

---

## Nesfatin-1 Monoclonal Antibody [NESFATY1]

(Catalog No. A-0621)

---

### Background

Nesfatin-1 is metabolic polypeptide derived from the posttranslational processing of nucleobindin-2 (NUCB2). Originally recognized as a hypothalamic neuropeptide, it is also found in other parts of the brain, the pancreas, gastric endocrine cells and adipocytes. Involved in the regulation of hunger and fat storage, an increase in the Nesfatin-1 in the hypothalamus has been shown to suppress appetite and promote energy expenditure, independent of leptin signaling.

### Description

Mouse monoclonal antibody to Nesfatin-1, Clone NESFATY1

### Formulation

Liquid. In PBS containing 0.02% sodium azide

### Immunogen

Recombinant rat nesfatin-1 (aa 25-106)

### Specificity

Mouse, Rat

### Isotype

IgG1

### Purification

Protein G- affinity purified

### Storage

Store at +4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

### Application

ELISA, IP (1:100), WB (1:1000)

### Ordering Information

| Products                                  | Size   | Cat. No.   |
|---|--------|------------|
| Nesfatin-1 Monoclonal Antibody [NESFATY1] | 100 µg | A-0621-100 |

*This product is for research purposes only. Not intended for use in diagnostic procedures.*