
HA-TAG Tag Antibody

(Catalog No. A26004)

Description

HA-Tag Tag Antibody is specific for the HA (Haemagglutinin) epitope tag peptide sequence YPYDVPDYA, when fused to either the amino- or carboxyl- terminus of target proteins, including expression in many frequently used expression vectors. HA-TAG Tag Antibody has been tested against both the immunogen and recombinant proteins containing the HA sequence in ELISA and Western blotting. In Western blotting of bacterial extracts this antibody has been shown not to cross-react with any endogenous proteins.

Antibody Type

Rabbit Polyclonal Antibody

Purification

Purified IgG - liquid

Immunogen

Keyhole limpet hemocyanin conjugated epitope tag peptide YPYDVPDYA (114-122) from haemagglutinin influenza. A cysteine residue was used to facilitate coupling at the C-terminal end.

Isotype

Polyclonal IgG

Size

50 µg

Formulation

Phosphate buffered saline. 0.09% Sodium Azide (NaN₃)

Specificity

HA

Storage

Store at +4°C or at -20°C if preferred. Storage in frost-free freezers is not recommended. This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody.

Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

ELISA; Immunohistology - Paraffin (This product requires antigen retrieval using heat treatment prior to staining of paraffin embedded sections. Citrate buffer is recommended for this purpose); Western Blotting

Ordering Information:

Products	Size	Cat. No.
HA-TAG Tag Antibody	50 µg	A26004-1

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