



CERTIFICATE OF ANALYSIS

Product **BIOTINYLATED ANTI-MALTOSE BINDING PROTEIN**
AFFINITY PURIFIED

Catalog BA-0701

Amount 0.25 mg biotinylated IgG

Lot No. L1113

Concentration 1.0 mg/ml

Produced in goat

Solution 10 mM phosphate, 0.15 M NaCl, pH 7.8, 0.08% sodium azide.
Immunohistochemical grade bovine serum albumin has been added for stabilization.

Storage Refrigerate

Immunoglobulin biotinylated 99%

This biotinylated antibody is to be used in western blot detection of maltose binding protein (MBP) or MBP-containing fusion (chimeric) proteins. This product has been optimized for use with the VECTASTAIN[®] ABC-AmP[™] Kit (Cat. No. AK-6000). The recommended dilution for use is 1:1000 in 1x casein solution (10x Casein Solution; Cat. No. SP-5020).