

**Purified Mouse Anti-species independent
Nitrotyrosine Antibody *EM-30,
monoclonal***Catalog number: V1031855
Unit size: 0.1 mg**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Species independent
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Nitrotyrosine
Clone	EM-30

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, IHC(P), IHC(F)

Applications

3-Nitro-L-tyrosine is a 3-nitrotyrosine comprising L-tyrosine having a nitro group at the 3-position on the phenyl ring. It has a variety of applications in the study of important biological activities such as oxidase activity, intrinsic activity and dismutase activity, especially in cell-lines like plant cells, host cells and cultured cells. 3-Nitrotyrosine has been widely used in the examination of important cell-based pathways such as signaling pathway and serotonin pathway. It has a variety of applications in the examination of essential cellular processes like protein synthesis, SPase-catalyzed hydrolysis and chemical basis. It is essential for the examination of vital cellular enzymes like ceramide glycanases, tryptophan-specific permease and tyrosine dehalogenase