

**Purified Mouse Anti-human β -Endorphin
Antibody *B31.15, monoclonal***Catalog number: V103215
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	β -Endorphin
Clone	B31.15

Biological Properties

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

Applications

beta-Endorphin has shown use in a multitude of applications, specifically in the field of biochemistry and biology. It is vital to the examination of vital biological activities like neuronal activity, modifying activity and physiological activity, particularly in cell-lines such as intercalated cells, epidermal keratinocytes and SH-SY5Y cells. It is essential for the research of vital biological receptors such as ghrelin receptors, opioid receptors and glucocorticoid receptors. Moreover, beta-Endorphin is essential for the application of critical cellular pathways like signaling pathways and pioglitazone pathways, and it is essential for the application of vital biological enzymes like superoxide dismutase.