



Development and Validation of a 2'3'-cGAMP ELISA

Daniel Tew, Tatyana Maxey, and Yogan Khatri

Cayman Chemical

— Developed in collaboration with **BIOLOG** - LIFE SCIENCE INSTITUTE -



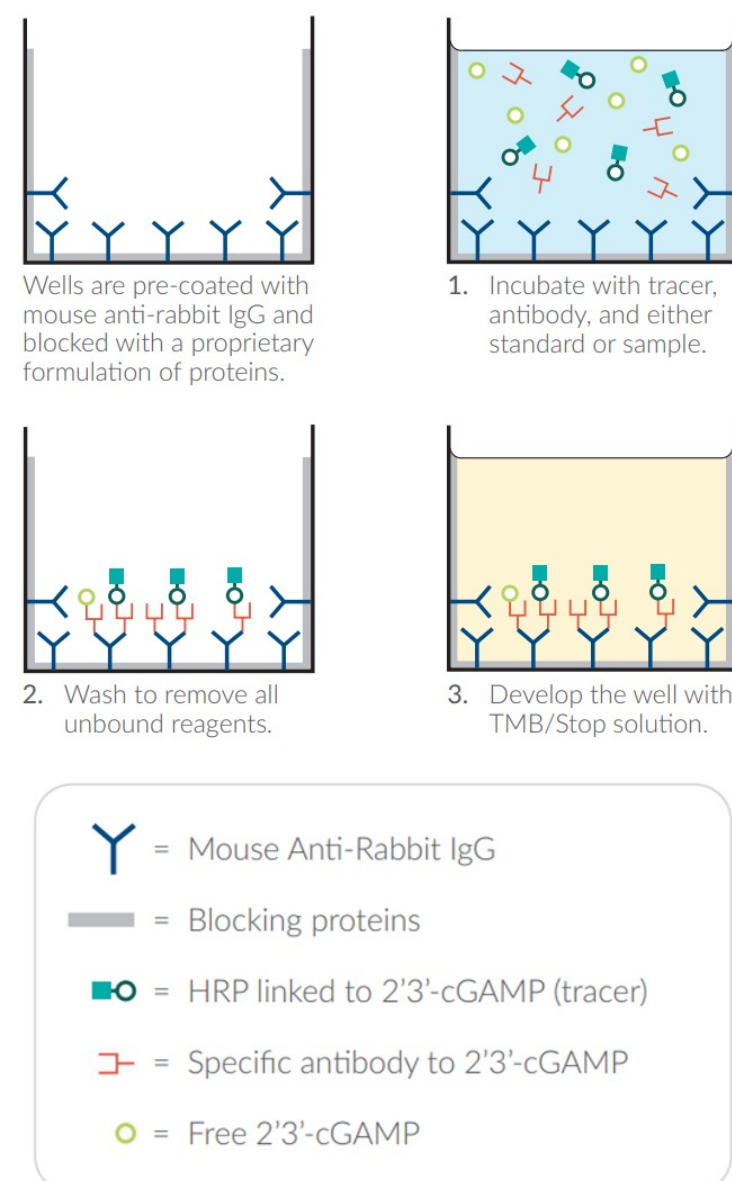
Scan to hear full ePoster presentation

WWW.CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

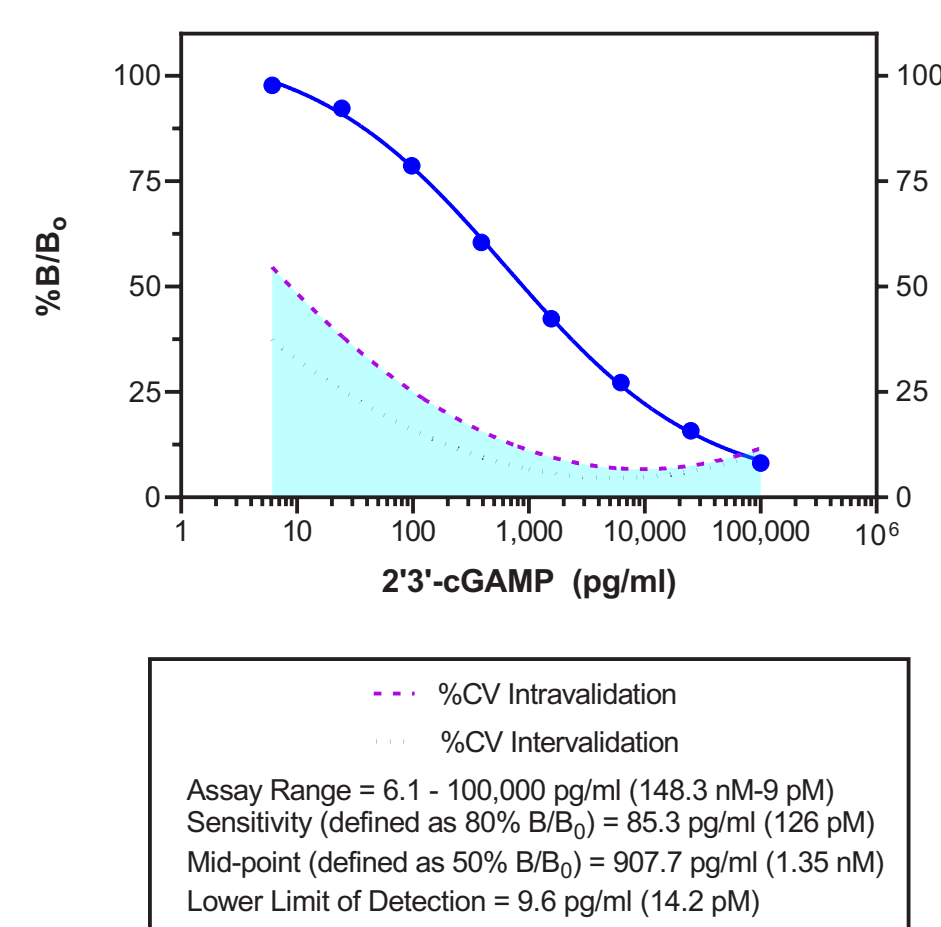
ELISA schematic



SPECIFICITY CHARACTERISTICS	
Compound	Cross Reactivity
2'3'-cGAMP	100%
2'2'-cGAMP	0.8%
3'3'-cGAMP	<0.01%
cyclic di-AMP	<0.01%
cyclic di-GMP	<0.01%
cGMP	<0.01%
cAMP	<0.01%
ATP	<0.01%
GTP	<0.01%

WWW.CAYMANCHEM.COM

2'3'-cGAMP ELISA Validation & Performance Characteristics



Intra-assay precision was determined by analyzing 24 replicates of three matrix controls (spiked M-PER™ samples) in a single assay.

Matrix Control (pg/ml)	%CV
7,165	8.4
828	14.5
102	21.3

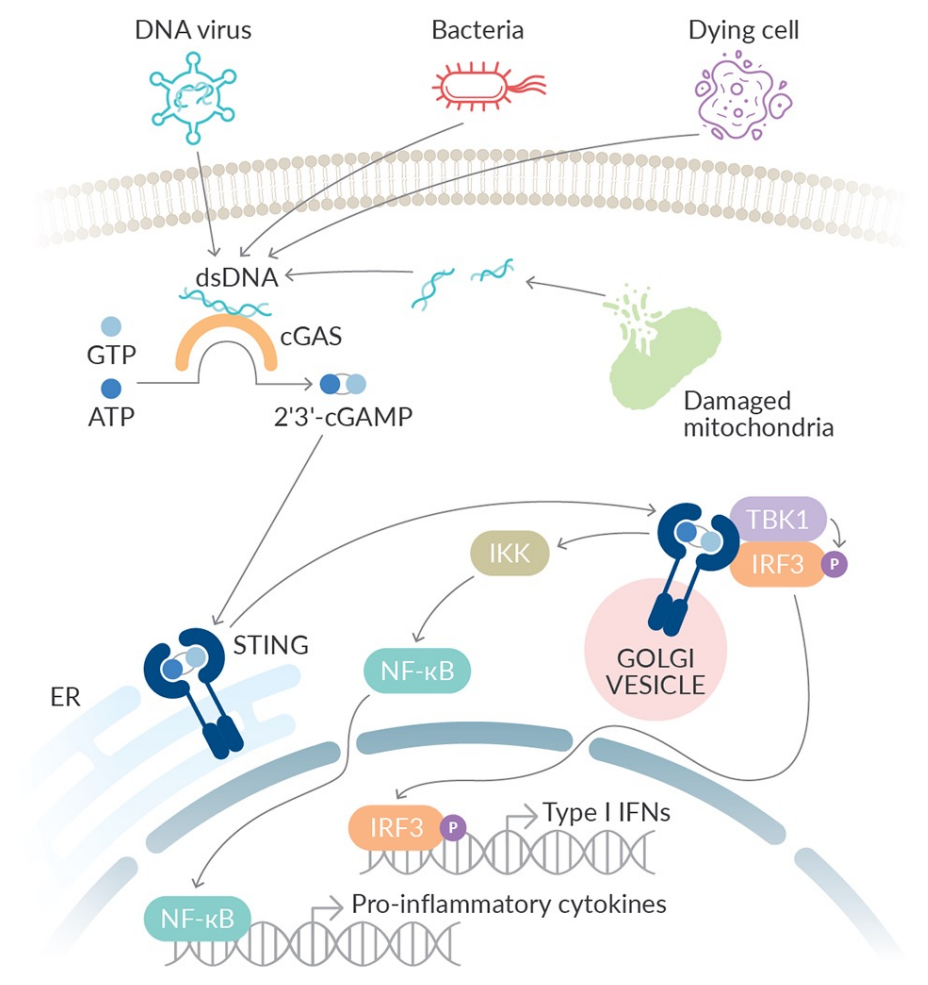
Inter-assay precision was determined by analyzing replicates of three matrix controls (spiked M-PER™ samples) in separate assays over several days.

Matrix Control (pg/ml)	%CV
7,530	9.6
811	9.6
72.2	18.7

WWW.CAYMANCHEM.COM

cGAS Pathway

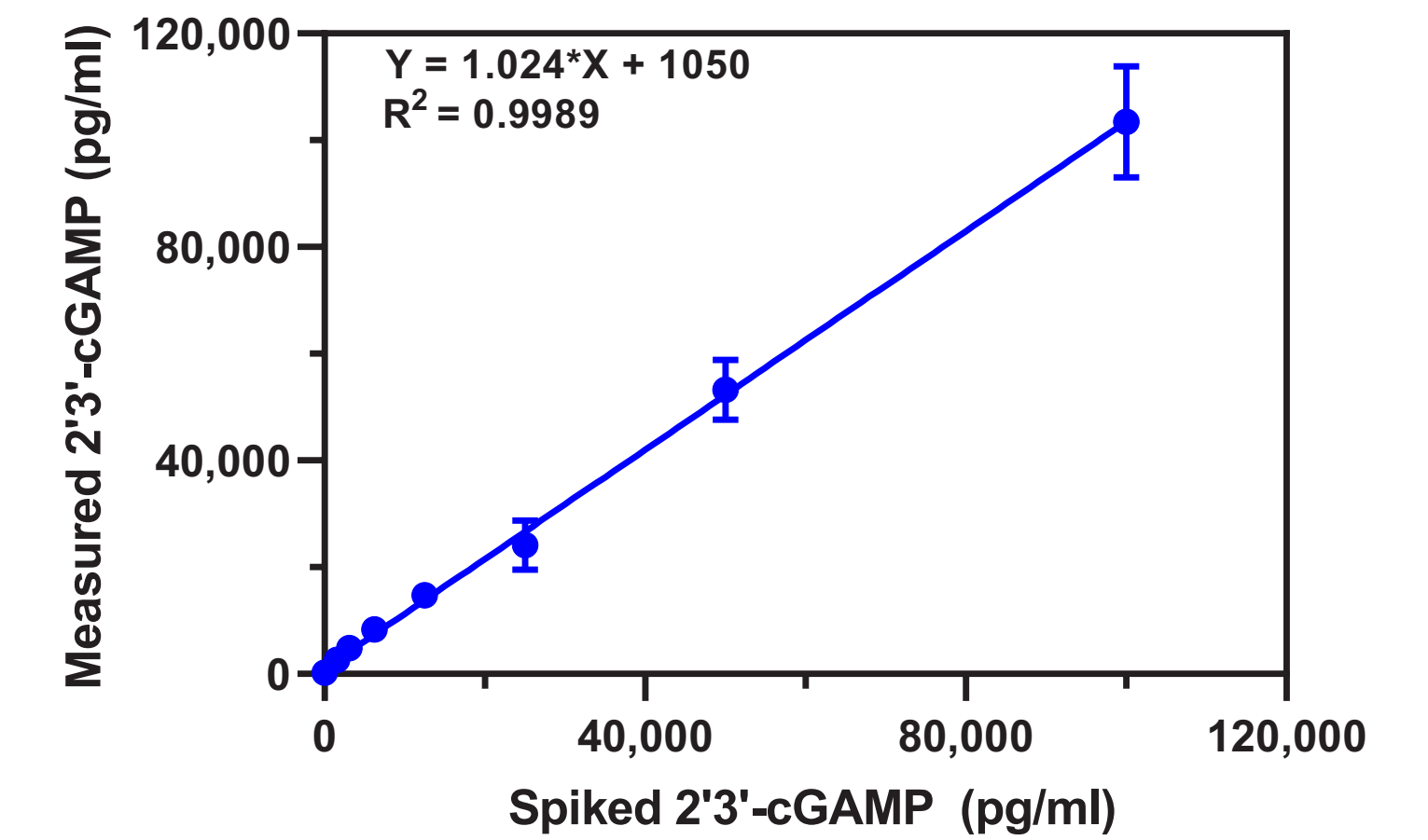
The cGAS-STING pathway is activated by an exogenous (DNA/RNA virus or bacteria/dead cells/tumor cell) or an endogenous (mtDNA released by stressed mitochondria) agonistic DNA. cGAS, consisting of a two-lobed catalytic domain and an extended N-terminal domain, synthesizes the noncanonically linked dinucleotide 2'3'-cGAMP in which DNA stands sandwiched between the two cGAS promoters recognizing their own DNA thread during its active and stable confirmation. The second messenger 2'3'-cGAMP binds to stimulator of interferon genes (STING) dimers residing at the endoplasmic reticulum (ER) and translocates to ER-Golgi intermediate compartments, leading to STING activation.



Upon activation, TANK-binding kinase-1 (TBK1) phosphorylates the CTT of STING and recruits interferon regulatory factors (IRFs) for phosphorylation, IRF dimerization, and nuclear translocation. This leads to the activation of downstream signaling and the transcription of target genes, including type I IFNs and ISGs (IFN-stimulated genes), in addition to NF-κB transcription factors.

WWW.CAYMANCHEM.COM

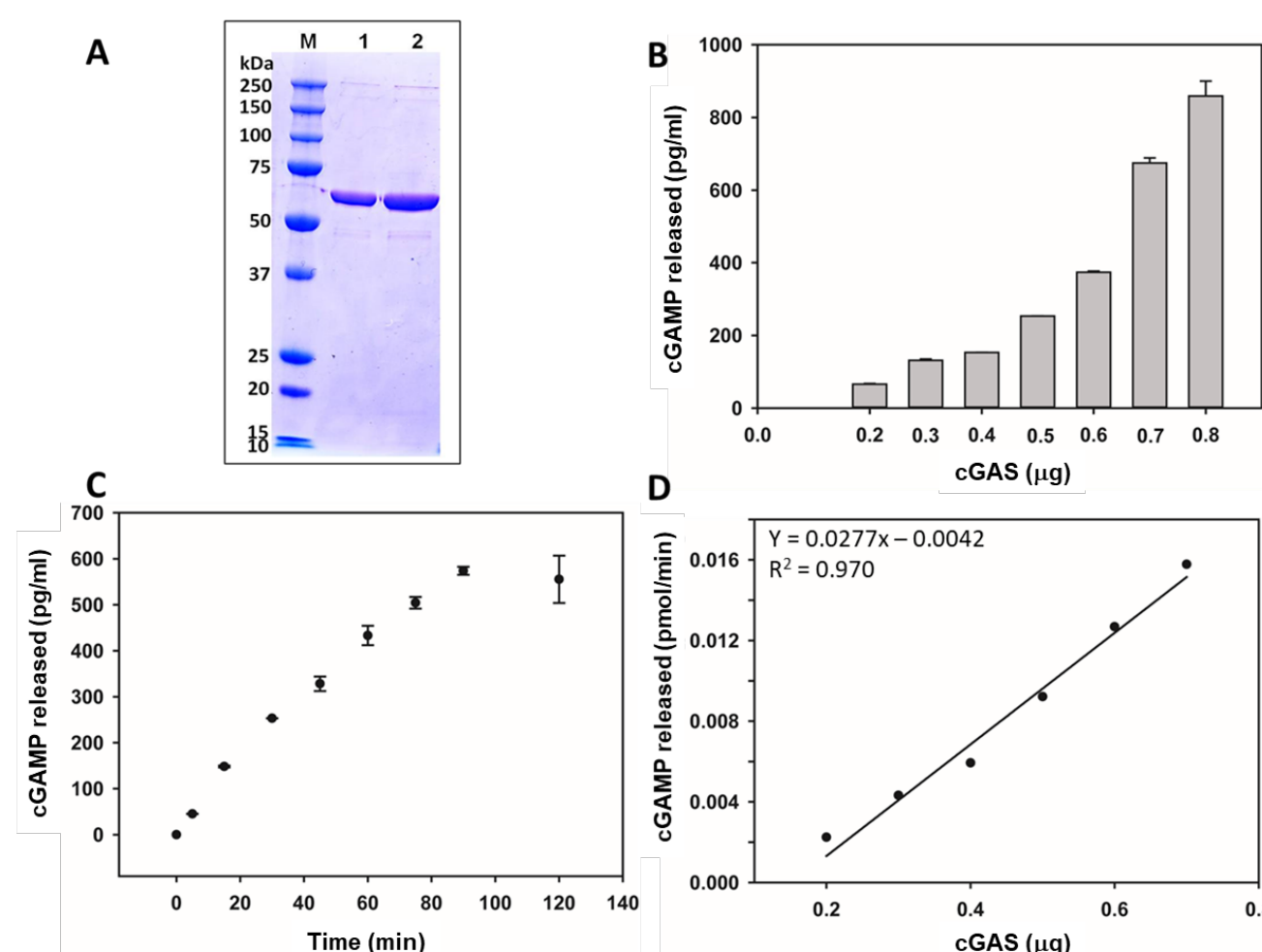
Spike & Recovery of 2'3'-cGAMP in THP-1 Cell Lysates



THP-1 cell lysate in M-PER™ (Thermo Scientific) was spiked with different amounts of 2'3'-cGAMP, serially diluted with Immunoassay Buffer C, and analyzed using the 2'3'-cGAMP ELISA. The error bars represent standard deviations from multiple dilutions of each sample.

WWW.CAYMANCHEM.COM

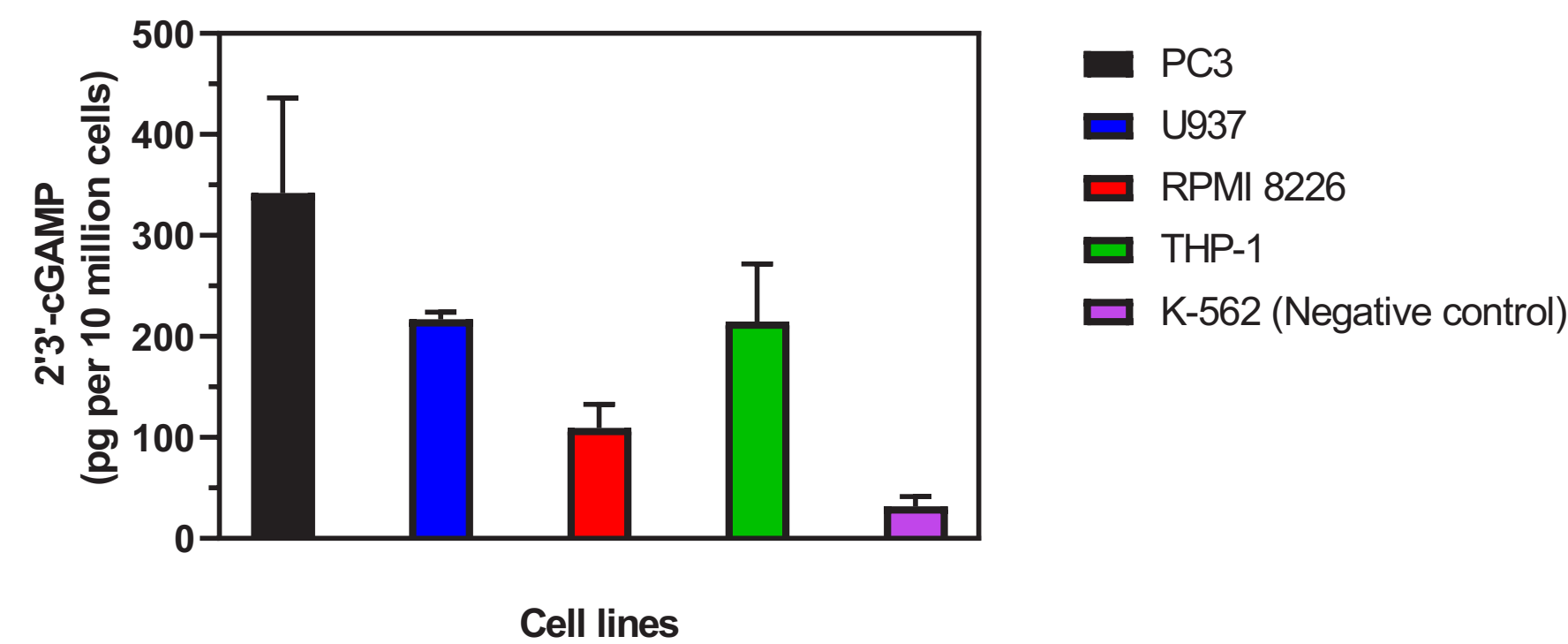
cGAS Characterization



Analysis of active human cGAS enzyme (Item No 22810). (A) Recombinant N-terminal His-tagged cGAS (full length, 2-522) expressed in *E. coli*. Lane 1- Marker, 2 and 3 - 2 and 4 μg of purified cGAS, respectively. Measurement of cGAMP released by (B) increasing concentration of cGAS enzyme assayed for 30 min and (C) increasing time period of cGAMP using 0.5 μM cGAS (D) The specific activity of cGAS was measured by plotting the rate of released cGAMP vs individual concentration of cGAS using 2'3'-cGAMP ELISA Kit (Item 501700). One unit of cGAS produces 1 pmole of cGAMP per minute at 37°C in 80 mM Tris-HCl, pH 7.5, containing 200 mM NaCl, 20 μM ZnCl₂, and 20 mM MgCl₂ with 4 μg circular DNA (3623 bp) and 0.25 mM each of GTP and ATP

WWW.CAYMANCHEM.COM

Quantification of 2'3'-cGAMP in Cell Lysates by ELISA



PC3 cells were cultured in F-12K media; RPMI 8226, U937, and THP-1 cells were cultured in RPMI 1640 media; K562 cells were cultured in IMDM media. All media contained 10% fetal bovine serum and 1% polysaccharide. Cells were lysed with M-PER™ (Thermo Scientific) at a concentration of 25 million cells per 2 ml lysis buffer. Clarified cell lysates were quantified using the 2'3'-cGAMP ELISA (Item No. 501700).

WWW.CAYMANCHEM.COM

Conclusions

The 2'3'-cGAMP ELISA is a sensitive and selective assay for measuring a wide range of 2'3'-cGAMP concentrations in cell-based and *in vitro* experiments, allowing researchers to identify cGAS modulating compounds. This assay is a valuable tool to monitor 2'3'-cGAMP released from the cGAS-STING pathway.

Acknowledgements

A special thanks to Laura Kostrzewa (Protein Core, Cayman Chemical Company) & Bill Ho (Cell Biology, Cayman Chemical Company) for supplying the cGAS protein and cells.

References

- 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).
- Hall, J., Braut, A., Vincent, F., *et al.* Discovery of PF-06928215 as a high affinity inhibitor of cGAS enabled by novel fluorescence polarization assay. *PLoS One* **12**(9), e0184843 (2017).
- O'Neill, L.A.J. Sensing the dark side of DNA. *Science* **339**(6121), 763-764 (2013).
- Cai, X., Chiu, Y.-H., and Chen, Z.J. The cGAS-cGAMP-STING pathway of cytosolic DNA sensing and signaling. *Mol. Cell* **54**(2), 289-296 (2014).
- Hopfner, K.-P. and Hornung, V. Molecular mechanisms and cellular functions of cGAS-STING signaling. *Nat. Rev. Mol. Cell Biol.* **21**(9), 501-521 (2020).

WWW.CAYMANCHEM.COM