
C-MYC Tag Antibody

(Catalog No. A26001)

Description

Mouse Anti C-MYC

Antibody Type

Monoclonal Antibody

Purification

Purified IgG - liquid

Immunogen

Synthetic peptide sequence AEEQKLISEEDLL

Isotype

IgG1

Size

100 µg

Formulation

Phosphate buffered saline. 0.09% Sodium Azide

Specificity

Human

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: 1/100 - 1/500; Flow Cytometry: (Use 10ul of the suggested working dilution to label 1x10⁶ cells in 100ul) Neat - 1/10; Immunohistology - Frozen; Immunohistology - Paraffin: (This product does not require protein digestion pre-treatment of paraffin sections prior to staining. This product does not require antigen retrieval using heat treatment prior to staining of paraffin sections). Where this antibody has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the antibody for use in their own systems using appropriate negative/positive controls.

Ordering Information:

Products	Size	Cat. No.
C-MYC Tag Antibody	100 µg	A26001-1

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