
blaNDM-1 Polyclonal Antibody, FITC Conjugated

(Catalog A55776)

Background

Confers resistance to many beta-lactam antibiotics, including some carbapenems. Does not confer resistance to the polymyxin colistin or the fluoroquinolone ciprofloxacin.

Description

blaNDM-1 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Klebsiella pneumoniae

Isotype

IgG

Uniprot ID

C7C422

Purification

Protein G purified

Immunogen

Recombinant Klebsiella pneumoniae Metallo-beta-lactamase type 2 protein (29-270AA)

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

NDM-1, Metallo-beta-lactamase NDM-1, blaNDM-1

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
