
IL-33 Monoclonal Antibody [NESSY1], Biotin Conjugated

(Cat. No. A-0613)

Background

The protein encoded by this gene is a cytokine that binds to the IL1RL1/ST2 receptor. The encoded protein is involved in the maturation of Th2 cells and the activation of mast cells, basophils, eosinophils and natural killer cells. Several transcript variants encoding different isoforms have been found for this gene.

Description

Mouse monoclonal antibody to IL-33, Clone NESSY1 (biotin conjugated)

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Immunogen

Recombinant human IL-33 (aa 112-270)

Specificity

Human, Mouse, Rat

Uniprot ID:

O95760

Isotype

IgG1

Purification

Ascites purified.

Storage

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

Alternative Names

Interleukin-33, IL-1F11, NF-HEV

Application

ELISA

Ordering Information

Products	Size	Cat. No.
IL-33 Monoclonal Antibody [NESSY1], Biotin Conjugated	100 µg	A-0613-100

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