING R232 variant (N242A mutant) (2 µg)
Lane 3: STING R232 variant (N242A mutant) (4 µg)

SDS-PAGE Analysis of STING R232 variant (N242A mutant).
This protein has an apparent molecular weight of 28.7 kDa.



Binding activity of STING N242A variant. STING N242A variant (5 µg) was incubated with serial dilutions of 3'3' cGAMP (Item No. 17966) and SPYRO™ 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, 05/29/2024

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

Stimulator of interferon genes (STING) is a component of the innate immune response that recognizes and binds to cyclic dinucleotides (CDNs), which either originate from bacteria or are signals of intracellular stress, leading to activation of NF- κ B and transcription of immunomodulatory genes, including type I interferon (IFN).¹⁻⁵ The R232 variant is the most common variant in the human population, found at a frequency of 57.9% in the 1000 Genome Project.⁶ STING is composed of four transmembrane domains at the N-terminus, a helix α 1 domain involved in protein dimerization and ligand sensing, and a cytoplasmic C-terminal domain containing the cyclic dinucleotide-binding domain, as well as the TBK1/IRF1-binding site, TBK1 phosphorylation site, and IRF3 docking site.⁷ Various mutations in STING either reduce or increase its activity or binding affinity.^{6,8-10} The asparagine-to-alanine substitution at position 242 (N242A) reduces c-di-GMP and 2'3'-cGAMP binding to STING.^{11,12} Cayman's STING R232 variant (N242A mutant; human, recombinant) protein can be used for ELISA, enzyme activity assay, and Western blot applications.

References

1. 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).
2. 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).
3. 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).
4. 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).
5. Hopfner, K.-P. and Hornung, V. Molecular mechanisms and cellular functions of cGAS-STING signalling. *Nat. Rev. Mol. Cell. Biol.* **21**(9), 501-521 (2020).
6. 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).
7. Li, Y., Wilson, H.L., and Kiss-Toth, E. Regulating STING in health and disease. *J. Inflamm. (Lond)* **14**, 11 (2017).
8. Jeremiah, N., Neven, B., Gentili, M., *et al.* Inherited STING-activating mutation underlies a familial inflammatory syndrome with lupus-like manifestations. *J. Clin. Invest.* **124**(12), 5516-5520 (2014).
9. Ni, G., Konno, H. and Barber, G.N. Ubiquitination of STING at lysine 224 controls IRF3 activation. *Sci. Immunol.* **2**(11), eaah7119 (2017).
10. Sauer, J.D., Sotelo-Troha, K., von Moltke, J., *et al.* The N-ethyl-N-nitrosourea-induced *Goldenticket* mouse mutant reveals an essential function of Sting in the *in vivo* interferon response to *Listeria monocytogenes* and cyclic dinucleotides. *Infect. Immun.* **79**(2), 688-694 (2011).
11. Shu, C., Yi, G., Watts, T., *et al.* Structure of STING bound to cyclic di-GMP reveals the mechanism of cyclic dinucleotide recognition by the immune system. *Nat. Struct. Mol. Biol.* **19**(7), 722-724 (2012).
12. 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).