

His-Tag Polyclonal Tag Antibody, HRP Conjugated

(Catalog # A26404)

Background

A polyhistidine-tag is an amino acid motif in proteins that consists of at least six histidine (His) residues, often at the N- or C-terminus of the protein. Polyhistidine-tags are often used for affinity purification of polyhistidine-tagged recombinant proteins that are expressed in Escherichia coli or other prokaryotic expression systems.

Description

Rabbit polyclonal antibody to His-tag, HRP conjugated

Purification

Affinity purified

Specificity

Species Independent

Isotype

IgG

Immunogen

This antibody is produced by immunizing rabbits with a synthetic peptide CHHHHHH(KLH-coupled).

Formulation

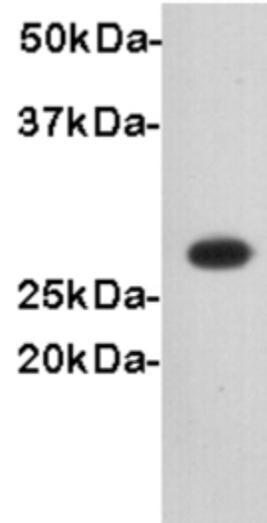
Liquid. 1*TBS (pH7.4), 0.5% BSA, 25% Glycerol. Preservative: 0.05% Sodium Azide.

Storage

Store at 4°C after thawing. Aliquot and store at -20°C or -80°C. Avoid repeated freeze/thaw cycles.

Application

WB: 1:2,000



▲ Western blot analysis on His-Tag recombinant protein (26kDa) using His-Tag Polyclonal Tag Antibody, HRP Conjugated.

Products

His-Tag Polyclonal Tag Antibody, HRP Conjugated

Size

100 µl

Cat. No.

A26404-100

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