
Amatoxin Polyclonal Antibody

(Catalog # A64220)

Background

Major toxin from Amanita phalloideae. Acts by binding non-competitively to RNA polymerase II and greatly slowing the elongation of transcripts from target promoters.

Description

Amatoxin Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Concentration

1 mg/ml

Dilution

1:500 to 1:2000 with PBS + 0.1% tween-20. Diluted antibody should be used within the assay day

Specificity

Amanita phalloides

Isotype

IgG

Uniprot ID

P85421

Purification

Protein G purified

Immunogen

Peptide sequence from Amanita phalloides Alpha-amanitin proprotein

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

Amatoxin, Alpha-amanitin, Gamma-amanitin

Application

ELISA
