
MRJP1 Polyclonal Antibody, HRP Conjugated

(Catalog # A53729)

Background

Major royal jelly protein 1: induces the differentiation of honeybee larvae into queens through an Egfr-mediated signaling pathway. Promotes body size increase by activating p70 S6 kinase, stimulates ovary development by augmenting the titer of vitellogenin (Vg) and juvenile hormone, and reduces developmental time by increasing the activity of mitogen-activated protein kinase and inducing the 20-hydroxyecdysone protein (20E). Most abundant protein found in the royal jelly which is the food of the queen honey bee larva. The royal jelly determines the development of the young larvae and is responsible for the high reproductive ability of the honeybee queen

Description

MRJP1 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Apis mellifera

Isotype

IgG

Uniprot ID

O18330

Purification

>95%, Protein G purified

Immunogen

Recombinant Apis mellifera Major royal jelly protein 1 protein (20-432AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

56-kDa protein 4 Bee-milk protein Royalactin MRJP1

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
