

PRODUCT INFORMATION



cGAS (161-522) (human recombinant)

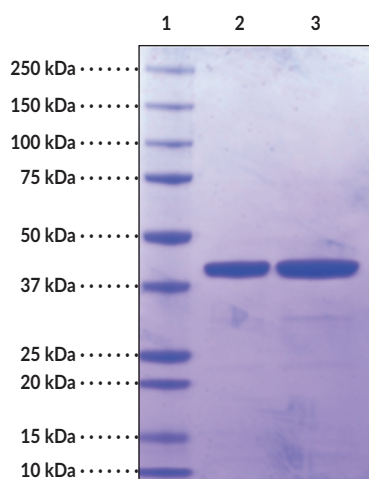
Item No. 25001

Overview and Properties

Synonyms: C6orf150, 2'3'-cGAMP Synthase, cGAMP Synthase, h-cGAS, Cyclic GMP-AMP Synthase, Mab-21 domain-containing protein 1, MB21D1
Source: N-terminal histidine-tagged human cGAS protein (truncated), purified from *E. coli*
Amino Acids: 161-522
Uniprot No.: Q8N884
Molecular Weight: 44.6 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: *batch specific* (≥90% 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

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

Image



Lane 1: MW Markers
Lane 2: cGAS (161-522) (2 µg)
Lane 3: cGAS (161-522) (4 µg)

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

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, 07/31/2018

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

Cyclic GMP-AMP (cGAMP) synthase (cGAS) (161-522) is a truncated form of cGAS (Item No. 22810) that contains the nucleotidyltransferase and Mab21 domains.¹ cGAS is a nucleotidyltransferase located in the cytosol that acts as a cytosolic DNA sensor to detect foreign DNA from microbial pathogens as part of the innate immune response.^{2,3} Upon binding to cytosolic DNA, cGAS produces the cyclic dinucleotide second messenger cGAMP, which activates stimulator of interferon genes (STING), leading to activation of the type I interferon (IFN) pathway.²⁻⁴ *In vitro*, fibroblasts, macrophages, and dendritic cells isolated from cGAS knockout (cGAS^{-/-}) mice do not produce type I IFNs following DNA transfection or DNA virus infection.⁵ Similarly, cells containing a frame-shift mutation in the cGAS locus fail to mount an immune response to HIV and other retroviruses.⁶ *In vivo*, cGAS^{-/-} mice infected with herpes simplex virus 1 (HSV-1) have lower levels of IFN- α and IFN- β , shorter survival times, and higher post-mortem levels of HSV-1 in the brain.⁵

References

1. Tao, J., Zhang, X.-W., Jin, J., *et al.* Nonspecific DNA Binding of cGAS N Terminus Promotes cGAS Activation. *J. Immunol.* **198**(9), 3627-3636 (2017).
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. Ablasser, A., Goldeck, M., Cavlar, T., *et al.* cGAS produces a 2'-5'-linked cyclic dinucleotide second messenger that activates STING. *Nature* **498**(7454), 380-384 (2013).
5. Li, X.-D., Wu, J., Gao, D., *et al.* Pivotal roles of cGAS-cGAMP signaling in antiviral defense and immune adjuvant effects. *Science* **341**(6152), 1390-1394 (2013).
6. Gao, D., Wu, J., Wu, Y.-T., *et al.* Cyclic GMP-AMP synthase is an innate immune sensor of HIV and other retroviruses. *Science* **341**(6148), 903-906 (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