

# Bacterial Signaling & Sensing

Cayman offers a wide range of products to study bacterial signaling and sensing by cyclic dinucleotide (CDN) second messengers and multiple quorum-sensing systems.



## Cyclic Dinucleotides (CDNs)

Bacteria use CDN second messengers, including 3'3'-cGAMP, 3'2'-cGAMP, cyclic di-GMP, and cyclic di-AMP, to regulate exopolysaccharide (EPS) synthesis, biofilm formation, antiphage defense, potassium homeostasis, cellular osmolarity, motility, and more. CDNs from pathogenic bacteria can be detected by the host protein stimulator of interferon genes (STING) during infection, alerting host cells to the presence of a pathogen and leading to activation of innate immune responses.

### ELISAs to Quantify Bacterial CDNs

Powered by **bioLog**  LIFE SCIENCE INSTITUTE

#### Cyclic di-GMP ELISA Kit

Item No. 501780

- Measure cyclic di-GMP in bacterial cell lysates
- LLOD: 7.7 pM (5.3 pg/ml)

#### Cyclic di-AMP ELISA Kit

Item No. 501960

- Measure cyclic di-AMP in bacterial cell lysates
- LLOD: 31.4 pM (20.7 pg/ml)

#### 3'3'-cGAMP ELISA Kit

Item No. 502130

- Measure 3'3'-cGAMP in bacterial and mammalian cell lysates and cell supernatants
- LLOD: 26 pM (17.5 pg/ml)

#### 3'2'-cGAMP ELISA Kit

Item No. 502340

- Measure 3'2'-cGAMP in fruit fly homogenates and bacterial lysates
- LLOD: 4.0 pM (2.7 pg/ml)

## Quantification of Cyclic Dinucleotides During Growth Phase in Three Bacterial Species Using ELISAs

Learn how Cayman's ELISA kits were used to measure and compare the levels of 3'3'-cGAMP, cyclic di-GMP, and cyclic di-AMP in bacterial lysates and supernatants.



Download the application note at [www.caymanchem.com/bacterialCDNs](http://www.caymanchem.com/bacterialCDNs)

## CDNs, Metabolites, & Negative Controls

Item No.	Product Name
17966	3'3'-cGAMP (sodium salt)
17753	Cyclic di-AMP (sodium salt)
17144	Cyclic di-GMP (sodium salt)
22485	Cyclic di-IMP (sodium salt)
33890	Cyclic di-UMP (sodium salt)
33886	5'-pApA (sodium salt)
33889	5'-pGpG (sodium salt)
36245	<sup>13</sup> C <sub>20</sub> , <sup>15</sup> N <sub>10</sub> -Cyclic di-GMP (sodium salt)

## Microbial Evasion of STING Signaling

In *Mycobacterium tuberculosis*, cyclic di-AMP is hydrolyzed by cyclic di-NMP phosphodiesterase (CdnP) to prevent activation of host STING and induction of innate immune responses. Cayman offers purified recombinant CdnP as well as the CdnP inhibitor C82 to study this enzyme.

Item No.	Product Name
22809	CdnP ( <i>Mycobacterium tuberculosis</i> strain ATCC 25618/H37Rv recombinant)
33563	C82

See all STING signaling products at [www.caymanchem.com](http://www.caymanchem.com)

# Quorum Sensing

Bacteria utilize quorum-sensing systems to control population-dependent expression of a variety of genes, including those involved in bioluminescence and virulence factor production. Quorum sensing signaling molecules, also called autoinducers, include acyl-homoserine lactones in Gram-negative bacteria, the *Pseudomonas* quinolone signal (PQS) 2-heptyl-3-hydroxy-4(1H)-quinolone and other quinolones in *P. aeruginosa*, and a class of cyclic dipeptides called diketopiperazines.

## Acyl-homoserine Lactones (AHLs)

Item No.	Product Name
10011207	N-( $\beta$ -ketocaproyl)-L-Homoserine lactone
9001147	N-3-hydroxydecanoyl-L-Homoserine lactone
10007895	N-3-oxo-dodecanoyl-L-Homoserine lactone
10011206	N-3-oxo-octanoyl-L-Homoserine lactone
10007898	N-butyryl-L-Homoserine lactone
10011201	N-decanoyl-L-Homoserine lactone
10011203	N-dodecanoyl-L-Homoserine lactone
10007896	N-hexanoyl-L-Homoserine lactone
10011199	N-octanoyl-L-Homoserine lactone
10011200	N-tetradecanoyl-L-Homoserine lactone

See all homoserine lactones at [www.caymanchem.com](http://www.caymanchem.com)

## Quinolones

Item No.	Product Name
29186	2-heptyl-3-hydroxy-4(1H)-Quinolone
9002699	2-nonyl-3-hydroxy-4-Quinolone
9003969	2-nonyl-3-hydroxy-4-Quinolone-d <sub>4</sub>

## Diketopiperazines

Item No.	Product Name
25739	Cyclo( $\Delta$ -Ala-L-Val)
34758	Cyclo(Phe-Pro)
27961	Cyclo(L-Pro-L-Val)
24943	Cyclo(L-Phe-L-Val)

See all microbial diketopiperazines at [www.caymanchem.com](http://www.caymanchem.com)

## Quorum Sensing Inhibitors

Item No.	Product Name	Description
27483	Ambuic Acid	A cyclohexanone that inhibits the biosynthesis of cyclic peptides involved in quorum sensing, including GBAP in <i>E. faecalis</i> , AIP-I in <i>S. aureus</i> , and LsrD698 and LsrD826 in <i>L. innocua</i>
31777	Amicoumacin B	Inhibits violacein production and downregulates the <i>viaA</i> , <i>viaB</i> , <i>viaD</i> , and <i>viaE</i> genes in <i>C. violaceum</i>
14198	10Z-Hymenialdisine	A natural marine sponge alkaloid that inhibits bacterial quorum sensing at micromolar concentrations
14016	Iberin	Inhibits AHL signaling in <i>P. aeruginosa</i> without affecting growth (IC <sub>50</sub> = 31-62 $\mu$ M)
28407	10-Methyldodec-2-en-4-olide	A bacterial volatile lactone that inhibits induction of an AHL-responsive reporter by the AHL N-dodecanoyl-L-homoserine lactone in a <i>P. putida</i> AHL biosensor strain
18163	Nifuroxazide	A nitrofurantoin antibiotic that inhibits quorum sensing and virulence factor production in <i>P. aeruginosa</i>
11440	Penicillic Acid	A mycotoxin that inhibits quorum sensing in <i>P. aeruginosa</i> by selectively repressing various virulence factor and other quorum sensing-regulated genes
36609	Qstatin	Inhibits activation of the quorum-sensing regulator LuxR in a $\beta$ -galactosidase reporter assay (25 $\mu$ M)

View a complete list of quorum sensing products  
at [www.caymanchem.com](http://www.caymanchem.com)

## Discover Our Latest Bacterial Signaling & Sensing News

Read about the synthesis and degradation of bacterial CDNs, their role in biofilm formation, and available research tools and assay kits at

[www.caymanchem.com/biofilms](http://www.caymanchem.com/biofilms)

