
rodA Polyclonal Antibody, HRP Conjugated

(Catalog #A55803)

Background

Cell wall protein regularly arranged in interwoven fascicles of clustered proteinaceous microfibrils, or rodlets, to form the outer spore coat protein. It is involved in resistance to environmental stress and may well be associated with conidial hydrophobicity. It is important in the morphogenesis of the dispersible conidia.

Description

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

Formulation

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

Specificity

Neosartorya fumigata

Isotype

IgG

Uniprot ID

P41746

Purification

>95%, Protein G purified

Immunogen

Recombinant Neosartorya fumigata Hydrophobin protein (19-159AA)

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

Rodlet protein, rodA, hyp1, AFUA_5G09580

Application

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