



CERTIFICATE OF ANALYSIS

Product ANTI-BIOTIN M
(MONOCLONAL ANTIBODY)

Catalog No. MB-9100

Amount 1 ml

Lot No. ZB0505

Produced in mouse

Clone Hyb-8

Class IgG1

Solution Reconstitute by adding 1 ml water. The resulting culture supernatant contains 0.08% sodium azide.

Storage Refrigerate lyophilized powder. After reconstitution, solution should be stored frozen.

This product is recommended for *in situ* hybridization or immunohistochemical applications. For these applications the concentrated stock solution should be diluted at least 1:200 for a standard 30 minute room-temperature incubation. Optimal dilution of the anti-biotin should be established for each application. This reagent can be detected with a VECTASTAIN[®] ABC System or the ImmPRESS[™] Anti-Mouse IgG Polymer Detection System (Cat. No. MP-7402).