
c-Myc Monoclonal Antibody [9E10]

(Catalog No. A-0642)

Description

Mouse monoclonal antibody to c-Myc, Clone 9E10

Size

0.5 mg

Formulation

Liquid. In PBS.

Immunogen

Synthetic peptide corresponding to aa 408-438 of human pp62 c-myc.

Specificity

Human

Isotype

IgG1

Purification

Purified

Storage

Store at -80°C. Avoid multiple freeze/thaw cycles.

Application

IHC, IP, WB

Ordering Information

Products	Size	Cat. No.
c-Myc Monoclonal Antibody [9E10]	0.5 mg	A-0642-500

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