

# PRODUCT INFORMATION



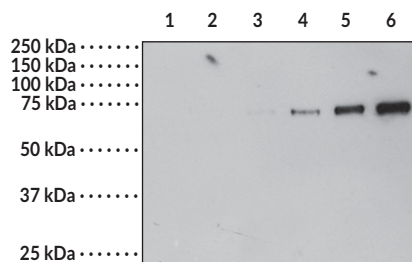
## Nitrotyrosine Monoclonal Antibody (Clone CC.22.8C7.3)

Item No. 189542

### Overview and Properties

<b>Contents:</b>	This vial contains 50 µg or 200 µg of protein G-purified IgG.
<b>Synonym:</b>	3-Nitrotyrosine, NT
<b>Immunogen:</b>	Peroxynitrite-treated KLH
<b>Cross Reactivity:</b>	(+) Chlorotyrosine (weakly)
<b>Species Reactivity:</b>	Species Independent
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Stability:</b>	≥3 years
<b>Clone:</b>	CC.22.8C7.3
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG2b
<b>Applications:</b>	ELISA and Western blot (WB); the recommended starting concentration for ELISA is 2 µg - 15 ng/ml and 2 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Lane 1: BSA (1 µg)  
Lane 2: Nitrotyrosine BSA (0.01 µg)  
Lane 3: Nitrotyrosine BSA (0.025 µg)  
Lane 4: Nitrotyrosine BSA (0.05 µg)  
Lane 5: Nitrotyrosine BSA (0.1 µg)  
Lane 6: Nitrotyrosine BSA (0.2 µg)

**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, 08/16/2024

**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

---

Nitrotyrosine is a post-translational modification that is formed by the nitration of tyrosine.<sup>1</sup> Under conditions of oxidative stress, tyrosine is oxidized by reactive oxygen species (ROS) or, in the presence of hydrogen peroxide and nitrite, by myeloperoxidase (MPO) to yield a tyrosine radical that reacts with reactive nitrogen species (RNS), such as nitric oxide or peroxynitrite, to form nitrotyrosine. It exists in a free or protein-bound form and is commonly used as a marker of nitrosative or oxidative stress.<sup>2</sup> Nitrotyrosine residues have been found in a variety of proteins, including LDL, surfactant protein A, angiotensin II, and human and bovine serum albumin.<sup>3</sup> Free nitrotyrosine production induced by peroxynitrite is inhibited by the polyphenols epicatechin gallate, gallic acid, catechin, or epicatechin in cell-free assays, as well as in aortic rings isolated from normotensive and spontaneously hypertensive rats administered the antioxidant N-acetyl-cysteine (NAC; Item No. 20261).<sup>4,5</sup> Nitrotyrosine levels are increased in the affected tissues of numerous pathological conditions, including atherosclerosis, cancer, ulcerative colitis, Alzheimer's disease, and Parkinson's disease.<sup>3</sup> Autoantibodies that recognize nitrotyrosinated proteins are increased in the synovium of patients with rheumatoid arthritis and are positively correlated with joint and tendon inflammation.<sup>1</sup> Cayman's Nitrotyrosine Monoclonal Antibody (Clone CC.22.8C7.3) can be used for ELISA and Western blot (WB) applications.

## References

---

1. Smallwood, M.J., Nissim, A., Knight, A.R., *et al.* Oxidative stress in autoimmune rheumatic diseases. *Free Radic. Biol. Med.* **125**, 3-14 (2018).
2. Teixeira, D., Fernandes, R., Prudêncio, C., *et al.* 3-Nitrotyrosine quantification methods: Current concepts and future challenges. *Biochimie* **125**, 1-11 (2016).
3. Oldreive, C. and Rice-Evans, C. The mechanisms for nitration and nitrotyrosine formation *in vitro* and *in vivo*: Impact of diet. *Free Radic. Res.* **35(3)**, 215-231 (2001).
4. Pannala, A.S., Rice-Evans, C.A., Halliwell, B., *et al.* Inhibition of peroxynitrite-mediated tyrosine nitration by catechin polyphenols. *Biochem. Biophys. Res. Commun.* **232(1)**, 164-168 (1997).
5. Cabassi, A., Dumont, E.C., Girouard, H., *et al.* Effects of chronic N-acetylcysteine treatment on the actions of peroxynitrite on aortic vascular reactivity in hypertensive rats. *J. Hypertens.* **19(7)**, 1233-1244 (2001).

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