

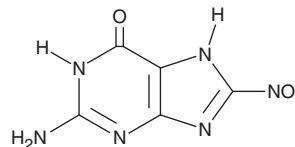
# PRODUCT INFORMATION



## 8-Nitroguanine

Item No. 89295

CAS Registry No.: 168701-80-2  
Formal Name: 2-amino-1,9-dihydro-8-nitro-6H-purin-6-one  
MF:  $C_5H_4N_6O_3$   
FW: 196.1  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{\max}$ : 397 nm  
Supplied as: A solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

8-Nitroguanine is supplied as a solid. A stock solution may be made by dissolving the 8-nitroguanine in the solvent of choice. 8-Nitroguanine is soluble in 1 M sodium hydroxide at a concentration of approximately 50 mg/ml.

### Description

8-Nitroguanine is a nitrative guanine derivative formed by oxidative damage to the guanine base in DNA by reactive nitrogen species (RNS) during inflammation and *in vitro* by reaction of DNA with peroxynitrite and other RNS reagents.<sup>1-3</sup> It is mutagenic and induces G:C to T:A transversion in DNA. Incorporation of 8-nitroguanine as an 8-nitroG:anti-G base pair into a primer template stalls human DNA polymerase  $\beta$  and induces a 2:1 preference for deoxyadenosine (dA) insertion over deoxycytosine (dC).<sup>3</sup> 8-Nitroguanine levels are increased in the lung tissue of mice with conditional expression of mutant K-Ras(G12V) that developed lung adenocarcinoma and in the lung tissue and peripheral lymphocyte DNA of cigarette smoke-exposed rats.<sup>4,5</sup> It is also increased in the lung tissue of influenza- or Sendai virus-infected mice and the colon epithelial cells of mice in an inflammatory bowel disease (IBD) model.<sup>6,7</sup>

### References

1. Hiraku, Y. Formation of 8-nitroguanine, a nitrative DNA lesion, in inflammation-related carcinogenesis and its significance. *Environ. Health Prev. Med.* **15**(2), 63-72 (2010).
2. Ohshima, H., Sawa, T., and Akaike, T. 8-nitroguanine, a product of nitrative DNA damage caused by reactive nitrogen species: Formation, occurrence, and implications in inflammation and carcinogenesis. *Antioxid. Redox Signal.* **8**(5-6), 1033-1045 (2006).
3. Bhamra, I., Compagnone-Post, P., O'Neil, I.A., et al. Base-pairing preferences, physicochemical properties and mutational behaviour of the DNA lesion 8-nitroguanine. *Nucleic Acids Res.* **40**(21), 11126-11138 (2012).
4. Ohnishi, S., Saito, H., Suzuki, N., et al. Nitrative and oxidative DNA damage caused by K-ras mutation in mice. *Biochem. Biophys. Res. Commun.* **413**(2), 236-240 (2011).
5. Hsieh, Y.-S., Chen, B.-C., Shiow, S.-J., et al. Formation of 8-nitroguanine in tobacco cigarette smokers and in tobacco smoke-exposed Wistar rats. *Chem. Biol. Interact.* **140**(1), 67-80 (2002).
6. Akaike, T., Okamoto, S., Sawa, T., et al. 8-nitroguanosine formation in viral pneumonia and its implication for pathogenesis. *Proc. Natl. Acad. Sci. U.S.A.* **100**(2), 685-690 (2003).
7. Ding, X., Hiraku, Y., Ma, N., et al. Inducible nitric oxide synthase-dependent DNA damage in mouse model of inflammatory bowel disease. *Cancer Sci.* **96**(3), 157-163 (2005).

#### 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, 11/09/2022

#### 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